





Cytobank **簡易マニュアル**

MAPSS-LS-202312-13

内	容
	_

1 クラウド型サイトメトリー解析ソフトウエア Cytobank について	7
1-1 Cytobankとは	8
1-2 必要要件	8
1-3 データセキュリティ	8
1-4 クリニカルデータの保存についての注意点	8
1-5 アカウント・ライセンス	9
1-5-1 購入ライセンスの種類	9
1-5-2 ライセンス終了後	9
1-5-3 アカウントアクセスレベル(ライセンスなし)でできること	9
1-5-4 アカウントのデータ容量	
1-6 利用規約とプライバシーポリシー	
1-7 お問い合わせ先	
1 - 8 Cytobank 日本語関連資料・ガイドビデオ	
1-9 Cytobank 英語資料	12
2 基本機能	13
2-1 アカウント作成(Premium cytobank)	14
2-2 ログイン画面(Experiment Manager)	15
2-2-1 ログイン方法	15
2-2-2 Experiment Manager	15
2-2-3 Experiment を開く	17
2-2-4 Labelを付与・解除する	17
2-2-5 Experiment のアーカイブ化	18
2-2-6 Project の作成とその管理	19
2-3 Experiment の新規作成・データアップロード	22
2-3-1 Experiment 画面	25
2-3-2 機能ツール一覧	25
2-4 Actions:	26
2-4-1 View Summary	26
2-4-2 Clone	
2-4-3 upload	31
2-4-4 Export	
2-4-5 Link to parent: Experimentのリンク付けと解除	
2-4-6 Reset:Experimentを初期化	
2-4-7 Delete: Experiment を削除	
2-5 スケール調整	

2-5-1 スケールの種類	38
2 - 5 - 2 スケール調整画面: Scale Editor	39
2-6 コンペンセーション(蛍光補正)	40
2-6-1 Compensation Editor	40
2-6-2 新規にマトリックスを作成する	40
2-6-3 マトリックスをインポートする	41
2-6-4 自動コンペンセーション	42
2 - 7 データ QC : PeacoQC	43
2-7-1 自動データ QC ツール: PeacoQC とは	43
2 - 7 - 2 PeacoQC 設定·実行	43
2 - 7 - 3 PeacoQC 結果の評価	46
2-7-4 実行上限の目安	49
2-8 ゲーティング	51
2 - 8 - 1 Cytobank でのゲーティングとポピュレーションの定義	51
2-8-2 サーバー保存のタイミングと Experiment gate 情報としての保存	51
2-8-3 Gating Editor:ゲーティング画面の各機能	51
2-8-4 プロットに表示されるイベント数が多すぎる場合:Time Slice	53
2-8-5 測定中の異常イベントのクリーンアップ	54
2-8-6 ゲーティング後のコンペンセーション/チャネル変更:Hidden gate	54
2-8-7 Boolean expression : ポピュレーションの定義と編集	56
2-8-8 Check gates:ゲート位置の一括確認	60
2-8-9 Gate tailoring : ファイルや親集団でのゲート位置調整	60
2-8-10 ポピュレーションの階層構造の確認と編集	66
2 - 8 - 1 1 ゲーティングストラテジーの自動表示: Gating hierarchy	69
2-8-12 ゲート情報のインポート	69
2-9 Sample tags:サンプルタグ	72
2-9-1 サンプルタグとは	72
2-9-2 サンプルタグのつけ方	72
2 - 9 - 3 他の Experiment からのサンプルタグのインポート	73
2-9-4 サンプルタグの一括編集・追加	74
2-9-5 チャネルとパネル設定: Assign panels	75
2-9-6 ファイル属性の管理	77
2 - 1 0 Figure 作成(Illustrations; Event level plot)	79
2 - 1 0 - 1 新規 illustrator の作成	79
2 - 1 0 - 2 Plots 設定	79
2 - 1 0 - 3 Layout 設定	85

2-10-4 ファイル統合プロットの表示: Virtual concatenation	
2 - 1 0 - 5 Save 設定	91
2 - 1 0 - 6 Export 設定	92
2-10-7 Templates ボタン	93
2-10-8 印刷ボタン	93
2-10-9 ロックボタン	93
2-10-10 プロット上でのレイアウト変更操作	94
2 - 1 0 - 1 1 Statistic table : プロット下の統計表	94
2-11 統計グラフと統計的有意差検定(Illustrations; summary plot)	96
2-11-1 統計グラフ設定の基本	96
2-11-2 Heatmap : ヒートマップ	
2-11-3 Box Plot:箱ひげ図	102
2-11-4 Violin plot:バイオリンプロット	103
2-11-5 Bar chart : 棒グラフ	105
2-11-6 Line chart & Stain index:線グラフとステインインデックス	107
2-11-7 Summary dots : サマリードット	111
2-11-8 統計的有意差検定機能(共通)	113
2-11-9 統計グラフのエクスポート	121
2-1 2 共有・プロジェクト・権限	121
2 - 1 2 - 1 ユーザーコネクション(Premium cytobank のみ)	121
2-12-2 データセキュリティ・デフォルトのアクセス権限について	122
2-12-3 Experimentの共有	122
2-12-4 アクセス権限を付加した共有ができるプロジェクトの作成と管理	123
3 Advanced analyses(機械学習解析)	126
3-1 機械学習による解析	127
3-2 自動ゲーティング	128
3-2-1 自動ゲーティングとは	128
3-2-2 トレーニングデータの準備	128
3-2-3 モデルトレーニングの設定・実行	129
3-2-4 新規データへのモデルの適用方法	132
3-2-5 自動ゲーティング(モデル作成/適用)の結果の評価	133
3-2-6 自動ゲーティング後の解析例	139
3-2-7 トレーニング上限と計算時間の目安、トラブルシューティング	141
3-3 次元削減アルゴリズム(tSNE-CUDA、UMAP、opt-SNE、viSNE)	145
3-3-1 軸スケールの調整(前処理)	145
3-3-2 ゲーティングによる解析対象の絞り込み(前処理)	146

3-3-3 設定と実行	146
3-3-4 複数のアルゴリズム結果の比較表示	158
3-3-5 結果のマップ上でのゲーティング	159
3-3-6 結果の良い/悪いの見方	161
3-3-7 結果の illustrations での表示と比較	162
3-4 FlowSOM	167
3-4-1 FlowSOM とは	167
3-4-2 tSNE/UMAP マップ上に FlowSOM メタクラスターを表示する場合	167
3-4-3 FlowSOM 設定・実行	167
3-4-4 FlowSOM 結果の出力ファイル	170
3-4-5 FlowSOM 結果の Cytobank への出力	172
3 - 5 SPADE	179
3-5-1 SPADEとは	179
3-5-2 SPADE 設定・実行	179
3-5-3 SPADE 結果の表示と解析	180
3-5-4 バブルの作成	182
3-5-5 選択したノードのプロット表示	182
3-5-6 バブル内の細胞(イベント)の出力	182
3-6 CITRUS	
3-6-1 CITRUSとは	
3-6-2 CITRUS設定·実行	
3-6-3 CITRUS 結果の確認	186
3-6-4 CITRUS 結果ファイル	187
3-6-5 CITRUS 結果ファイルの見方	190
3-6-6 CITRUS 結果から見出した候補クラスターの詳細解析・グラフ化	190
4 各種機能	192
4-1 ダウンロード・エクスポート	193
4 - 1 - 1 Experiment アーカイブ (acs) のダウンロード	193
4-1-2 統計計算結果・サンプルタブ・ゲート情報・Experiment内 FCS ファ	イル&添付ファイ
ル・Illustrations で表示している統計表・Illustrations で表示している細胞情	青報の FCS/text
ファイルのエクスポート	193
4-2 Experiment の複製	
4-2-1 Experimentの削除	
4-2-2 ゴミ箱に入れた Experiment を戻す・完全消去する	
4-3 FCS ファイル詳細情報	195
4-4 FCS ファイル以外のデータを Cytobank ヘインポート(DROP 機能)	195

	4-5 アカウントの確認・設定変更	196
	4-5-1 パスワードの変更	196
	4-5-2 使用データ容量と内訳の確認	197
	4-5-3 ユーザーコネクションのリクエストとリストの確認	197
	4-5-4 登録 E-mail アドレスの変更	198
	4-5-5 その他細かいアカウント設定の変更	198
	4-6 ログアウト	199
5	バージョン対応表	200

1 クラウド型サイトメトリー解析ソフトウエア Cytobank について

1-1 Cytobank とは

シングルセルデータ解析のためのクラウドプラットフォームのソフトウエアです。

非常に高いセキュリティのクラウド上に、データや解析結果、関連ファイルを保存および一元管理することが できます。

フローサイトメトリーのハイパラメータデータやその他シングルセルデータを、機械学習アルゴリズムをベースにした高次元解析ツールで簡単に解析することができます。

データや解析を世界中の共同研究者と、ハイセキュリティなクラウド上で安全にシェアすることができます。

1-2必要要件

・ウェブブラウザ(Chrome, Firefox, Safari 最新バージョン推奨)

・E-mail アドレス

・インターネット (通信速度 50 MBPS download/ 10 Mbps upload 以上推奨) Speedtest.net にて速度環境を調べることができます。

※Microsoft Edge には対応しておりません。

※ブラウザの動作するデバイスであれば端末のスペックは要求しませんが、操作性からパソコンでの使用を おすすめします。

※複数のブラウザやデバイスから重複してログインするとサーバーセキュリティが不審者からのアクセスと判断 し、アカウントを一時停止することがあります。

1-3データセキュリティ

Cytobank 上のデータは完全に安全な状態で保存されています。データと解析は暗号化され、定期的 にバックアップされています。

セキュリティの詳細について記載された文書が必要な方はお気軽にお問合せください。

1-4クリニカルデータの保存についての注意点

Cytobank は研究用途のみにお使いいただけ、診断にはご使用になれません。 クリニカルデータをアップロードする際には、匿名化する必要があります。 21 CFR Part 11 には対応しておりません。

1-5アカウント・ライセンス

Cytobank へのアクセスレベルは2つのレベルになっており、レベル1はアカウントアクセス、レベル2はサブ スクリプションアクセスです。アカウントアクセスレベルでは閲覧とエクスポートのみが可能、サブスクリプション アクセスではデータのアップロードや解析等すべての機能が可能です。アクセスレベルを上げるためには、トラ イアル/Enterprise/Premium のいずれかのライセンスが必要となります。

1-5-1購入ライセンスの種類

EnterpriseとPremiumの2種類のライセンスがあります。

Enterprise

- 独立した専用サーバー+ライセンス(10 ライセンス以上)
- サーバー内アカウント数は制限なしで、管理者権限によるライセンス管理(ライセンスの移動)が可能
- デフォルトデータ容量は5 TB/サーバー
- データ共有は同一サーバー内のみで、他のサーバー(Premium Cytobank 含む)とは共有・ア クセスできません。
- ラボ / 企業単位向け

■ Premium

- Premium Cytobank サーバー内のアカウント単位ライセンス
- 最小単位:1ライセンス
- デフォルトデータ容量は 500 GB/アカウント
- Premium Cytobank 内アカウントでのデータ共有が可能(Enterprise アカウントとは不可)

1-5-2ライセンス終了後

Enterprise ではサーバー内でライセンスが移動しただけですので、アカウントアクセスはデータ・解析結果 とともにサーバー内に維持されます。

Premium ではライセンス終了後 90 日間はアカウントとデータ・解析結果は維持されますが、複数回の メール通知の後、データは完全削除されます。

データが削除される前にデータやアーカイブファイル(.acs)などをエクスポート・ダウンロードください。

1-5-3アカウントアクセスレベル(ライセンスなし)でできること

サブスクリプションアクセスレベルの時に作成した Illustrations や figure を表示したり、生データやアーカ イブファイル (.acs) をダウンロードしたりすることができます。データを追加したり、解析したりすることはで きません。

1-5-4アカウントのデータ容量

Enterprise では契約時に購入したサーバーの容量(デフォルト 5 TB)を全アカウントで共有します。 Premium では 500 GB/アカウントです。

容量が不足しそうな場合、容量のみの追加購入も可能ですのでお気軽にお問合せください。

1-6利用規約とプライバシーポリシー

Cytobank の利用規約(Terms of Use)とプライバシーポリシー(Privacy Policy)はお使いの Cytobank ページ下部からいつでもご確認いただけます。



1-7お問い合わせ先

ベックマン・コールター株式会社 ■Cytobank 専用サポートメール cytobank_support@beckman.com

■電話(お客様サポートセンター) 0120-566-730 または 03-6745-4704 9:00~17:30 (土・日・祝日・その他休業日を除く)

1-8 Cytobank 日本語関連資料・ガイドビデオ

■ Cytobank サーバーの日本語サポート Experiment

Cytobank サーバーの日本語サポート Experiment にアクセスすると最新の日本語マニュアルなどの情報を取得いただけます。 公開 Experiment なのでどなたでもアクセスできます。

Cytobank ログイン> Experiment manager の検索ボックスに"日本語"と入力> 日本語サポート Experiment にアクセス

Experiment Manager	9、日本語		Group linked			
• New experiment	Q Searc	h results				Settings
Q All 1170		Name	<i>V</i>	PR	Created	Updated ▼
Mine 190	0	診日本語サポート_Cytobank support for Japanese	1	Cytobank staff	Apr 12	Apr 13
& Shared with me 812						
Cytobank curated 27						
Public 143						

最新版マニュアルは Experiment summary の Attachments セクションからダウンロードいただけま

す	þ											
	日本語サポー = Actions	►_Cytobank supp Y Sample tags	ort for Japanese	🗘 Gates	តំរិ Advanced anal	yses 🚺	Illustrations		🐨 Scales	A Compensation	Sharin;	g lic
	? Experim	nent summary	Clone									
0 0 1	-	Name: Labels:	日本語サポート_C None + イ	Cytobank suppo	ort for Japanese		P Create Update	R: Cytobank st d: 6:24 pm, Apr 12 d: 06:10 PM	aff 2023			
1	 Notes Purpose: Cytobank日本ユーザー様向けサポート Comments: ■最新版マニュアル: Attachmentsセクションからダウンロードください。 ご質問等はsupport_cytobank@beckman.comまでご連絡ください。 											
	Lillust This exp	rations eriment does not conta eate an illustration	in any illustrations yet.									
	@ Attac	hments										
	File N	lame				Download	Date U	ploaded By	Size	md5sum		
	MAPS	SS-LS-202304-26_I	Premium_Cytobank	k_簡易マニュフ	7リレ_V4.000.pdf	B	Apr 13	Cytobank staff	17.2 MB	98f738e		

■ Cytobank 最初のガイドビデオ

Cytobank お使い始めのための4つのガイドビデオ

https://www.beckman.jp/flow-cytometry/software/cytobank-premium/learningcenter/basics

■ Cytobank ラーニングセンター

機能ごとの案内ページで、特徴の説明やガイド動画も豊富に用意されています。 https://www.beckman.jp/flow-cytometry/software/cytobank-premium/learning-<u>center</u>

■Cytobank 資料まとめページ

日本語ガイド動画やアプリケーションノート、テクニカルノートなど Cytobank の情報をまとめてご提供しております。

https://www.beckman.jp/flow-cytometry/software/cytobank-premium/jp-helpfullinks?web=1&wdLOR=cA8BE831A-4B24-41F4-9BB9-C1B18BB56EBC

1-9 Cytobank 英語資料

Cytobank support portal https://support.cytobank.org/hc/en-us

*本簡易マニュアルと製品マニュアル(英語版)の間に矛盾もしくは不一致が生じた場合は英語版が 優先するものとします。

2基本機能

2-1アカウント作成 (Premium cytobank)

ベックマン・コールターウェブサイトの申し込みフォーム(<u>https://www.beckman.jp/flow-</u> <u>cytometry/software/cytobank-premium/free-trial</u>)に登録をした後、Premium Cytobank サーバー(<u>https://premium.cytobank.org/cytobank/signup)</u>にてアカウントを 作成します。

フリートライアル申し込みフォーム	
	Register Now for Cytobank Premium
	Free and full-featured for 30 days - No credit card
	required!
	Username
17ドレス•	
	First name
26*	Last name
S.•	Contact email
	Confirm email
## F *	Use the address you want to receive Cytobank email (e.g. jane@companyx.com or john@institution.org)
(*	Privacy 🗹 Do not display my email address to other users
	Password:
	Confirm password:
用門•	Passwords should be 8 to 40 characters, with at least
a・サービス情報やWEBセミナー、展示会などのベックマン・コールターからのから	one letter and one number.
畑らせを希望しますか?	Pulsubmitting this form I confirm that I have reviewed and agree with th
ail : •	Privacy Policy and the Cytobank Terms of Use. I also understand my
参望する 参望しない	privacy choices as they pertain to my personal data as provided in the
*************************************	Privacy Policy Under Your Privacy Choices.
ere and - All Sector All And All Mark All Ma	
- 19 (A	Register

セキュリティのため、アカウントと申し込みフォームの情報が照合がされ、米国オペレーションチームによって認証されるとアカウントが登録され、認証メール(英文)が登録メールアドレスに届きます。認証メールのリンクをクリックすると、自動的に 30 日間の Cytobank premium トライアル(すべての機能を使用可能)が開始されます。

※認証作業は米国時間で行われるため、時差や祝日によって遅れる場合があります。3営業日を経過 しても認証メールが届かない場合は、メールソフトのゴミ箱や迷惑メールに入っていないかをご確認の上、 cytobank_support@beckman.comまでご連絡ください。

※Enterprise ではご所属のサーバーでのアカウント作成が必要となりますので、ご施設の管理者にお問 合せください。

2-2ログイン画面(Experiment Manager)

2-2-1ログイン方法

アカウント作成後は、下記 URL または Google で Cytobank と検索し、Cytobank ウェブサイトの左 上 Login> Premium からログインページに入ることができます。

Premium Cytobank : <u>https://premium.cytobank.org/</u>



アカウント登録した User name または Email と、Password を入力し Login ボタンを押します。 パスワードを忘れた場合は Login ボタン下の Forget password をクリックしてガイドに従います。

2 - 2 - 2 Experiment Manager

Cytobank では解析の単位(データセット)を Experiment として管理します。

ログインすると Experiment Manager 画面が表示されます。Experiment Manager はメール管理 ソフトのような構成で、アクセスできる Experiment の管理をすることができます。

アクセス可能な Experiment は、作成したものの他に共有しているものや公開されているものも含まれます。

	(1) Cytobank Prenium Experiment Mana	Exper	(4) iments	Pro	© ojects € © ⊈ Group link	ed			(12) Help	ynakane
		t 474	60	8	nk curated _{Name}	¥	PR	Size	Created	₀ [©] Settings Updated ▼
	Mine Mine	169			PBMC Experiment (fluorescence)	11	Qianjun Zhang	401.2 MB	Mar 7, 2013	
	A Shared with me	163			Dillustration Editor Example Data	8	Angela Landrigan	45.3 MB	Sep 2, 2020	L.
	Cytobank curated Public	20		1	© Healthy Human PBMC w/ 26 Surface Markers from Fluidigm	1	Angela Landrigan	182.0 MB	Jun 7, 2017	Jan 24
	Archived	3		2	© Fluidigm Basic Human PBMC Panel	1	Angela Landrigan	73.3 MB	Feb 13, 2013	Jan 21
\bigcirc	> Projects			2	© Science, Bendall et al (Mass Cytometry)	3	Angela Landrigan	12.3 GB	Jan 11	Jan 12
3	PRs				D U937 Experiment (Fluorescence)	15	Angela Landrigan	15.6 MB	May 21, 2014	Jan 2
_	Advanced analyses			1	© Cytek Aurora - 23 color immunophenotyping (FlowSOM demo)	3	Angela Landrigan	518.9 MB	Nov 9, 2018	Dec 28, 2020
	FlowSOM	26		4	D Fluidigm Cytobank Human PBMC Analysis Guide	2	Chris Ciccolella	670.8 MB	Sep 17, 2013	Dec 11, 2020
	visne	126	3	1	DITRUS Demo: BCR-XL stimulation of human PBMCs	16	Angela Landrigan	254.3 MB	Oct 18, 2016	Dec 2, 2020
	SPADE	27	A	1	~					
	🕨 🖤 Labels		(9)	Y	10					
	- Demo	16	C	N						
	V & A	16 4								

- Cytobank ロゴマーク:どの操作画面からでも Experiment Manager へ行くことができます。Ctrl キー(Windows)や Cmd(Mac)キーを押しながらクリックすると、ブラウザの機能により別タブで 開くことができます。
- ② New Experiment ボタン: <u>新規 Experiment 作成</u>を開始します。
- ③ カテゴリーセクション: Experiment を所有者ごとや Project ごと、Project Reader ごと、解析種類ごと、Label ごとに自動的または任意に分類して呼び出すことができます。
 各カテゴリー名の右の数字はそのカテゴリーに含まれる Experiment 数を示しています。数字部分をクリックすると Star this filter と hide this filter が表れます。
 Star this filter : 選択するとそのカテゴリーをリストの上部 All の下に移動します。
 Hide this filter : 選択するとリストの一番下に作成される Hidden filters へ格納します。
 - ・ All: アクセスできるすべての Experiment を中央のリストに表示します。
 - ・ Mine:ご自身が作成した(Clone を含む) Experiment のみを表示します。
 - Shared with me:他のアカウントの Experiment で、アクセス許可をしてもらった共有 Experiment のみを表示します。
 - ・ Cytobank curated: Cytobank が提供するデモデータ Experiment のみを表示します。
 - ・ Public: 公開されている Experiment のみを表示します。
 - ・ Trash: ゴミ箱に入れた Experiment のみを表示します。
 - Project: クリックすると作成・共有している Project が表示され各々の Experiment のみを 表示することができます。
 - PRs: クリックするとアクセスできる Experiment の Project Reader リストが表示され各々の Experiment のみを表示することができます。
 - Advanced analyses: クリックすると各種の解析が表示され、各々の解析結果
 Experiment のみを表示することができます。
 - Labels: クリックすると Label のリストが表示され、各ラベルを付与された Experiment だけ を表示することができます。
- ④ Experiments ボタン:新規作成や最近開いた Experiment へのリンクが開きます。
- S Project ボタン: Project の新規作成や全 Project の表示、最近使用した Project のリンクが開きます。
- ⑥ フィルターボックス:入力した文字でフィルターをかけ、目的の Experiment を見つけやすくします。
- ⑦ Group linked ボタン: チェックを入れると、解析結果の Experiment やクローンなどの関連
 Experiment をまとめて表示します。Linked Experiment も一覧に表示したい場合はチェックを 外します。
- ⑧ 選択チェックボックス: Experiment を選択すると画面上部にアクションボタンが表示されます。 Apply label ボタンで<u>ラベル付与</u>、Archive ボタンで<u>アーカイブ化</u>、Delete ボタンで<u>削除</u>をまとめて することができます。Unselect ボタンでは選択を解除することができます。

Cytobank Expe	riments Projects
Experiment Manager	Actions for 51 selected experiments 🖉 Apply label 🖀 Unarchive 🛱 Delete 🛞 Unselect
① New experiment	🖥 My experiments
Q All 474	☑ Name

- ⑨ 水色丸: Linked clone experiment (関連するクローンや Link to parent で紐づけられた Experiment) がある場合、その数が表示されています。
- ① Settings ボタン:中央の Experiment リストに表示する項目を設定することができます。
- 12 Helpボタン:サポートサイト(英語)のリンクなどが表示されます。
- ③ My profile ボタン:アカウント設定やログアウトができます。

2-2-3 Experiment を開く

- ① 開きたい Experiment 名の部分をクリックします。
- ② 右側にウィンドウが表れ、その Experiment の概要を確認することができます。
- ③ 上部水色の帯部分のメニューバーから使用したい機能を選択します(特にない場合は Experiment summary を開きます)。または、概要部分を下にスクロールして、Linked experiments や各 Advanced analysesの解析結果をクリックし、それらを直接開くこともできま す。



2-2-4 Label を付与・解除する

Experiment に任意のラベルをいくつでも付けることができ、様々な分類をすることができます。

- ① 対象の Experiment のチェックボックスにチェックを入れます。
- ② 上部にアクションボタンが表れますので、Apply label をクリックします。



③ 新規の Label を付与する場合は、Label name ボックスに名前を入力し、Create label and apply ボタンを押します。



- ④ Label が作成され付与されます。
- ⑤ Label の名前部分をクリックまたはカテゴリーセクションの Label をクリックすると、その Label が付与 された Experiment のみが表示されます。
- ⑥ 上部の Rename label ボタンで、Label の名前を変更することができます。
- ⑦ Edit color ボタンで、色を変更できます。
- ⑧ Delete ボタンで、その Label 自体を削除することができます。
- (9) Label の名前の右に X マークをクリックすると、その Experiment の Label を解除することができます。

Cytobank	Expe	riments Projects
Experiment Manage	er	Q label:demo
Onew experiment		'Demo' labeled Rename label Edit color Delete
Q AII	474	□ Name
🖬 Mine	169	4 1 0- Demo Gating illustrations Demo 🛞 < 🌖
🤽 Shared with me	163	
Cytobank curated	20	
Public	120	6 15 1-SPADE demo Demo
Archived	3	7 ElowSOM from viSNE recult (Clone) Demo
🛱 Trash	2	
🕨 🚞 Projects		
🕨 👤 PRs		
・ 品 Advanced analyses		
> @ Labels (5)		
🥏 Demo 🏏	16	

2-2-5 Experiment のアーカイブ化

しばらく使用しない Experiment をアーカイブフォルダに格納することができます。

① 対象の Experiment のチェックボックスにチェックを入れます。

- ② 上部にアクションボタンが表れますので、Archive をクリックします。
- ③ カテゴリーセクションの Public の下に Archived が作成され、これに格納されます。

2-2-6 Project の作成とその管理

複数の Experiment を Project としてまとめて管理することができます。 プロジェクトチームでの共同作 業にとても有用です。

Project は単に Experiment をまとめるだけでなく、ユーザーのアクセス権限も管理することができます。

- 2-2-6-1 Project の新規作成
- ① Experiment manager 画面上部の Projects ボタンをクリックし、New を選択します。



- ② 名前 Name (と必要であれば説明 Description)を入力します。
- ③ Project のアクセス権限を設定します。
 - Sharing level: illustrations 機能への権限を View Only(閲覧のみ)か
 illustrations(ユーザーが各々illustrations を作成できる)から選択します。
 - Cloning level: Clone (Experiment の複製)機能への権限を、Disabled (不可)、
 Clone Files Only (生データのみダウンロードや複製可能)、Full Clone (フルクローンが
 可能)から選択します。
- ④ Create project をクリックします。

૾ૢૺ	Cytobank	Experime	ents Projects		
<	New pro	oject			
		Name	TEST		
		Description			
	As you like				
					.11
		Sharing level	View only	~	
		Cloning level	Disabled	~	
			Create Proj	ect	

- ⑤ Edit project '名前'の画面が表示されます。
- ⑥ Edit Details 欄は前画面の設定が表示されます。 Manager はいつでも編集可能です。
- ⑦ Edit Project Members 欄で Project メンバーを追加します。メンバーの役割は Managers, Leaders と Members の 3 種類があります。 Managers は、メンバーや Experiment の追加、 Project のアクセス権限の設定変更などすべてのことができます。 Leaders と Members は名称は 異なりますが権限は同じで、 Project の設定(アクセス権限やメンバー)や Project 内の

Experiment (のゲートやスケール、コンペンセーションなどの設定を含む)を変更することはできません。また、Leaders と Members は③で設定されたアクセス権限 (Sharing level と Cloning level)の制限を受けます。

 ⑧ Add 役割名ボックスに Username またはアカウント登録されている E-mail アドレスを入力すると 候補ユーザーが表示されるので、選択して追加します。
 ※候補ユーザーは user connections リストが対象となります。 User connections に登録され ていないユーザーを追加したい場合は、ボックスに表示される Connect with a user をクリックし、 ユーザーコネクションリクエストを行ってください。



⑨ メンバーを削除する場合は名前の右側に表示される X マークをクリックします。

dit Project Members Managers										
Managers can add and remove project members, leaders, managers, Type and then click the account that appears										
Add Manager: darth vader 🗡										
Leaders Can access Add Leader: Darth Vader g to the sharing and										
Members										
Members can access project experiments according to the sharing an										
Geoff Kraker 2 Click to remove										
☑ Invite a new user										

- Project から Experiment を削除する場合は、左のボックスで Experiment を選択し、ボックス下の Remove Experiment(S)ボタンを押します。※Project から削除されるだけで、この操作で Experiment 自体は削除されません。

Add experiments current unassigned to projects:	Remove experiments from this project:
pbmc	Filter displayed:
(12100) - PBMC (Proper scales) (Clone) [12100) - PBMC (Proper scales) (Clone) [12307] - PBMC (Proper scales) (Clone) [12313] - PBMC (Proper scales) (Clone) [12313] - PBMC (Proper scales) (Clone) [12429] - PBMC (Proper scales) (Clone) [12433] - PBMC (Proper scales) (Clone) [12430] - PBMC (Proper scales) (Clone)	[3790] PBMC (Proper scales) (Clone)
[10000] DBMC (Dranas agains) (Clana) 20 files (synar	t otot

- 2-2-6-2 Project の編集・削除
- Experiment manager 画面上部の Projects ボタンをクリックし、 View all projects を選択し ます。
- ② 編集したい Project の Actions 欄の : ボタンをクリックし、閲覧したい場合は view を、編集したい

場合は Edit を、削除したい場合は Delete を選択します。

③ 編集する場合、設定については上述の Project の作成を参照ください。

2-3 Experiment の新規作成・データアップロード

新規の解析を行う場合は Experiment を作成しデータをアップロードします。

- ① Experiment Manager 左上の 🕣 New Experiment をクリックします。
- ② Experiment Name、Purpose を入力して(その他は任意)、Create Experiment ボタン(2項目入力するとアクティブになる)を押します。

Cytobank	Experiments	Projects	Help	🖸 ynakane
? Create N	New Experiment	Create via ACS File		
	* Experiment Name			
	Project	None		
	* Primary Researcher	Yuko Nakane		
	* Principal Investigator	Yuko Nakane		
	,	☑ Invite a new user Allow Principal Investigator to have full access to experiment		
	Source(s)	Nothing selected		
	* Purpose			
	Comments			
		Create Experiment		

※ACS ファイル経由で Experiment をアップロードをしたい場合は、Create New Experiment 画面の Create via ACS File ボタンから行うことができます。

③ Add Files ボタンをクリックしてアップロードファイルを選択するか、アップロードしたいファイルを直接 枠の中へドラッグ&ドロップします。

※データセットのサイズが大きい場合は Zip 圧縮し、枠上部の Upload File Using a Zip file ボタンからもアップロードでき、時間を短縮することができます。

Cyt	obank	Experiments Pro	ojects						Help	ynakane
tes	t 5									Sharing
≡	Upload	f Experiment Files								Private
?	D. A	dd More								
		Name			φ	Size	φ	Туре	\$	
		Sample 1(Control).fcs				34.71 KB		FCS		
		Sample 2.fcs				34.71 KB		FCS		
		Sample 3.fcs				34.71 KB		FCS		
		Sample 4.fcs				34.71 KB		FCS		
		Sample 5.fcs				34.71 KB		FCS		
		Sample 6.fcs				34.71 KB		FCS		
		Sample 7.fcs				34.71 KB		FCS		
		Sample 8.fcs				34.71 KB		FCS		
	You	u can also still drag and drop	files 🎽			0.27 MB selected	1	8 files		
				🕆 Upload files						

④ ファイルを選択後、Upload files をクリックするとアップロードが開始します。

※インターネット通信速度が不安定な環境の場合、速度が低下した時にアップロードが停止してエラーが 表示されたり、アップロードをやり直して重複してしまうことがあります。これを回避したい場合は多数のデー タファイルを一度にアップロードせず、数から十数ファイルずつアップロードすると良いです。 ※ファイルアップロードのやり直し等が発生し、Experiment内にファイルが重複してしまった場合は、 <u>Clone Experiment</u>で、重複したファイルのサイズの小さい方(または不要なファイル)を FCS files to clone の選択から外して Clone を作成します。

- ⑤ アップロードが完了すると Set up new experiment ページが表示されます。
- Set up new experiment では、データの QC・前処理の 4 つのステップ(1.スケール調整、2.コンペンセーション、3.サンプルタグ、4.パネル・チャネル名)が表示されており、各ステップをクリックすると、機能ボタンにポップアップが現れてガイドします。

Cytobank	Experi	ments Projects					🔧 Admin	Help	💽 Cyto	obankSuppor
PBMC Experim Actions	ent (Fluores	cence) ags 🏠 Gates	है Advanced analyses	L. Illustration	s		10 Scales	∆∆ Com	pensation	Sharing
Set up ne	ew experin	ient				Click here to a	djust your exp	eriment sc	ales.	
C							Learn more			
Succ	ess: rou up	oaded 10 files to t	nis experiment. Here are i	the next steps:						
-	1. A	djust your scales ?			(i)					
	<u>A</u> 2. S	et the experiment o	ompensation ?		(i)					
	🌱 3. A	nnotate your files w	ith sample tags ?		i					
	97 4. F	leview or adjust you	r panel assignments or char	nnel names ?	í					
	5. C	one with setup and	ready to start exploring yo	ur experiment da	ta ?					

- ガイドされた機能で調整が完了したら、機能ウィンドウを Close するかメニューバーに表示される
 Set up experiment ボタンで戻ります。
- ⑧ ステップが完了すると緑色のチェックマークが付きます。
- ④ 4ステップが完了したら、5. Done with setup and ready to start exploring your experiment data?をクリックします。

(10) 解析とレポートのボックスが表示されますので、次に行いたい作業をクリックします。

0	Cytobank		Experiments	Projects								🔧 Admin	Help	💽 Cyto	bankSupport
	PBMC Experin	ment (Fl 🌱 Sa	luorescence) ample tags	🗘 Gates	ភ្លំ Adv	anced analyse	25	L Illustration	5			E Scales	🕰 Cor	npensation	Sharing
<	? Set up n	iew exp	periment								Configure uplo	oad landing page	• (j)		
	Plea	ase revi	iew the follow	wing steps to	set up y	our experime	ent:								
		10 ²	1. Adjust y	our scales ?					i	•	✓				
		2. Set the experiment compensation ?			on ?			i	•	~					
		۷	3. Annotati	e your files wi	th sample	tags ?			i	•	*				
		æ	4. Review of	or adjust your	panel ass	ignments or o	channel	l names ?	i	•	✓				
	Set	up coi	mplete! Nov	v you can sta	art explo	ring your ex	perim	ent data:							
			An	alysis			Re	eporting							
	☆ Gating Editor ④				e illustration	í									
	state Run an advanced analysis ①					ort statistics	í								

※各ステップは任意にスキップすることもできます。調整済みのファイル等で不要な場合、5. Done with setup and ready to start exploring your experiment data?をクリックし、すぐに先の作 業へ進むこともできます。

2-3-1 Experiment 画面

Experiment 画面では、上部に Experiment 名と機能ツールが用意されています。

Cytobank	Experiments	Projects					Help	
PBMC (Fluoresc Actions	ence) 🎙 Sample tags	ル- Data QC	🗘 Gates	ដាំ Advanced analyses	L Illustrations	™ ™ Scales	A Compensation	Sharing Private

2-3-2機能ツール一覧

- <u>Actions</u>: <u>Experiment サマリー</u>へのアクセスや<u>複製</u>、各種 <u>Upport/Export</u>機能
- Sample Tags: サンプルタグ機能
- <u>Data QC</u>:測定時の異常イベントを自動検出・除去
- Gates : ゲート作成・確認・編集機能
- <u>Advanced Analysis</u>: 高度な解析ツール(クリックすると以下が選択できます)
 - <u>Dimensionality Reduction</u>:次元削減アルゴリズム(tSNE-CUDA, UMAP, opt-SNE, viSNE)
 - ・ Clustering: クラスタリングアルゴリズム (FlowSOM, SPADE)
 - <u>CITRUS</u>
 - Automatic gating
- Illustrations: <u>Figure の作成や統計グラフ、統計的有意差検定機能</u>
- <u>Scales</u>:スケール設定
- <u>Compensation</u>:コンペンセーション設定
- Sharing Private: <u>Experimentの共有機能</u>

2-4 Actions:

Experiment サマリーの表示や、複製、データの Export/Import などを行うことができます。



2-4-1 View Summary

Experiment の詳細情報、関連する Experiment との相関、各種解析結果、fcs ファイル以外のファ イル、fcs ファイル情報などすべての情報がまとめられています。

- ・ Notes: Experimentの目的やコメントを記載できます(クリックで編集可)。
- Linked experiments: 関連する Experiment との相関図が表示されます。クリックするとその Experiment へ移動できます。下図 Linked experiments 例は Demo test という Experiment のデータセットで FlowSOM、SPADE、viSNE の解析を行い、viSNE 解析結果を 複製(クローン)したときのものです。

्रे	Cytobank Premium	Experiments	Projects		
[Demo test Actions	L Illustrations	🌱 Sample tags	តិ Advanced analyses	🗘 Gating
<	? Experime	nt Summary	Clone		
^					
_		Name: D	emo test		
		Project: N	one		
1		Labels: N	one		
1					
4		PR: MYuko Naka	ne		
1		PI: 🛛 Yuko Naka	ne		
	Uploa	der: 🔃 Yuko Naka	ne		
1	Creat	ted: 11:22 pm, Jan 2	22 2020		
Ø	Upda	ted: 11:05 PM			
0					
	Notes				
5		Purpose: de	emo		
		Comments: N	one		
	@ Linker	evneriments			
	<i>⊘</i> LIIIKet	experiments			
			De	mo test	
			*		1
	e	FlowSOM demo	🌩 SP/	ADE demo	VISNE demo
					↓ SNE demo

 Illustrations:作成中と保存された illustration のリストが表示されています。名前をクリックする とその illustration に移動します。Experiment を共有していて、他のユーザーが illustration を 作成している場合は Author のユーザー名で判断できます。自分が作成した illustration につい ては、右端のゴミ箱アイコンをクリックすると削除でき、Delete all your illustrations ボタンで全て 削除もできます。

Illustrations Onew					
Name	Author	Created	Updated		
L CD19 histogram		May 2017	Jul 16	8	
🔏 Heatmap (Raw values)		May 2017	May 2017		
🛃 Density dot plots		May 2017	Jul 16	8	Ô
😹 Histogram overlay		May 2017	May 2017	8	
				Delete all y	our Illustrations

- ・ Clones: その Experiment の複製が作成された場合、リストで表示されます。
- ・ 各種 Analyses : 解析の結果や実行ステータスがリストとして表示されます。実施したものだけが表示されます。不要な結果は右端のゴミ箱アイコンをクリックすると削除できます。

ፋ Dim	Dimensionality reduction analyses 💿 🔞 New											
	Name	Settings	Size	Author	Status	Created	Updated	Delete				
1	Cytobank's tSNE-CUDA analysis (200i) 🧷	o ^O Settings	24.6 MB		Completed	Jul 19	09:52 AM	â				
14	Cytobank tSNE-CUDA analysis (400i) 🧷	° Settings	24.6 MB		Completed	Jul 19	09:51 AM	â				
14	Cytobank t5NE-CUDA analysis 🥒	° Settings	24.6 MB		Completed	Jul 16	09:48 AM	8				
U	Cytobank UMAP analysis 🧷	₀ ^O Settings	24.6 MB		Completed	Jul 16	Jul 20					
U	Cytobank UMAP analysis (copy) 🖉	° Settings			New	Jul 20	Jul 20					

 Attachments: Attach a File ボタンで fcs 以外のファイルを添付できます。fcs ファイルも添付フ ァイルとしたい場合は、Zip フォルダに格納することで添付できます。Word、Excel、PowerPoint はもちろん、顕微鏡の画像ファイルなど関連するファイルを添付し、Experiment 内にまとめて管理 できます。PNG、JPEG、TIFF ファイルの場合は位置情報(EXIF データ)はファイルのメタデータか ら削除されます。FCS と PDF ファイルのメタデータは削除されません。

6	Attachments						
	File Name	Download / View	Date	Uploaded By	Size	md5sum	
	experiment_12_May-03-2017_03-08-PM.acs Original ACS File that this experiment was created from.	B	Jul 16	the second second	663 MB	60dcb4b	
	Upload an attachment						

- ・ Protocols:実験や測定のプロトコルファイルをファイルアップロード、URL、手入力から選択して添付できます。
- FCS Files: Experiment に含まれている fcs ファイル数とその名前、細胞数(Events)、ファイ ルサイズや Sample tags のリストを確認することができます。また、Panel は各 FCS ファイルのチャ ネルパラメータパネルを表示しており、同一のファイルには同一の番号のパネルが割り当てられていま す。Panel 部分をクリックすると、パネルの内容(チャネル数、名前、ファイル名)を確認・編集でき る <u>Assign panels</u>機能へ移動します。

FCS files — 3					
File Name	Sample Name	Sample Tags	Panel	Events	Size
Marrow1_01_Basal1.fcs @ details		Basal, basal, Plate 1	Panel 1	351,619	55 MB
Marrow1_06_BCR.fcs @ details		BCR, bcr, Plate 1	Panel 1	472,082	73.8 MB
Marrow1_12_IL7.fcs @ details		IL7, II7, Plate 1	Panel 1	507,989	79.5 MB
			Total:	1,331,690	208 MB

※fcs ファイルの詳細について

FCS file summary: Marrow1_01_Basal1.fcs		
 Marrow1_01_Basal1.fcs - FCS File Information 		
FCS File Summary		
Version:FCS3.0Sample Name:Marrow1_01_Basal1.fcsFile Size:57668673md5sum:e499bceecf6a33c16f858b8c74c4627bEvents:[] 351619Channel Count:41Tube Name:File Size:Plate Name:File Size:Well ID:File Size:Cytometer:2021-07-16 20:58:16 UTC	Cell Length	
* FCS File Diagnostics		
We have not detected any problems with this FCS file.		
▼ FCS File Keywords		
Required Keywords (FCS3.0)	Optional Keywords	
\$BEGINANALYSIS 0 \$BTIM Acc	quisition Begin Time:	
\$BEGINDATA 3149 \$COM Con	mment:	

アップロードした fcs ファイルはデータ取り扱いの信頼性の観点から、削除できない仕様になっています。特定のファイルを除いた Experiment にしたい場合は、Clone 機能の FCS files to clone で除きたいファイル以外のファイルのみを含む Experiment を複製することが可能です。

2 - 4 - 2 Clone

Experiment を複製することが出来ます。

Experimentを共有した後に複製することで元の Experiment を上書きすることなく、共有者も再解 析することができるようになります。

また、option 設定によって、元 Experiment の一部の情報を反映したものとしてや、特定のゲートポピ ュレーションのみでの複製を作成することもできます。

① Clone ボタンをクリックします。Create Clone of Experiment ウィンドウが表示されます。

	PBMC Experiment (Fluore	scence) tags 🖒	7 Gates	ភ្លំ Advanced analyses	L Illustrations
	Experiment actions View summary 		Clone	🥒 Edit detail	s
	🗞 Clone	-			
54.	1. Upload	e: PBM	C Experime	nt (Fluorescence)	
6	🛃 Export	t: None	C Experime	nt (Fluorescence)	
8 1	🔗 Link to parent	s: None			
Q	🗊 Reset				
2	💼 Delete				
1.21					

- ② Experiment name:に複製した Experiment 名を入力します。デフォルトでは元の名前+ (Clone)となります。
- ③ Sharing:で共有アクセスを選択できます。
 - Clone individual user access:元の Experiment を共有した共有者に、Clone した Experiment へのフルアクセスも可とします。
 - Clone project membership: Project 内のメンバーのアクセス権限が Clone にも反
 映される設定です。デフォルトではチェックが入っています。
 - Link to originating experiment: チェックを入れると作成した Clone は元の Experiment に紐づけられます。デフォルトでチェックが入っています。

S Create clos	ne of experiment			×
Experiment name Purpose	Sample Experiment		5	
Owners	Principal Investigator Cytobank Support	Primary Researcher Cytobank Support		
Sharing	Allow PI full experiment access Clone individual user access Link clone to this experiment	 ▲ Ø Clone project membership Ø 		

- ④ Clone options: で Clone に含める (反映する) 内容をすべてまたは個別に選択できます。
 - Select all : すべて
 - ・ Clone gates : ゲート
 - ・ Clone compensation : コンペンセーション
 - ・ Clone sample tag : サンプルタグ
 - ・ Clone file panels:マーカーパネルファイル
 - ・ Clone illustrations:保存した illustrations
 - ・ Clone attachments and protocols:添付ファイルとプロトコル
 - Clone reagents: 蛍光色素情報

※gateとcompensationなど、一緒である必要がある項目は、Cytobankが自動で組み 合わせます。

 ⑤ FCS files to clone : Clone に含めるファイルを選択することができます。Select files…ボタン で選択するか、下のリストで Clone に含めるファイルのみを選択します。

S Create Clo	ne of Experiment			
Experiment name	Sample Experiment (Clone)			
Purpose	A purpose goes here			
Owners	Principal Investigator Cytobank Support		Primary Researcher Cytobank Support	
Sharing	Clone sharing access	Ł	Clone project membership	-
Clone options	Se	ect all	(Full clone)	
	Cione gates	$\hat{\Omega}$	Clone compensations	M
Choose what cloning options to include	Clone sample tags	*	Clone file panels	
within the new experiment	Cione illustrations		Clone attachments and protocols	0
	Clone reagents	*		
FCS files to clone	Select files Q. Filter displayed		1	
11 files selected	All displayed			
Choose what files	None displayed			
to clone based on these options	Only experimental files			
	PE - comp_pe.fcs PE_Cy7 - comp_pe_cy7.fcs PacBlu - comp_pacblu.fcs PerCP_Cy55 - comp_percp_cy55.fcs			

 Split Files by population:元 Experiment 内の特定のゲートポピュレーション (populations)を各 FCS ファイルに分割して Clone を作成することができます。Select population ボタンをクリックするとゲートリストがポップアップしますので、チェックを入れて選択するこ とができます。

Experiment name	PBMC Experiment (Fluorescence) (QJ) (Clone)			IE.	Select Populations to split files by	Go to Gating Editor	×		
Purpose	Stimulate PBMC with IL6, IL10, and LPS. Measu	re phosphorylation of Stat3 and p38 after	5 /						
Owners	Principal Investigator Cytobank Support Allow Pf full experiment access Cone individual user access Link to originating experiment P	Clone		All None Inverse All None Inverse I intact cells CD33+ monocytes CD33- lymphocytes	e I ^A Z]			
Clone options	Select all (Full clone)	Δ		CD20+ B cells CD3+ T cells CD3+CD4+ T cells		Ĵ		
	Cione sample tags			CD3+CD4-T cells					
	Cione illustrations	Cione attachments and protocols	@ FCS files	C 0	Displaying all 8 populations to split files by. I of 8 are selected.	С	lose		
FCS files to clone 11 files selected	Select Hirs)	11 file	selecter					
Split files by populations	Select pop	ulations	Spl	it files by	S Select populations				
~ Marx as Clone	Yes	Cancel Su	Mari	k as Clon	e Yes 🗸				

2-4-2-1 不要な FCS ファイルの Experiment からの削除

アップロードした fcs ファイルはデータ取り扱いの信頼性の観点から、削除はできない仕様になっています。 上述の手順⑤で不要なファイルを対象から外した Clone を作成することで、必要なファイルのみの Experiment とすることができます。

2-4-2-2 特定のポピュレーションをファイルに分けて Clone Experiment を作成 上述の手順⑥で FCS ファイルとして分けたいポピュレーションを選択します。 例えば、元 Experiment に 3ファイルがあり、 Split files by populations で popA と popBの2つのポピュレーションを選択する と、 popA の細胞のファイル 3 つと popBの3 つを含む Clone Experiment が作成されます。

2-4-3 upload

PBMC Experiment (Fluorescence) ■ Actions ▼ Sample tags 🗘 Gates R Advanced analyses Experiment actions nt View summary 🗞 Clone eriment: 狩 Upload sample tags ± Export ① Upload gates from gating-ML Upload more files 🔗 Link to parent 🗊 Reset the experiment compensation ? 📅 Delete otate your files with sample tags ? 🔧 Edit reagents [admin only] 4. Review or adjust your panel assignments or char

Experiment に下記 3 種類のファイルをローカルからアップロードができます。

2-4-3-1 Upload Sample Tags : サンプルタグファイルのアップロード <u>エクスポート後編集したサンプルタグファイル</u>をアップロードして Sample tags に適用します。

2-4-3-2 Upload Gates from Gating-ML: <u>Gating-ML ファイル</u>のアップロード ※Cytobank が使用している Gating-ML のフォーマットは the International Society for Advancement of Cytometry (ISAC)によって規格された Gating-ML 2.0 であり、その他の規格 ファイルはエラーとなります。

Spidlen, Josef *et al.* "ISAC's Gating-ML 2.0 data exchange standard for gating description." Cytometry. Part A : the journal of the International Society for Analytical Cytology vol. 87,7 (2015): 683-7. doi:10.1002/cyto.a.22690

2-4-3-3 Upload More Files: FCS ファイルの追加アップロード

2-4-4 Export

Experimentの様々な情報を Export できます。



2-4-4-1 Export Statistics:統計計算結果の出力

様々な統計値やサンプルタグ、ファイルキーワード情報を任意に選択し、全ファイルの一覧表を csv ファイ ルとしてエクスポートします。

- ① Actions> Export> Export Statistics をクリックします。
- ② 新規に作成する場合は Define or new を選択します。テンプレートがあり、呼び出す場合は Load a saved template リストから選択します。



- ③ (必要な場合)Export Optionsを開きファイルの設定をします。
 - Filename: ファイル名
 - Orientation one FCS file per: FCS ファイルを項目として行(Column)と列 (Row)のどちらに配置するかを選択します。
 - Compensation: どのコンペンセーションを適用するかを選択します。

- FCS Files to include:出力対象のファイルを選択します。Experiment にインポートされている File type が設定されているファイルから、出力するファイルの種類を以下のパターンで選択できます。
 - サンプルファイルのみ (Only experimental files)
 - ・ サンプルとコントロールファイル (Experimental and other control files)
 - サンプルとコンペンセーション単染色ファイル (Experimental and compensation files)
 - ・ サンプルとコントロールとコンペンセーションファイル
- Delimiter type : 出力ファイル形式の選択
- Append time stamp to filename : ファイル名にタイムスタンプを入れます。
- Include descriptive file header : ファイルのヘッダーを含めます。
- Include Gating ML file with archive: アーカイブと Gating ML ファイルを含めます。
- ④ Define what to export で、統計計算、サンプルタグ、ファイルキーワードから選択します。
- ⑤ 統計計算を選択するとメニューが現れ、以下の順番に対象を指定していきます。
 - Pick a statistic type:使用する統計
 - ・ Median:中央値
 - ・ Mean:平均値
 - Coefficient of variation:変動係数
 - Percent of population : population A /population B ※BはAの上流ポピュ レーションという前提。計算は厳密には (A⊆B)/B で算出される。
 - Standard deviation:標準偏差
 - Variance:分散
 - Minimum:最小值
 - ・ Maximum:最大値
 - Event count:細胞数
 - Event count ratio: Population A 細胞数 /population B 細胞数※ポピュレーション階層考慮なし。
 - ・ Channel range: チャネルの範囲
 - Geometric mean:幾何平均
 - ・ Percent in gate:指定したゲート内イベントが何%含まれるか
 - Pick a population : 対象の細胞集団
 - Pick a channel : 使用するチャネル(複数選択可)※使用する統計によっては非表示
 - Enter column header name, separated by comma:指定された対象がコンマ区切りで表示されます。

 6 Add column ボタンを押すと Column defined の表に追加できます。表では、Column name をクリックすると名前を変更できます。また、左の×ボタンで表から削除できます。

PBMC E	Experiment	t (Fluorescence Illustrations	e) 🌱 Sample Tags	🔶 SPADE	🗳 visne	() CITRUS	🟠 Gating		Scales 🕅 Comps
К Е	xport ex	periment s	tatistics	Settings Ter	mplates			Export statistics file	preadsheet here
, D	Export C Define what	options	Click to exp	pand configurat	tion options	cted:	Pick population Y:	Enter column header names, separated by	
	Add	id Statistic Sample Tag	Medians Means Coefficients of Percent of Pop Standard Devi Variances	Variation outation (1) ations	Ungated intact cells CD33+ monocyl CD33- lymphoc CD3+ T cells CD20+ B cells	tes ytes	Ungated Intact cells CD33+ monocytes CD33- lymphocytes CD3+ T cells CD20+ B cells	 commas: % of CD3+ monocytes in intact cells, % of CD20+ B cells in intact cells, % of CD3+CD4+ T cells in intact cells, % of CD3+CD4+ T cells in intact cells 	*
Template bi different ph to export c	uilder: click hases to cre Columns d	through the ate a statistic efined	Minimums Maximums Event Counts Event Count R	ratios (i) 🔻	CD3+CD4- T ce CD3+CD4+ T ce DSTAT3+ Monor All	ells cvtes	CD3+CD4+ T cells CD3+CD4+ T cells DSTAT3+ Monocytes All None	Add columns	
			Statistic definitio	ons appear as ro	ws in this table			Click to add the statistics for export	A Delete all rows
	#		Column name	Y	Category	Туре	Populatio	on Channel	Delete statistics
	1	Ever	it Counts of Ungated		Statistic	Event Count	ungated	1	· •
	2	Event	Counts of intact cel	Is	Statistic	Event Count	intact cell	ls	0
	3	Event Co	unts of CD33+ mono	cytes	Statistic	Event Count	cD33+ mono	cytes	8

- ⑦ Export statistics file ボタンでエクスポートを開始します。
- 2-4-4-2 Export events:細胞情報を FCS または Text ファイルでエクスポート
 - ① Actions> Export> Export events をクリックします。
 - ② Export events ウィンドウの Settings で以下について設定します。
 - Filename: エクスポートするファイルの名前を入力します。
 - Compensation:適用するコンペンセーションを選択します。
 - Transformation:スケール設定の情報を反映するかを選択します。
 - File format: ファイル形式を FCS か Text か選択します。
 - 3 エクスポートの対象とする Populations と FCS Files、(Text 形式で出力する場合は Channels)を Choose ボタンを押して、選択します。
 - ④ Submit export ボタンを押します。

-					
Filename experiment_o	demo_exported_ev	vents			
Compensation File-Internal Com	npensation -				
Transformation Untransform Transform	med ed based on scale	esettings			
File format OFCS • Text					
Include header with FCS file	ename				
Populations	FCS	Files		Channels	
Populations	Choose FCS	Files	Choose	Channels	Choose
Populations Ungated	Choose IL10	Files	Choose	Channels FSC-A	Choose
Populations Ungated Unselected populations:	Choose 1L10 1L6	Files	Choose	Channels FSC-A FSC-W	Choose
Populations Ungated Unselected populations: CD3-	Choose IL10 IL6 LPS	Files	Choose	Channels FSC-A FSC-W SSC-A	Choose
Populations Ungated Unselected populations: CD3+ CD3+	Choose IL10 IL6 LPS Uns	Files) tim1	Choose	Channels FSC-A FSC-W SSC-A pStat3	Choose
Populations Ungated Unselected populations: CD3+ CD3+CD4+ CD3+CD4+ CD3+CD4+	Choose IL10 IL6 LPS Uns Uns	Files	Choose	Channels FSC-A FSC-W SSC-A pStat3 CD33	Choos
Populations Ungated Unselected populations: CD3- CD3-CD4- CD3-CD4- CD4- CD4- CD4-	Choose A IL10 IL6 LPS Uns Uns	Files	Choose	Channels FSC-A FSC-W SSC-A pStat3 CD33 DS-TR A	Choose

- ⑤ ステータスバーが表示されます。完了するとメール通知がくるので、Cytobankから離れても大 丈夫です。
- ・一度エクスポートしたファイルは Experiment summary ページの Exported files セクションに表示され、いつでもダウンロードできます。

2-4-4-3 Export Experiment to ACS: ACS 形式ファイルのエクスポート

Experiment に含まれるデータや illustrations、Attachment ファイルなどすべてをまとめてアーカイブ 化し、ACS ファイルとしてエクスポートします。エクスポートした ACS ファイルはダウンロードすることで、ローカ ルへ保存することができます。ファイルは Cytobank 独自フォーマットのため、他のソフトウエアでは機能し ませんが、サーバーの異なる Cytobank へはアップロードすることで移行できます。また、90 日以上 Cytobank のライセンスを停止しサーバーから Experiment が削除された後に、ローカルに保存しておい た ACS ファイルをアップロードすることで復元することも可能です。

- ① Actions> Export> Export ACS をクリックします。
- ② その Experiment に紐づいている Experiment によっては下記の3つから対象を選択します。
 - Express export:その Experiment だけを ACS としてエクスポートします。

🚍 Actions - 🛛 🦅 S	ample tags	Gates	Advanced analyses
Experiment actions Oview summary	ľ	Clone	🖉 Edit details
⊗ Clone		e: 1-SPADE demo	
🛓 Export	•	Export statistics	3
Link to parent		Export events	
S Reset		😨 Export experime	ent to ACS
Delete		Express export	~5.3 MB
		Export with advance	ed analyses ~165.2 MB
NOCES	Source(Export with linked	descendants ~208.4 MB

- Export with advanced analysis : その Experiment + Linked Experiment で紐づ いている Advanced analysis をまとめて ACS としてエクスポートします。
- Export with linked descendants: その Experiment + Linked Experiment で紐づ いているすべての下流 Experiment をまとめて ACS としてエクスポートします。

※情報量が多く出力に時間がかかるため、ファイル出力終了後にメールにて連絡が入ります。

2-4-4-4 Export Sample Tags: サンプルタグ情報をファイルとしてエクスポート Experiment の各 FCS ファイルにアノテーションした <u>Sample tags</u>の情報を <u>csv ファイルとしてエクスポ</u> <u>ートして、Excel で一覧表表示し、編集や tags を追加</u>することができます。編集後は表を.txt、.csv ま たは.tsv で保存し、<u>Upload sample tags</u>でアップロードして Experiment に適用することができま す。 2-4-4-5 Export Gating-ML: ゲート情報を Gating-ML ファイルとしてエクスポート Cytobank では、International Society for Advancement of Cytometry(ISAC)によって 開発された <u>Gating-ML 2.0</u>の規格を採用しています。Gating Editor で作成したゲート情報(最後 に Apply gates ボタンを押して保存した <u>Experiment gate</u>のバージョン)をファイルとしてエクスポート し、<u>ダウンロードして他の Experiment にインポート</u>したり、ローカルで保存や編集したりもできます。

2 - 4 - 4 - 6 Download Files : fcs ファイルや Attachment ファイルをダウンロード Experiment の FCS ファイルや Attachment ファイルを一括ダウンロードします。

>	Filename \$	÷	Description	÷	Size	÷	Туре 🔶	Events
* o B	pbmc_lrs005_il10.fcs		IL10		22.3 MB		FCS	32537
٥B	pbmc_lrs005_il6.fcs		IL6		22.4 MB		FCS	32683
C ti	pbmc_lrs005_lps.fcs		LPS		22.3 MB		FCS	3254
٥B	pbmc_lrs005_unstim1.fcs		Unstim1		22.4 MB		FCS	3266
o B	pbmc_lrs005_unstim2.fcs		Unstim2		22.3 MB		FCS	3246
οB	comp_alexa 488.fcs		Alexa 488		494 KB		FCS	6981
o B	comp_alexa 647.fcs		Alexa 647		472 KB		FCS	6663
οB	comp_pe.fcs		PE		455 KB		FCS	6428
	comp_pe_cy7.fcs		PE_Cy7		508 KB		FCS	7182
٥B	comp_pacblu.fcs		PacBlu		498 KB		FCS	7041
08	comp_percp_cy55.fcs		PerCP_Cy55		433 KB		FCS	6114

- ① ダウンロードしたいファイルにチェックボックスを入れます。一番上のボックスですべてのファイルを一括オン/オフできます。
- ② Filename 隣のダウンロードアイコンをクリックすると、直接ダウンロードが始まります(ブラウザ経由で OSの設定するフォルダへ保存)。
- ③ Zip&Download Files ボタンをクリックするとチェックボックスオンのファイルが一括ダウンロードの処理 に入ります。ダウンロードリンクの記載されたメールが登録メールアドレスに届きます。

※合計サイズが大きすぎる場合、エラーとなりメールも届きません。1日経過してもメールが届かない場合 はファイル数を減らすか、Export events にて FCS フォーマットでのエクスポートをご検討ください。 ※メールのリンクには有効期間(半日程度)があり、期限を過ぎるとエラー画面が表示されます。

2-4-5 Link to parent : Experiment のリンク付けと解除

任意の Experiment を関連する Experiment (親) として紐づけ (リンク) したり、解除したりするこ とができます。リンクされると <u>Experiment Summary</u> 画面の Linked experiments の相関図に親 Experiment として表示されます。

※下流の Experiment を残し、親または上流の Experiment を削除したい場合は、この機能で先に リンクを解除してから、削除を実行してください。

- 2-4-5-1 任意の Experiment を親としてリンク付けする
- ① リンクを解除したい Experiment の子(下流側) Experiment を開きます。
- ② Actions > Link to parent をクリックします。
- ③ Select an experiment to link as parent で親 Experiment を指定します。
 ※フィルターボックスに名前や Experiment ID 入力すると便利です。



- ④ Apply changes を押します。
- 2-4-5-2親 Experiment からリンクを解除
- ① リンクを解除したい Experiment の子(下流側) Experiment を開きます。
- ② Actions > Link to parent をクリックします。

P	Edit linked parent experiment	×
Curr	ently linked parent: [156745] - 1-SPADE demo	
Š	Unlink parent	
Seleo	t an experiment to link as parent:	
Sele	ect a parent experiment	•
	Cancel Apply changes	

③ OK を押します。

2-4-6 Reset: Experiment を初期化

2-4-7 Delete: Experiment を削除

Experimentは Trash(ゴミ箱)へ移動します。

2-5スケール調整

メニューバーの Scales で、スケール調整を行うことができます。

PBMC Experim	nent (Fluorescence)							Sharing
Actions	🌱 Sample tags	J- Data QC	🗘 Gates	n Advanced analyses	L Illustrations	Scales	A Compensation	Private
? Experime	ent summary	Clone	∥ Ec	dit details				

※Advanced analysis では、スケール調整が大きな影響を及ぼしますので、実行前に確認し、必要な場合調整を行ってください。

2-5-1スケールの種類

以下の3種類が選択できます。

- Linear
- Log
- Arcsinh: Biexponential, logicle, hyperlog 等と同じく0 近傍がリニア、それ以降が Log のハイブリッドなスケールの表示形式で、パラメータ Argument は0 近傍のリニア幅を調 整するパラメータです。

lick	to edit scale settings for a channel	(1) 🖸	R Filter o	displayed	Clear				
	Channel	Туре	Arg.	Minimum	Maximum		File/sample	name	(4
	Ax488-A, pStat3	Arcsinh	150	-200.0	262144.0	(2)	IL10 - pb	mc_lrs005_il10	•
	Ax647-A, pp38	Arcsinh	150	-200.0	262144.0		_	40 4 4	
	Ax700-A	Arcsinh	150	-200.0	262144.0		-	4	
	Ax750-A	Arcsinh	150	-200.0	262144.0	,			
	FSC-A	Linear	1	1.0	262144.0				
	FSC-W	Linear	1	1.0	262144.0				
	PacBlu-A, CD4	Arcsinh	150	-300.0	262144.0		:		
	PacOrange-A	Arcsinh	150	-200.0	262144.0				
	PE-A, CD33	Arcsinh	150	-500.0	262144.0				
	PE-Cy7-A, CD3	Arcsinh	150	-400.0	262144.0		FSC-A	and the time time tare t	-
	PE-TR-A	Arcsinh	150	-200.0	262144.0		Population	Ungated	
	PerCP-Cy55-A, CD20	Arcsinh	150	-300.0	262144.0		Compensation Plot type	comp-pok	
	Qdot525-A	Arcsinh	150	-200.0	262144.0		. iotijpe	at managram	
	Qdot605-A	Arcsinh	150	-200.0	262144.0				
	Qdot655-A	Arcsinh	150	-200.0	262144.0				
	Qdot705-A	Arcsinh	150	-200.0	262144.0				
	SSC-A	Linear	1	1.0	262144.0				
	Time	Linear	1	1.0	262144.0	-			
Scal	tolit scales for all selected and unfiltered eType Argument Minimu port scales from another experiment scl an experiment	(5)	Maxin	num	Apply				

2-5-2スケール調整画面: Scale Editor

- ① 入力する文字列でリストをフィルター表示することができます。
- Type、Arg.、Minimum、Maximumの各項目をクリックすると、選択または数値入力で変更することができます。
- ③ まとめて変更したい場合は、表左の□にチェックを入れ、Bulk edit scales for all selected and unfiltered channels で変更したい項目を変更し、Apply ボタンを押して適用します。
- ④ Channel 名にオーバーカーソルで表示される目玉マークをクリックすると、右側にそのチャネルのプロットが表示されます。
- ⑤ 他の Experiment のスケール設定を適用したい場合に、元となる Experiment を選択します。
- ⑥ Reset All Scales ボタンを押すとリセットできます。

2-6コンペンセーション(蛍光補正)

コンペンセーションはメニューバーの Compensation ボタンで設定ができます。

PBMC Experim	ent (Fluorescence)				Sh	aring
Actions	🌱 Sample tags	🗘 Gates	ភ្លំ Advanced analyses	L Illustrations	Scales 🖾 Compensation	Private
? Experime	nt summary	Clone			Go to Compensations	

※Cytobank ではコンペンセーションは機器ソフトウエアや Kaluza などで補正を済ませた FCS ファイルを インポートし、File-internal compensation でお使いになることを基本的には強くお奨めいたします。 ※コンペンセーションの再調整が必要なデータセットの場合は、解析を始める前にコンペンセーションを実 施ください。Advanced analysis では用いるコンペンセーションを変更できなくなっております。

2-6-1 Compensation Editor

Experiment で使用するコンペンセーションマトリックスを Experiment compensation:のプルダウンメ ニューで選択できます。FCS ファイルに記載されているコンペンセーションマトリックスを用いる場合は Fileinternal compensation を選択します。



2-6-2新規にマトリックスを作成する

① New matrix ボタンをクリックします。



- コンペンセーションに用いるパラメータをすべて選択し、Create Compensation ボタンを クリックします。
- ③ Name 部分にマトリックス名を入力し、ボックス右の Rename ボタンを押します。
- ④ マトリックスに値を入力します。
- ⑤ 作成したマトリックスを Experiment に適用するには、Experiment compensation:
 のプルダウンメニューで選択します。

2-6-3マトリックスをインポートする

Experiment 内のある FCS ファイルに記載されているマトリックスや他の Experiment で使用されてい るマトリックス、他のソフトウエアで作成した CSV 形式のマトリックスファイルをインポートすることができます。 同一のマトリックスが全ファイルに適用されます。

- ① import ボタンをクリックします。
- ② Import new compensation(s)ウィンドウが表示されます。
- ③ インポート元を選択します。
 - File-internal compensation: どのファイルに記載されているマトリックスをインポートする かを選択します。
 - Experiment: アクセス可能な Experiment がリスト表示されるので、Experiment 名
 や ID から選択します。
 - CSV File: Choose file for upload ボタンから CSV ファイルを選択します。



④ Import compensation ボタンを押してインポートします。

2-6-4 自動コンペンセーション

① Automatic ボタンをクリックします。

Actions	🌱 Sample tags	🗘 Gates	តំ Ad	vanced analyse	s 🛛 🚹 Illustrati	ons
^① <u>∕</u> Cor	mpensation Editor	(i)	Experime	nt compensatio	n: File-Internal Com	pensation 👻
C	reate new compensation	n: 💮 Ner	w matrix	웹 Import	° Automatic	

- Automatic compensation ウィンドウが表示されますので、コンペンセーションに用いるパラメータ を選択します。
- ③ 各パラメータチャネルのポジティブとネガティブコントロールのファイルを与えます。Cytobank で自動振り分けをしますが、正しく割り当てられているかご確認ください。
- ④ Calculate Compensation ボタンを押します。
- ⑤ 算出された値の入ったマトリックスとその下に各コントロールファイルのポジティブゲートとネガティブゲートが表示されます。



⑥ マトリックスの値は、セルをクリックして値を入力し、微調整することができます。その下の各コントロール ファイルのゲートは変更できません。

2-7データQC:PeacoQC

2-7-1 自動データ QC ツール: PeacoQC とは

Peak Extraction And Cleaning Oriented Quality Control (PeacoQC) アルゴリズムはフロ ー、マス、スペクトルデータ(スケール調整、コンペンセーション/アンミキシング済)でワークし、測定時のク ロッグ、一時的なインテンシティや流速の変化といった望ましくない状態のときのシグナルの変化をデータか ら除外することができます (Emmaneel et al., 2022)。このアルゴリズムでは、isolation tree (IT) と mean absolute deviation distance(MAD)よって除くべき異常値のイベント検出し、除去しま す。

シグナル強度の変化は誤って新規のポピュレーションに見えてしまうなど、データ解析者をしばしば翻弄する 原因となってしまっています。そのような異常なイベントを除いてから解析を始めることは、結果の質と信頼 性を高めます。Cytobankの PeacoQC: QC プロセスなら、自動的に異常イベントを除去した "Clean"ポピュレーションを作成してくれます。



FIGURE 1 Schematic overview of the proposed method. The raw fcs files should first be preprocessed according to the suggested workflow in the blue box, which can include the removal of margins, compensation, and transformation. The method itself starts with a preprocessed fcs file. First, the sample is binned and density peaks are determined, clustered, and filtered per bin. The cleaning part of the method starts with the usage of an isolation tree, then it will remove peaks based on their MAD distance and it ends by connecting disjointed removed regions. The method returns a cleaned fcs file [Color figure can be viewed at wileyonline[ibrary.com]

Emmaneel A, Quintelier K, Sichien D, Rybakowska P, Marañón C, Alarcón-Riquelme ME, Van Isterdael G, Van Gassen S, Saeys Y. PeacoQC: Peak-based selection of high quality cytometry data. Cytometry A. 2022 Apr;101(4):325-338. doi: 10.1002/cyto.a.24501. Epub 2021 Oct 3. PMID: 34549881; PMCID: PMC9293479.

<u>こちらのガイドビデオ</u>もご参照ください。

2-7-2 PeacoQC 設定·実行

① メニューバー> Data QC をクリックし、PeacoQC 名を入力します。



② FCS files ボックスの Choose ボタンをクリックし、ファイルを選択します。選択できるファイルは Time パラメータを有していて、かつ 500 MB 未満です。500MB 以上のファイルは Time スライス(横軸 Time でゲートを作成し、Clone 機能の Split files by populations で分割)が必要となりま す。800 万細胞以上 1500 細胞未満の細胞イベントのファイルは実行できません。

FCS files	Select FCS files	×
DSS		
LL	Q All None Inverse 1/2	
YSL		
	✓ DSS - DSS_QC	
	✓ LL - LL_QC	
	VSL - YSL_QC	

③ Channels ボックスの Choose ボタンをクリックします。 Time チャネルは自動的に選択されていま す。 少なくとも 2 チャネルを実行対象に選択します。

Channels	- Salact Channels	×
Choose		
Time		
Y585-PE-A	Q All None Inverse 1 ^A _Z	
Y610-mCHERRY-A		
Y710-PC5.5-A	□ Y675-PC5-A [FL6-A]	
Y763-PC7-A	□ Y710-PC5.5-H [FL7-H]	
R660-APC-A	✓ Y710-PC5.5-A [FL7-A]	
R712-APCA700-A	Y763-PC7-H [FL8-H]	
R763-APCA750-A	✓ Y763-PC7-A [FL8-A]	
Unselected channels:	R660-APC-H [FL9-H]	
FSC-A	R660-APC-A [FL9-A]	
FSC-H	R712-APCA700-H [FL10-H]	
SSC-A	R712-APCA700-A [FI 10-A]	
SSC-H	R763-APCA750-H [FI 11-H]	
8 of 49 selected	✓ R763-APCA750-A [FL11-A]	

④ 必要な場合 Advanced settings を設定します。

Remove Margins	
ON	
Use experiment scales	
O Use file-internal scales	
OFF	
Anomaly detection method 🚯	
Both (recommended)	
O MAD (Mean absolute deviation) distance	
O IT (isolation tree) position	
MAD parameter 🚯	
6	
IT limit parameter 🚯	
0.6	
Max bins 🗿	
500	
Consecutive bins 🕕	
6	

- Remove margins: ON にするとダイナミックレンジ外のすべてのイベントを除去します。サイト メトリーデータでは自動で ON に設定されます。スペクトルサイトメトリーデータではデフォルトでは OFF になっていて、必要な場合 ON に変更することもできます。マスサイトメトリーデータの場合 は自動的に OFF で、ON は選択できないようになっています。
- Anomaly detection method:異常イベントの検出と除去に用いる方法の選択を行います。デフォルトでは2つの方法の両方(Both)に設定されています。
 1つ目の方法である IT(isolation tree)は Time パラメータ上の各ビンについてイベントが、密度ピーク内にあるか外れているかで判断します。この方法は微妙な変化を検出できないことがあります。

2つ目の方法である MAD (mean absolute deviation) は、密度ピーク範囲内イベント の中央値を用いて算出された平均絶対偏差に対して設定値の倍数より大きいイベントをすべ て除去します。

- MAD parameter:デフォルトの設定値は6です。下げると基準を厳しくすることができ、上げると緩めて除去されるイベント数を減らすことができます。
- IT Limit parameter : isolation forest tree の構築時に、標準偏差についてどの程度 を外れ値とするかに影響します。デフォルトは 0.6 で値を下げるとより厳密になります。
- MAX bins: PeacoQC プロセスでは Time パラメータについて bins を作成します。各 bin について評価をして、上記の検出方法で異常があるとされた bin を丸ごと除去します。範囲は 40 から 1,000,000 で、デフォルトは 500 です。ほとんどのケースではデフォルト設定で良好で あるため変更することはおすすめしません。
- Consecutive bins:異常 bin と異常 bin の間の bin が設定値以下だった場合、間の bin も削除します。これはノイズ領域の bin が他の安定領域の bin と類似する場合にアルゴリ ズム厳密性が不十分になるためで、値を上げるとより多くのビンを除去します。デフォルトは5 で す。
- ⑤ Transformations でスケールとコンペンセーションが適切に適用されているかを確認します。Data scaling で File internal が設定されている場合は、各ファイルの \$ PnR を読み込んで最大値と 最小値を使用します。Compensation では Experiment compensation に設定されているコ ンペンセーションが使用されます。



2-7-3 PeacoQC 結果の評価

- 2-7-3-1 PeacoQC 結果画面ページでの評価
- PeacoQC が完了すると自動的に結果画面が表示されます。通知メールのリンクボタン、または Actions>View summary の Data QC セクションから、PeacoQC 結果をクリックします。

	🚍 Actions 💦 🌱 Sample ta	ags -M- Data QC		Actions	🌱 Sample tags	-∿- Data QC
<	Experiment actions View summary 		<	? Experim	nent summary	Clone
	🚳 Clone	Your Trial plan	^	-∿- Data	QC	• New
	土 Upload	2022 PeacoOC proc		Name	2	
	🛃 Export		5	11252	2022 PeacoQC proc	ess 🧷
	 Link to parent Reset Delete 	1 T	-/ 3 () 1	Jump to Data 11222 11222	QC 2022 PeacoQC proc 2022 PeacoQC proc	ess (copy) 🥒
		1	9		-	

 PeacoQC 結果画面には、PeacoQC report と Download run info and plots ボタンがあり ます。

Dor	acoOC cotup		Comusettio					P. Developed and
Pea	acouc setup		Copy setting	s			Analysis complete!	Download run
Pro	ocess Name: Cytob	oank's PeacoQC pro	ocess 🖉					
	PeacoQC rui	n information						
			Pe	acoQC rej	port			
		0.2	1.4	0.0	1.4	0.0	Tube1.fcs	
		0.1	2.8	NA	2.8	0.0	Tube2.fcs	
		0.2	0.2	NA	0.2	0.0	Tube3.fcs	
		0.2	2.1	0.0	2.1	0.0	Tube4.fcs	
		0.2	0.0	NA	0.0	0.0	Tube5.fcs	
		0.2	3.3	NA	3.3	0.0	Tube6.fcs	
		0.1	4.0	0.0	4.0	0.0	Tube7.fcs	
		0.2	0.0	0.0	0.0	0.0	Tube8.fcs	
	Consecutive bins IT limit MAD	Events per bin Remove margins	% Full analysis	% IT analysis	% MAD analysis	Consecutive cells	Incr/Decr	
	c	Consecutive b	insIT limit M	AD Event	s per bin	Removed p	ercentage	
		5	0.6	6 0 20	0 400 60	0 0 50 100		
		ncr/Decr Increasing cl No increasing	nannel g or decreasir	ng effect				

- ③ PeacoQC report には各 FCS ファイルについて、以下の情報が数値で表示されているとともに、結果を把握しやすいようにカラーヒートマップになっています。表示される項目は Advanced settings で選択した方法や値によります。
 - % Remove margins: margins によって除去された細胞イベントの%

- % Full analysis: すべての方法(IT・MAD の合計 + consecutive cells)によって除去 された細胞イベントの%
- % IT analysis: IT によって除去された細胞イベントの%
- % MAD analysis: MAD によって除去された細胞イベントの%
- % Consecutive cells : Consecutive cells によって除去された細胞イベントの%
- Incre/Decre:シグナル強度の上昇/下降傾向の有無

※Time に対するシグナルの単調な上昇/下降傾向を PeacoQC は検出し表示しますが、異常値 としては扱いません。ファイルに含まれる細胞数が少ない場合、この機能が過度に働く場合がありま すのでご注意ください。

- ④ より詳細な情報を確認したい場合は、Download run info and plots ボタンをクリックして、Zip フォルダをダウンロードします。
- ⑤ フォルダを開くと以下のようなファイル類が保存されています。
 - 📕 plots
 - 202-217-Cytobank's-PeacoQC-process-heatmap.pdf
 - 202-217-Cytobank's-PeacoQC-process-log.txt
 - 202-217-Cytobank's-PeacoQC-process-report.tsv
- ⑥ このフォルダには上述の PeacoQC report の PDF ファイルとその数値情報の tsv ファイル、run information と Log のファイルと plots フォルダが保存されています。 Plots フォルダには PeacoQC にかけた FCS ファイルごとの png ファイルが保存されており、設定で選択したすべてのチャネルについて X 軸 Time の密度ピークが含まれています。

密度ピークの背景が白の領域は異常なしで、ピンク色の領域は IT、紫色は MAD、薄ピンク色は consecutive cells によって異常が検出され、除去された領域となります。



- 2-7-3-2 PeacoQC 結果が作成した FCS ファイルを用いた Cytobank での評価
- ① PeacoQC が完了すると、「オリジナルの FCS ファイル名_QC.fcs」という名前のファイルが自動的に 作成されます。このファイルには anomaly という名前のパラメータが追加されており、オリジナルの

		Barton	and and files		
Assign	paneis	Configure o	control files		
Panel 1 (Or	iginal pre-QC files)	Uisible	Panel 1 (Cle	aned files) 🥖	
			20 channels	5	Q
19 channel	S	Q	FL4-A	HLA-DR-PC5.5 B488	8-690BP-A
Name	Reagent		FL5-A	PD-1-PC7 B488-780	BP-A
FSC-A	FSC-A		FL6-A	CCR7-AF647 R638-0	560BP-A
FSC-H	FSC-H		FL7-A	TNFa-AF700 R638-7	12BP-A
SSC-A	SSC-A		FL8-A	CD4-APC-A750 R63	8-780BP-A
SSC-H	SSC-H		FL9-A	CD45-PB V405-450	BP-A
TIME	TIME		FL10-A	CD8-KO V405-525B	P-A
FL1-A	CD3-FITC		FL11-A	CD16-BV-605 V405	-610BP-A
FL2-A	IFNg-PE		FL12-A	CD14-BV-650 V405-	-660BP-A
FL3-A	CD137-ECD		FL13-A	CD45RA-BV-785 V4	05-780BP-A
FL4-A	HLA-DR-PC5.	5	FSC-Width	FSC-Width	
FL5-A	PD-1-PC7		Anomaly	Anomaly	
FL6-A	CCR7-AF647				
C O files			□ 8 files		٩
o mes		Q	_ Tube1 - T	ube1_QC	
□ Tube1			🗆 Tube2 - T	ube2_QC	
Tube2			Tube3 - T	ube3 OC	

FCS ファイルとは異なるパネルに割り当てられています。メニューバーの Sample tags> <u>Assign</u> panels で確認することができます。オリジナルの FCS ファイルは Original pre-QC files Panel に、作成された FCS ファイルは Cleaned files panel に割り当てられています。 Visible にチェック を入れると Experiment で扱うことができるようになります。 デフォルトでは Cleaned files panel だけが Visible に設定されています。

② Gating Editor を開くと Gates に Anomaly と Clean のゲートとポピュレーションが自動作成されているのを確認できます。Clean ゲートをクリックすると、PeacoQC が自動作成した Anomaly パラメータが X 軸のプロットが表示され、%を確認することができます。異常イベントが除去された Cleanポピュレーションから、マニュアルゲーティングや機械学習解析を開始すると良いです。



- メニューバー> <u>illustrations</u>> all saved を開くと、2つの illustration が自動作成されているの を確認できます。
 - PeacoQC Quality Check Event count ration heatmap を開くと、Clean 前 (Ungated) に対して後(Clean)の細胞数比(Fold of event count) がヒートマップ で表示されます。PeacoQC によって細胞数がどのくらい減ったかを把握できます。



 PeacoQC Quality Check-Cleaned and Anomaly events comparison は、Time パラメータが X 軸、最初のチャネルパラメータが Y 軸のドットプロットが表示されます。列に FCS ファイルが、行に Ungated、Anomaly、Clean ポピュレーションが並び、異常イベントがどのく らい、どのように発生しているかを把握できます。



2-7-4実行上限の目安

他の解析アルゴリズムと同様、サーバーの計算処理リソースを超えるデータセットの場合、PeacoQC が完 了せず Failed となります。 データセットの合計のサイズ上限は明らかとなっていませんが、以下のようなファイルが1つでも含まれる場合は実行できないようになっています。

- 500 MB 以上 /ファイル
- 500 チャネル以上 /ファイル
- 800 万細胞以上 /ファイル
- 1500 細胞未満 /ファイル
- 細胞イベント数とチャネル数の組み合わせが下表を超えている場合

Range of the number of events (Per file) *M=million	Max number of channels allowed (Per file)		
1.5k-0.2M	500		
0.2M-0.4M	400		
0.4M-0.6M	300		
0.6M-0.8M	200		
0.8M-1M	100		
1m-1.5M	80		
1.5M-2M	55		
2M-3M	43		
3M-4M	35		
4M-5M	29		
5M-6M	23		
6M-7M	17		
7M-8M	10		

PeacoQC が Failed になるのには、チャネル数とイベント数(1ファイルあたりの)が大きく影響します。 細胞数とチャネル数の下図に示すとおり、緑色は緑色が問題なく完了できる範囲、黄色が Failed となる 可能性のある範囲、赤色がメモリオーバーで Failed になる範囲です。



黄色の領域の設定がされている場合は、メモリ上限のため Failed になる可能性があるという Warning メッセージが表示されます。この時点でイベント数や対象ポピュレーションを削減することも可能です。 赤色の領域の設定がされている場合、設定したチャネル数での上限細胞イベント数を伝えるメッセージが 表示されます。このアドバイスに従って選択対象を減らすことで、実行可能となります。

2-8ゲーティング

2-8-1 Cytobank でのゲーティングとポピュレーションの定義

Cytobank ではゲートを作成すると同時に同名のポピュレーションを自動作成し、作成されたポピュレーションはポピュレーションとして扱います。

ゲート内容(位置やサイズの変更など)とポピュレーションは連動しますが、ゲート自体を削除しても<u>ポピ ュレーションは削除</u>されなかったり、ゲート名を<u>変更してもポピュレーション名</u>はそのままだったりします。 ゲートとポピュレーションを分けて扱えるようにすることで、あるゲーティングを後から上流に挿入したり、ある ゲートの NOT や OR を各々別名のポピュレーションとして設定したりすることができます。

メニューバーの Gating ボタンをクリックすると、Gating Editor と Gating hierarchy が選択できます。

■ Actions	🗘 Gates	ាំង Advanced an	
? Experiment summary	🏠 Gating Editor		
	🗧 🔄 Gating 🛛	nierarchy	

2-8-2サーバー保存のタイミングと Experiment gate 情報としての保存

Cytobank では Gating Editor の変更を常にサーバーで保存しています。これに加えて Apply gates ボタンを押した時の情報を Experiment gate 情報としてバージョン保存も行います。Gating Editor 以外の機能では、保存された最新の Experiment gate 情報のみを使用します。過去のバージョンの Experiment gate 情報は Cytobank からはアクセスできません。

Apply gates で Experiment gate 情報を保存しない状態で他の機能へ移動しようとするとメッセージが表示されます。

2-8-3 Gating Editor:ゲーティング画面の各機能



① Apply gates ボタン

Gating Editor の状態を Experiment 全体に適用(Apply)するボタンです。 Cytobank では Gating Editor の変更を都度サーバーで保存しています。これに加えて Apply gates ボタンを押した時の情報を Experiment gate 情報としてバージョン保存も行います。 Gating Editor 以外の機能では、保存された最新の Experiment gate 情報のみを使用しま す。過去のバージョンの Experiment gate 情報は Cytobank からはアクセスできません。

- ⑦ ゲーティングツール
 ⑦ プロットにゲートをかけるためのツールです。左から四角形・円形・ポリゴン・クアドラント(四分割)・
 スプリッド(二分割)・レンジゲートです。
- Import gates ボタン
 他の Experiment のゲート情報や Gating-ML をインポートします。
- ④ <u>Cluster gates</u>ボタン クラスターチャネルを用いてクラスターゲートを自動作成します。
- ⑤ Reset all ボタン ゲートやポピュレーションのすべての情報をリセットします。 Undo はできません。
- 6 <u>Compensation</u>

適用されているコンペンセーションマトリックス名が表示されています。鉛筆アイコンをクリックすると Compensation Editor に移動し、Experiment compensation を変更することができます。

⑦ Gates リスト

作成されたゲートのリストが表示されます。Cytobank ではゲートは囲いの XY 座標情報であるだけで、階層構造の情報はここには反映されません。ゲートを組み合わせて使用して定義した階層構造はポピュレーションとして扱います。

Gates 名部分にオーバーカーソルするとメニューが表示され、左から、ゲート名の変更(鉛筆マー

- ク)、<u>Gate tailoring</u>(Tマーク)、<u>Check gates</u>(五角形と四角アイコン)、ロック(錠前マー
- ク)、削除(ゴミ箱アイコン)ができるようになっています。

Filter displayed ボックスで文字列の一致する Gates だけを表示することもできます。

Gating Editor	(i) 🔲 Apply gate
Gates	Q Filter displayed
CD3+	🖉 T 🏠 🗗 🟛
CD4vsCD8	
<u> </u>	

8 Population

プロットに表示する対象のポピュレーションを示しています。クリックすると作成されたポピュレーションの リストが表示され、選択して変更することができます。

9 File

プロット表示しているファイルを示しています。左右の矢印をクリックして表示ファイルを変更することが できます。

10 Plot settings

クリックするとプロット設定を変更できます。



Plot type : プロット種類

Gate label : ゲートとともに表示する情報 Plot colors : カラー表示のパレット Kernel smoothing : スムージングの強度 Percent per contour : 等高線幅の% Outliers start at : 何%以下の密度をドット表示とするか

- Population tree ポピュレーションの階層構造
- Boolean expressions (旧 Population manager 機能を含む) 作成されたポピュレーションの一覧と編集・新規作成ができます。
- <u>Sunburst view</u> ポピュレーションの階層構造と%分布を Sunburst プロットで表示します。

2-8-4プロットに表示されるイベント数が多すぎる場合: Time Slice

FCS ファイルに含まれる細胞イベント数が多すぎてそのままではゲーティングが困難な場合、Time slice をして表示イベント数を減らして下流のゲーティングを行い、作成したゲートを Time slice なしの元ポピュ レーションに適用することができます。

- ① X軸を Time にしたプロットを表示します。
- ② 適当なイベント数になるように Time slice ゲートをかけます。
- プロットの上にある Population メニューで Time slice を選択し、X・Y 軸をゲーティングしたい設定 に変更します。



④ 少なくなったイベント数のプロットでゲーティングを行います。

⑤ Time slice の下に作成されたポピュレーションについて Boolean expression の expression で、設定されている population を Time slice から元ポピュレーションに変更することで、元ポピュレーションでの展開とすることができます。

2-8-5測定中の異常イベントのクリーンアップ

測定時のクロッグ発生や、機器トラブルで途中のデータが正常でないようなデータの場合、安定して測定できている領域にだけ Time slice をかけることで、解析の信頼性を向上させます。



自動検出・除去をしてくれる<u>データ QC:PeacoQC</u>もおすすめです。

2-8-6ゲーティング後のコンペンセーション/チャネル変更: Hidden gate

Cytobank ではゲーティング後にコンペンセーションの変更や修正、<u>チャネルの変更</u>を行うとそのゲートは隠 れゲート(Hidden gate)となり、ゲートが表示されなくなります。 隠れゲートとなったものはゲーティング画面のゲートリストのゲート名に斜線の目玉マークが付きます。コンペンセーションやチャネルをオリジナル(ゲートを作成した時の設定)に戻すと解除されます。

③ Gating Editor	Apply gates		2 + + 2	} €	ን Import / clu	ster gat
Gates	Q Filter displayed	Population	Ungated	-	^ v	
Scatter Ø		File	IL10 - pbmc_lrs005_il10	•	< >	
		260k - 240k -				*
ating Editor	pplygates		Import / cluster gates	Reset all Comp	pensation: File-Int	ternal 🥒
Gates Q. Filter display	File	IL10 - pbmc_lrs005_ii10	< >	Q. Filter displayed		
Warning: Selected gate compe pok' differs from experiment comp	260k - 240k - 240k - 220k - 220k - 200k - 20		•	 Ungated (325,37 Scatter (91.3%) 	74) 5 297,153)	
	180k – 160k –					

Hidden gate をクリックすると、理由がコンペンセーションの場合は赤い Warning メッセージが表示されますので、右上の compensation から、適用する Experiment compensation をオリジナルに戻してください。

理由がチャネルの場合は、X または Y 軸パラメータが Undefined for current panel と表示されます。



2-8-7 Boolean expression:ポピュレーションの定義と編集

Cytobank ではポピュレーションを(複数の)ゲートやポピュレーションの Boolean 式で定義するものと しています。例えばゲート A の集団(ポピュレーション a)の内のゲート B をポピュレーション b とする場合 は、ポピュレーション b =ポピュレーション a AND ゲート B となります。



使用できる演算子はAND、NOT、ORで、カッコを用いた式を作成し複雑なポピュレーションを定義することもできます。

- 2-8-7-1 設定方法
- Gating Editor 画面の右側 Population tree タブと Sunburst view タブの近くにある Boolean expression タブを開きます。
- ② 作成されたポピュレーションとそのポピュレーションを定義する式(Expression)のリストが表示されます。デフォルトではゲート名と同名のポピュレーションが作成されています。Population欄のポピュレーション名で鉛筆マークをクリックすると名前を変更することができます。
- ③ 各ポピュレーションは通常、ポピュレーション・Boolean 演算子・ゲートの組み合わせの式で定義されています。Expression 欄では、ポピュレーションは水色、ゲートは灰色のボックスで表示されています。Expression 欄のポピュレーションをマウスオーバーすると、そのポピュレーションを定義している式が吹き出し表示されます。式に入れられる要素は最大 50 です。



 ④ ポピュレーションの定義を編集するには、Expression欄の各要素をクリックするか右端の+マークを クリックして Expression builder を表示させ、変更または追加したい項目(カッコ、演算子、ポピ ュレーション、ゲート)を選択して式を編集します。

λF	Filter displayed 🕀 Ner	v population		
	Population	Expression		
	CD20+ B cells	CD33- lymphocytes	AND CD	20+ B cells
	CD3+ T cells	CD33- lymphocytes	CD3+ T cells	Expression Builder
	CD3+CD4-T cells	CD3+T cells CD3+CE	D4+ T cells	
	CD3+CD4+ T cells	CD3+T cells CD3+CE	D4- T cells	Parenthesis
	CD33- lymphocytes	Intact cells CD33- lym	nphocytes	()
	CD33+ monocytes	Intact cells CD33+ mo	onocytes	Operator
	Intact cells	Intact cells		AND OR
				AND NOT
				Population
				CD3+ T cells
				CD3+CD4+T cells
				CD3+CD4-T cells
				CD33- lymphocytes
				CD33+ monocytes
				Intact cells .

⑤ リスト上にある New population ボタンをクリックすると新規ポピュレーションを作成して、式を作ることもできます。

Q	-ilter displayed	• New population		-	
	Population	Expression	n		
	Population 1	Đ			
	CD20+ B cells	Expression Builder	× cytes	CD20+ B cells	
	CD3+T cells		cytes	CD3+ T cells	
	CD3+CD4-T cells	Parenthesis	^ CD3+	CD4+ T cells	
	CD3+CD4+ T cells	()	CD3+	CD4- T cells	
	CD33- lymphocytes	Operator	:D33-	lymphocytes	
	CD33+ monocytes	AND OR	:D33+	monocytes	
	Intact cells	AND NOT			
		CD3+ T cells CD3+CD4- T cells CD3+CD4+ T cells CD20+ B cells CD33- lymphocytes CD33+ monocytes	•		

- ⑥ ポピュレーションがたくさんある場合は New population ボタン左の Filter displayed ボックスで、
 文字列の一致するものだけを表示することもできます。
- ⑦ Population 欄のポピュレーション名の鉛筆アイコンをクリックすると名前を変更できます。

Рор	ulation tree (i) Boole	an expression	s Sunburst view		
Q	Filter displayed	• New po	opulation		
	Population		Expression		
	Population 1	I	CD33- lymphocytes	CD20+ B cells	\oplus
	CD20+ B cells	E	CD33- lymphocytes	CD20+ B cells	

⑧ リストの左端にあるボックスにチェックを入れると、複数のポピュレーションにまとめて Prefix:接頭に追加、Suffix:末尾に追加、Copy:コピー、Delete:削除をすることができます。

_	Рор	ulation tree	(i) Boolea	an expressio	ns Sunburst view	
	⊗ (Jnselect NA	Prefix	Ak Suffix	Copy 🛱 De	elete
		Population		*	Expression	
	V	Population 1	L		CD33- lymphocytes	AND CD20+B cells
		CD20+Bce	lls		CD33- lymphocytes	CD20+ B cells

2-8-7-2AND 演算子

階層的なゲーティングストラテジーを作成するには、AND 演算子を用いて親ポピュレーションと新規ゲート を組み合わせます。 典型的な例としては、CD3+ゲートを作成した後にその下に CD4+ゲートを作成した 場合は、CD3+ポピュレーション AND CD4+ゲートの新規のポピュレーション(CD3+CD4+)が作 成されることになります。



また、AND 演算子を用いて複数のゲートの重複する領域に位置する細胞イベントを同定することもでき ます。例えば CD33+ポピュレーションについて、pSTAT3+と pp38+の 2 つをゲーティングして、 CD33+pSTAT3+と CD33+pp38+の 2 つのポピュレーションが作成された後に、CD33+pSTAT3+ ポピュレーションに AND 演算子で pp38+ゲートを追加すると、トリプルポジティブのポピュレーションを作 成することができます。



2-8-7-3NOT 演算子

NOT 演算子はその後ろのポピュレーションに含まれる細胞イベントを除外します。例えば CD33+pSTAT3+とCD33+pp38+の2つのポピュレーションを作成し、



その後、ポピュレーション pSTAT3+ NOT pp38+を作成すると、pSTAT3 だけが陽性の集団とすることができます。



2-8-7-4OR 演算子

OR 演算子はすべてのポピュレーションの細胞イベントを含みます。 例えば CD33+pSTAT3+と CD33+pp38+の2つのポピュレーションを作成し、



その後、ポピュレーション p STAT3+ OR pp38+を作成すると、どちらかまたは両方に属する細胞イベン トすべてを含みます。



2-8-7-5カッコ()を用いた組み合わせ

カッコを用いることで、ひとつの演算子を用いた式だけでなく複雑な式を作成することできます。式に使用で きる要素は 50 個以下です。

2-8-8Check gates:ゲート位置の一括確認

Gate tailoring で Per File (TPF) をしている時など、設定したゲートがすべてのファイルに適切にかかっているかを一括で確認する方法です。

- ① Gates リストでチェックしたいゲートをオーバーカーソルすると、ゲートメニューが表示されます。
- ② メニューの左から3つめの Check gates ボタンをクリックします。



③ 別タブでウィンドウが開きすべてのファイルに適切にゲートがかかっているかを確認することができます。
 ※最後に Apply gates ボタンを押して保存した Experiment gate 情報が読み込まれますので、表示がおかしい場合は Apply gates ボタンを押してから再度確認ください。

2-8-9 Gate tailoring: ファイルや親集団でのゲート位置調整

ゲートを作成すると、デフォルトでは自動でそのゲートをすべてのファイルに適用します(Global mode)。特定のファイルや親ポピュレーションでのゲート位置だけを変更したい場合は Gate tailoring を使用します。

ファイルによってゲート位置を調整したい場合は <u>Per File</u>を、親ポピュレーションによってゲート位置を調整 したい場合は Per Population を使用することでゲート位置を調整することができます。

- 2-8-9-1ファイルによってゲート位置を調整したい場合: Per File (TPF)
- ① 表示するファイルを調整したいファイルに変更します。
- ② Gates リストでゲートをオーバーカーソルし、表示されるメニューの左から2つめTのボタンをクリックし、Per File を Global から Tailored にします。

Gating Editor	i Apply gate
Gates	Gate tailoring
CD4vsCD8 Cells Lympho T Singlets	Gate tailoring () Per File Global Tailored
	Per Population Global Tailored
	Apply tailored gate

- ③ プロット上でゲート位置を変更します。
- ④ 変更後のゲート位置を複数のファイルに適用したい時は Apply tailored gate ボタンを押して、適用したいファイルを選択し、Apply ボタンを押します。
- ⑤ Tailored を解除したい場合は、戻したいゲートのファイルを表示した後、Per File を Global に戻し ます。

2-8-9-2親ポピュレーションによってゲートの位置を調整したい場合: Per Population (TPP) 例えば IL10 刺激前後で pSTAT3 反応する細胞%を CD33+monocytes と CD4+T cells につい て解析したい場合、Global mode では片方に合わせるともう片方が合わなくなってしまいます。



これを解決するため、pSTAT3+ゲートを親ポピュレーションである CD33+monocytes と CD4+T cells とで異なるゲート位置に調整する必要があります。

このような場合に Gate tailoring の Per Population が有用です。

- ① ゲートを作成します。
- ② Gates リストでゲートをオーバーカーソルし、表示されるメニューの左から2つめTのボタンをクリックし、Per Population を Global から Tailored にします。



- ③ Population を調整したい親ポピュレーションに変更します。
- ④ プロット上でゲート位置を変更します(変更しないと正常に機能しません)。
- ⑤ 変更後のゲート位置を他の親ポピュレーションに適用したい時は Apply tailored gate ボタンを押して、適用したい親集団を選択し、Apply ボタンを押します。

pStat3+ T	Î	Т 🗘 🗹 🖬	Ô			
	Gate tailoring (i)					
	Per File					
	Global	Tailored				
	Per Population	on				
	Global	Tailored				
	Apply tailored gate					
Apply Tailored Gate						
Selected Gate: pStat	3+					
Tailored to Population: Select Population(s) to apply the tai Filter displayed:	CD33+ monocytes	All	None			
CD3+CD4+ T cells Intact cells Ungated			*			
0 selected	Cancel	v				

⑥ 各親ポピュレーションの下に TPP ポピュレーションが作成されます。 Population tree や Boolean expression タブで確認することができます。
 ※TPP は Boolean expression で編集することはできません。編集が必要な場合は一度 Global に戻す必要があります。

i Population tree	Boolean expressions	Sunburst view
Q, Filter displayed		
 Ungated (326,6 	23)	
 Intact cells (8 	8.6% 289,278)	
 CD3+CD4+ 	T cells (38.2% 110,53	9)
pStat3+ (CD3+CD4+ T cells) (0.0	033% 36)
 CD33+ mor 	nocytes (12.8% 37,114	ł)
pStat3+ (CD33+ monocytes) (0.3	16% 61)

⑦ Illustrations 機能で並べて表示することもできます。



2-8-9-3ファイルと親ポピュレーションの両方によってゲート位置を調整する:TPF&TPP 厳密なゲーティングが必要な場合は、ファイルと親ポピュレーションの両方によって異なるゲート位置が必要 になります。そのような場合は、TPFとTPPを組み合わせで使用すると良いです。

 Gates リストでゲートをオーバーカーソル> Gate tailoring ボタンを選択> Per File と Per Population の両方を Tailored にします。

pStat3+ T	0				
	Gate tailoring (i) Per File				
	Global	Tailored			
	Per Population				
	Global	Tailored			
	Apply tail	pred gate			

- この後変更したすべてのゲートは、そのファイル・親ポピュレーションの場合での位置となります。設定 原理は TPF、TPP と同じです。
- ③ 位置を調整したゲートを複数のファイル・親ポピュレーションの場合に使用したい時は、Apply tailored gate ボタンを押します。



④ ファイルと親ポピュレーションのリストからゲートを適用したいものを選択して Apply ボタンをクリックしま す。

Apply Tailored Gate	
Selected Gate: pStat3+ Tailored to Fcs File: Unstim1 - pbmc_lrs005_unstim1 Select Fcs File() to apply the tailored gate to: Filter displayed:	All None
IL10 - pbmc_Irs005_II10 IL6 - pbmc_Irs005_II6 LPS - pbmc_Irs005_Ips Unstim1 - pbmc_Irs005_unstim1 Unstim2 - pbmc_Irs005_unstim2	*
2 selected Tailored to Population: CD3+CD4+ T cells Select Population(2) to apply the tailored gate to: Filter displayed:	All None
CD3+CD4+ T cells CD33+ monocytes Intact cells Ungated	A V
1 selected Cancel Apply	

2-8-9-4複数の子ポピュレーションをまとめて他の親ポピュレーションの下にも展開する

定量的な比較を行う場合、ある親ポピュレーションで展開した子ポピュレーション解析を他の親ポピュレーションでも展開したいことがよくあります。

例えば、CD19+B cells 下の 3 つの機能性マーカーの展開を、CD4+T、CD8T、CD33+Mono でも 展開し、%ポピュレーションを調べたい場合などです。

i Population tree	Boolean expressions	Sunburst view			
Q Filter displayed					
 Ungated (351,61 	19)				
 Singlets (56.69) 	% 198,998)				
▼ CD45+ (65.	4% 130,086)				
 CD3+ T cells (56.4% 73,411) 					
CD4+ T	cells (50.5% 37,083)				
CD8+ T	cells (33.5% 24,573)				
(• CD19+B	cells (10.9% 14,222)				
PERK1/	2+ (1.65% 235)				
pSTAT3	+ (3.6% 512)				
pSTAT5	+ (3.56% 506)				
CD33+ M	onocytes (21% 27,304	1)			

Boolean expression で、3つの子ポピュレーションをコピーして編集する方法もありますが、TPPと Apply tailored gate 機能での以下の方法もあります。

 子ポピュレーション(例えば pSTAT5+)のゲートを Gates リストで選択し、Gate tailoring の Per Population を Tailored にします。

- Apply tailored gate 機能で CD4+T、CD8+T、CD33+Mono を TPP の対象親ポピュレーションにして Apply ボタンを押します。
- ③ すべての親ポピュレーションの下に pSTAT5+が作成されたことを確認します。



④ 残りの2つの子ポピュレーションについても同様にして作成します。



⑤ Cytobankのゲートとポピュレーションの定義のとおり、TPPで自動作成されたポピュレーションはその後独立するため、もし子ポピュレーションの位置を親ポピュレーションによって微調整せず同じにしたい場合は、適用したいゲート位置のものをプロットに表示した後に、Per Populationの TailoredをGlobal に戻すことですべての親ポピュレーションで同じゲート位置になります。

※TPP が適用されているポピュレーションはその下流も含めて Boolean expression で編集ができなくなる仕様です。

Pop	ulation tree () Boolean expressions Sunburst view	
Q	Filter displayed New population	
	Population *	Expression
	CD19+ B cells	CD45+ CD19+ B cells
	CD3+T cells	CD45+ CD3+T cells
	CD33+ Monocytes	CD45+ CD33+ Monocytes
	CD4+ T cells	CD3+T cells CD4+T cells
	CD45+	Singlets CD45+
	CD8+T cells	CD3+T cells CD8+T cells
	pERK1/2+ (CD19+ B cells)	CD19+ B cells pERK1/2+
	pERK1/2+ (CD33+ Monocytes)	CD33+ Monocytes pERK1/2+
	pERK1/2+ (CD4+ T cells)	CD4+T cells pERK1/2+
	pERK1/2+ (CD8+ T cells)	CD8+T cells pERK1/2+
	pSTAT3+ (CD19+ B cells)	CD19+ B colls pSTAT3+
	pSTAT3+ (CD33+ Monocytes)	CD33+ Monocytes pSTAT3+
	pSTAT3+ (CD4+ T cells)	CD4+T cells pSTAT3+
	pSTAT3+ (CD8+ T cells)	CD8+T cells pSTAT3+
	pSTAT5+ (CD19+ B cells)	CD19+ 8 cells pSTAT5+
	pSTAT5+ (CD33+ Monocytes)	CD33+ Monocytes pSTAT5+
	pSTAT5+ (CD4+ T cells)	CD4+T cells pSTAT5+
	pSTAT5+ (CD8+ T cells)	CDB+T cells pSTAT5+
	Singlets	Singlets

2-8-10ポピュレーションの階層構造の確認と編集

Cytobank ではゲートを作成すると同時に同名のポピュレーションを作成し、作成されたポピュレーションは ポピュレーションとして扱います。ゲート内容(位置やサイズの変更など)とポピュレーションは連動します が、ゲート自体を削除しても<u>ポピュレーションは削除</u>されなかったり、ゲート名を<u>変更してもポピュレーション</u> 名はそのままだったりします。

この場合は、<u>Population tree</u>タブや <u>Boolean expression</u>タブでポピュレーションを選択し、Delete ボタンでポピュレーションを削除したり、ポピュレーション名の部分の鉛筆マークをクリックして名前を変更した りします。

2-8-10-1 Boolean expression: ポピュレーション一覧とその編集

Gating Editor 右の Boolean expression タブを開くと Population と Expression の表が表示されます。ここでは各ポピュレーションの削除・コピー・新規作成や名前の変更、ポピュレーションの定義(他のポピュレーション、ゲート、Boolean 演算子で構成される式)を編集することができます。

ポピュレーションがたくさんある場合は New population ボタン左の Filter displayed ボックスで、文字 列の一致するものだけを表示することもできます。

詳細は Boolean expression を参照ください。

Poj	Dulation tree (1) Boolean expression	Sunburst view	
٩	Filter displayed New	population	
	Population	Expression	
	CD20+ B cells	CD33- lymphocytes AND CD	20+ B cells
	CD3+ T cells	CD33- lymphocytes CD3+ T cells	Expression Builder 🛛 🗙
	CD3+CD4- T cells	CD3+T cells CD3+CD4+T cells	
	CD3+CD4+ T cells	CD3+T cells CD3+CD4-T cells	Parenthesis
	CD33- lymphocytes	Intact cells CD33- lymphocytes	(
	CD33+ monocytes	Intact cells CD33+ monocytes	Operator
	Intact cells	Intact cells	AND OR
			AND NOT Population CD3+ Teells CD3+CD4+ Teells CD3+CD4+ Teells CD33+ monocytes Lttatr cells

2-8-10-2 Population tree: ポピュレーション階層構造表示 Gating Editor でゲートを作成すると、同名のポピュレーションも自動的に作成され、その後独立に扱う ことができます。中央のプロットを挟んで左側がゲーティング、右側がポピュレーションとなっています。



ポピュレーション側はデフォルトでは Population tree タブが表示されており、ポピュレーションの階層構造 が表示されています。階層レベルに沿って名前、% of parent、細胞数が表示されています。 オーバーカーソルすると編集メニューが表示され、名前の変更やコピー、削除をすることもできます。 クリックで複数選択もでき、上部に表示されるメニュー(Prefix:接頭に追加、Suffix:末尾に追加、 Copy:コピー、Delete:削除)で一括編集もできます。



ここで表示されるポピュレーション名、% of parent、細胞数の一覧の数値情報が必要な場合は Export statisticsで csv ファイル出力できます。

2-8-10-3Sunburst 表示

Population Sunburst タブを開くとポピュレーション階層と統計情報を多重円グラフで表示します。

- ① オーバーカーソルでそのポピュレーションの数値情報を表示することができます。
- ポピュレーションをクリックすると、そのポピュレーションを360°で表し、下層の構成比率を感覚的に 把握しやすくします。
- ③ 上位層を360°にしたい場合は左上のバーをクリックします。



④ 左にある Palette:で表示するカラーパレットを変更できます。デフォルトは blue になっています。



- ⑤ Print view ボタンでは別タブで Sunburst だけを表示します。
- ⑥ Export to PDF ボタンでは、PDF ファイルとしてダウンロードできます。

Blues, Orang	es, & Greens 🔻
Display labels:	
Print view	Export to PDF

2-8-11ゲーティングストラテジーの自動表示: Gating hierarchy

表示したいファイルとポピュレーションを指定すると、そのポピュレーションがどのようなゲーティングストラテジー で作成されたのか、上流のゲーティングを自動的に一括表示することができます。

展開図を作成する手間を省くことができるだけでなく、途中のゲーティングを取り違えることがないため、確認だけでなく、レビュアーによる評価にも便利です。

- ① メニューバー> Gates> Gating hierarchy を選択します。
- ② FCS Files ボックスの Choose をクリックし、表示したいファイルを選択します。
- ③ Populations ボックスの Choose をクリックし、確認したいポピュレーションを選択します。
- ④ 選択したポピュレーションに至るゲーティングストラテジーがファイルごと、ポピュレーションごとにすべて表示されます。
- ⑤ 出力が必要な場合は Print view ボタンで別タブ表示後、ブラウザの印刷機能で PDF 出力もできます。



2-8-12ゲート情報のインポート

インポートできるゲート情報は <u>Experiment gate</u>情報です。 インポート元とコンペンセーションとチャネルの設定が異なる場合は Hidden gate となります。

- 2-8-1 2-1 他の Experiment のゲート情報をインポート
- Gating Editor メニュー> Import Gates> Import from another experiment をクリックします。

(i) Gating Editor	(i) 🔳 App	ply gates	口口公-	† -∔ ŀ·			☆ Import gates *	🏠 Cluste	er gate
Gates	C Filter displayed	Population	Singlets	•	^	~	☆ Import from another e ☆ Upload gating-ML	xperiment:	(i) es
		File	Sample 1 IM Basic	•	<	>			

 インポート元となる Experiment の名前か Experiment ID で検索し、選択して Import Gates ボタンを押します。

Import gates from experiment				;	
Selec	t an experime	nt to import fr	om:		
Ent	er name or sele	ict		-	

- ③ インポート結果が表示されます。
- ④ Gating Editor に移動するとインポートされたゲートが確認できます。
- ⑤ 問題なければ Apply gates ボタンを押して Experiment gate 情報として保存します。

※現時点では、ゲート情報はすべて Global としてインポートされ、Tailored gates は解除されます。

Gates リストの T マークも表示されなくなります。テンプレートとして使用して、Gate tailoring での微調 整や、<u>Boolean expression</u>での編集が必要な場合もあります。

Tailored gates もインポートで反映したい場合は、Gating-ML でインポートしてください。

2-8-1 2-2 Gating-ML のゲート情報をインポート

Cytobank では、International Society for Advancement of Cytometry (ISAC) による Gating-ML 2.0 の規格を採用しています。

 Gating Editor メニュー> Import Gates> Upload gating-ML または、Actions> Upload> Upload gating-ML をクリックします。

	😑 Actions 🗸 👌 🛉 Sample tag	s 🎶 Data QC 🏠 Gates 🚓 Ad	dvanced analyses 👘 🗽 Illustrations			
<	 Experiment actions View summary 	Apply gates	$\Box \circ \diamond + +$		C	â Import gates
	🚳 Clone	Population	intact cells	~	*	
	⊥ Upload	🎲 Upload sample tags 🛛 💡 File	IL10 - pbmc_lrs005_i110 •	<	>	
	📩 Export 🔋	1 Upload gates from gating-ML				
	P Link to parent	Upload more files				*
	🗊 Reset	220k-				
	💼 Delete	4 200k-				
	gate B	180k-				

- ② Gating-ML ファイルを指定してアップロードします。
- ③ インポート結果が表示されます。

④ Gating Editor に移動するとインポートされたゲートが確認できます。

⑤ 問題なければ Apply gates ボタンを押して Experiment gate 情報として保存します。 Gating-ML ファイルのエクスポート方法もご参照ください。その際、Experiment gate_の情報が Gating-ML としてバージョン保存されることにご留意ください。

※Gating-ML ファイルは text editor で開いて編集が可能です。編集が必要な場合はサポートまでご 相談ください。

2-9 Sample tags:サンプルタグ

2-9-1サンプルタグとは

サンプルファイルの背景(疾患状態や投与化合物、タイムポイントなどの実験変数)の情報をサンプルタ グとして付加することができます。FCS ファイルにサンプルタグを付加することで、illustrations のレイアウト 項目として使用することができるので、統計比較をより簡単に迅速に行うことができます。



Experiment 内のサンプルタグは、Advanced analysis の結果 Experiment や、クローンにも引き 継がれるのでデータをアップロードした段階でサンプルタグをつけておくことをお勧めします。

2-9-2サンプルタグのつけ方

※事前にファイルの情報にどのようなサンプルタグを付加するかを決定しておくことをおすすめします。

- ① メニューバーの Sample Tags を開き、Add experiment dimensions をクリックします。
- ② Add experiment dimensions ウィンドウには Doses、Individuals、Timepoints などがデフ オルトとして用意されています。その他任意のサンプルタグを Define a custom dimension で設 定することができます。左でアイコンを、中央にカテゴリー名を入力し、Submit ボタンを押します。


- ③ Add カテゴリー名欄に設定したい条件(例タイムポイントの場合、経過時間など)をサンプルタグと してコンマ区切りで入力します。
- ④ 各サンプルタグが作成されますので、該当するファイルをドラッグ&ドロップまたは Move file to で移動して分類します。※入力した条件名とファイル名が部分一致すると自動で振り分けられます。
- ⑤ ② ④を繰り返して複数のサンプルタグをファイルに付加し、Illustrationsのレイアウトに用いることができます。

Add Timepoints	0	
Enter a comma separated list of Timepoir	nts to add:	
15min,30min		Add Timepoints
All Timepoints 15min 30min C All Timepoints Tag files with Timepoints by dragging them to	o the box for the desired tag. You can also filt	er the files displayed and move them in bulk
Untagged	15min Tagged Files	30min Tagged Files
Filter Order Files to	Patient_female01-15min-dose01 Patient_female01-15min-dose02 Patient_male01-15min-dose01 Patient_male01-15min-dose02 All files that are affected by the 15 minute time point are tagged as such	Patient_female01-30min-dose01 Patient_female01-30min-dose02 Patient_male01-30min-dose01 Patient_male01-30min-dose02 All files that are affected by the 30 minute time point are tagged as such

2-9-3他の Experiment からのサンプルタグのインポート

- ① 元 Experiment での dimension と同一名の dimension を Sample tags> Add experiment dimensions で作成します。
- ② Import from another experiment をクリックし、元 Experiment を指定して適用します。

d Clusters	0	Cluster Actions
nter a comma separated list of Cluste	rs to add:	
Cluster 1, Cluster 2, Cluster 3, etc.	the Theorements of the second analyses is in Institutions in the Institutions is even the Clusters is even the Clusters is add: Type into the box to find an existing experiment Sters Clusters by Angeing the to the box for the desired tag. You can also there the files displayed and move them in bulk using Move files ta.: Editors by Angeing the to the box for the desired tag. You can also there the files displayed and move them in bulk using Move files ta.: Editors by Angeing the to the box for the desired tag. You can also there the files displayed and move them in bulk using Move files ta.: Editors by Angeing the to the box for the desired tag. You can also there the files displayed and move them in bulk using Move files ta.: Editors by Angeing the to the box for the desired tag. You can also there the files displayed and move them in bulk using Move files ta.: Editors by Angeing the to the box for the desired tag. You can also there the files displayed and move them in bulk using Move files ta.: Editors by Angeing the to the box for the desired tag. You can also there the files displayed and move them in bulk using Move files ta.: Editors by Angeing the to the box for the desired tag. You can also there the files displayed and move them in bulk using Move files ta.: Editors by Angeing the to the box for the desired tag. You can also there the files displayed and move them in bulk using Move files ta.: Editors by Angeing the to the box for the desired tag. You can also there the files displayed and move them in bulk using Move files tag.: Editors by Angeing tag.:	Import Cluster tags from an experiment
	(Fluorescence) Sample Lags	Enter name or select
All Clusters	Type into the box to find an existing experiment	Import
Untagged		pbmc_lrs005_il10.fcs •
Untagged Fite Move files to V Move PBMC_LRS005_JL101/sts (IL10) genc/rs005_J101/st PBMC_LRS005_JL6.fts (IL6) ster (rs005 files)		pbmc_Ins005_II10.fos • Duplicate
Untagged Filter [Move files Is		Dupficate
Untagged Fitter Move fites to		pbmc, Iva005, IVA.006 • Duplicate

2-9-4サンプルタグの一括編集・追加

ファイル数や追加するサンプルタグが多い場合、下記の手順でサンプルタブを一括編集・追加をすることができます。

 Actions> Export> Export sample tags でファイルとデフォルトサンプルタグのスプレッドシートを エクスポート、ダウンロードします。



 ダウンロードしたファイルをエクセル等で開き、追加したい dimension や tags を表に追加、編集 し、.txt か.csv、.tsv 形式で保存します。

_	A	B	C	D	E	F	G	Н		J	K	L	M
1	FCS Filename	Conditions	Doses	Doses Nume	Doses Units	Timepoints	Individuals	Plate Shape	Sample Type	Plate Row	Plate Column	Plate	FCS File Category
2	comp_alexa 488.fcs	-	-	-	-	-	-	-	-	-	-	Comp	Compensation Control
3	comp_alexa 647.fcs	-	-	-	-	-	-	-	-	-	-	Comp	Compensation Control
4	comp_pacblu.fcs	-	-	-	-	-	-	-	-	-	-	Comp	Compensation Control
5	comp_pe.fcs	-	-	-	-	-	-	-	-	-	-	Comp	Compensation Control
6	comp_pe_cy7.fcs	-	-	-	-	-	-	-	-	-	-	Comp	Compensation Control
7	comp_percp_cy55.fcs	-	-	-	-	-	-	-	-	-	-	Comp	Compensation Control
8	pbmc_lrs005_il10.fcs	-	-	-	-	-	-	-	-	-	-	PBMC_LRS005	Experiment Files
9	pbmc_lrs005_il6.fcs	-	-	-	-	-	-	-	-	-	-	PBMC_LRS005	Experiment Files
10	pbmc_lrs005_lps.fcs	-	-	-	-	-	-	-	-	-	-	PBMC_LRS005	Experiment Files
11	pbmc_lrs005_unstim1.fcs	-	-	-	-	-	-	-	-	-	-	PBMC_LRS005	Experiment Files
12	pbmc_lrs005_unstim2.fcs	-	-	-	-	-	-	-	-	-	-	PBMC_LRS005	Experiment Files
	A	В	С	D	E	F	G	Н	1	J	K	L	М
1	A FCS Filename	B Conditions	C Doses	D Doses Nume	E Doses Units	F Timepoints	G Individuals	H Plate Shape	l Sample Type	J Plate Row	K Plate Columr	L	M FCS File Category
1 2	A FCS Filename comp_alexa 488.fcs	B Conditions -	C Doses	D Doses Nume -	E Doses Units -	F Timepoints -	G Individuals -	H Plate Shape -	l Sample Type -	J Plate Row -	K Plate Columr -	L Plate Comp	M FCS File Category Compensation Control
1 2 3	A FCS Filename comp_alexa 488.fcs comp_alexa 647.fcs	B Conditions -	C Doses - -	D Doses Nume - -	E Doses Units - -	F Timepoints -	G Individuals - -	H Plate Shape - -	l Sample Type - -	J Plate Row - -	K Plate Columr - -	L Plate Comp Comp	M FCS File Category Compensation Control Compensation Control
1 2 3 4	A FCS Filename comp_alexa 488.fcs comp_alexa 647.fcs comp_pacblu.fcs	B Conditions - - -	C Doses - -	Doses Nume - -	E Doses Units - -	F Timepoints - -	G Individuals - -	H Plate Shape - -	l Sample Type - -	J Plate Row - -	K Plate Columr - -	L Plate Comp Comp Comp	M FCS File Category Compensation Control Compensation Control Compensation Control
1 2 3 4 5	A FCS Filename comp_alexa 488.fcs comp_pachu.fcs comp_pachu.fcs comp_pe.fcs	B Conditions - - -	C Doses - - -	Doses Nume - - -	E Doses Units - - - -	F Timepoints - - -	G Individuals - - -	H Plate Shape - - - -	I Sample Type - - -	J Plate Row - - -	K Plate Columr - - -	L Plate Comp Comp Comp Comp	KCS File Category Compensation Control Compensation Control Compensation Control Compensation Control
1 2 3 4 5 6	A FCS Filename comp_alexa 488.fcs comp_alexa 647.fcs comp_pacblu.fcs comp_pe.fcs comp_pe_cry7.fcs	B Conditions - - - -	C Doses - - - - -	D Doses Nume - - - -	E Doses Units - - - -	F Timepoints - - - -	G Individuals - - - - -	H Plate Shape - - - - -	I Sample Type - - - -	J Plate Row - - - -	K Plate Columr - - -	L Plate Comp Comp Comp Comp Comp	M FCS File Category Compensation Control Compensation Control Compensation Control Compensation Control
1 2 3 4 5 6 7	A FCS Filename comp_alexa 488.fcs comp_aelxa 647.fcs comp_petits comp_pe.fcs comp_pec.fcs comp_pec_cy7.fcs comp_percp_cy55.fcs	B Conditions - - - - -	C Doses - - - - -	Doses Nume - - - - -	E Doses Units - - - - -	F Timepoints - - - - -	G Individuals - - - - -	H Plate Shape - - - - - -	I Sample Type - - - - -	J Plate Row - - - - -	K Plate Columr - - - - -	L Plate Comp Comp Comp Comp Comp Comp	M FCS File Category Compensation Control Compensation Control Compensation Control Compensation Control
1 2 3 4 5 6 7 8	A FCS Filename comp_alexa 488.fcs comp_pachuf.fcs comp_pachU.fcs comp_percs comp_percs_v55.fcs phmc_trs05_110.fcs	B Conditions - - - - - IL10	C Doses - - - - -	Doses Nume - - - - - - -	E Doses Units - - - - - - -	F Timepoints - - - - - -	G Individuals - - - - - -	H Plate Shape - - - - - - - -	l Sample Type - - - - - -	J Plate Row - - - - - -	K Plate Columr - - - - - -	L Plate Comp Comp Comp Comp Comp Comp PBMC_LRS005	M FCS File Category Compensation Control Compensation Control Compensation Control Compensation Control Compensation Control Experiment Files
1 2 3 4 5 6 7 8 9	A FCS Filename comp_alexa 488.fcs comp_pacblu.fcs comp_pacblu.fcs comp_pe_cry7.fcs comp_pe_cry7.fcs comp_percp_cry55.fcs pbmc_trs005_ill0.fcs	B Conditions - - - - - - - - - - 1L10 IL6	C Doses - - - - - -	Doses Nume - - - - - - - - - - - - - -	E Doses Units - - - - - - - - - -	F Timepoints - - - - - - - -	G Individuals - - - - - - - - -	H Plate Shape - - - - - - - - - - - - - -	I Sample Type - - - - - - - - - - - - -	J Plate Row - - - - - - - - - - - -	K Plate Columr - - - - - - - - - - - - - - - -	L Plate Comp Comp Comp Comp Comp Comp PBMC_LRS005 PBMC_LRS005	M FCS File Category Compensation Control Compensation Control Compensation Control Compensation Control Compensation Control Experiment Files
1 2 3 4 5 6 7 8 9 9 10	A FCS Filename comp_alexa 488.fcs comp_pacblu.fcs comp_pachU.fcs comp_pe.cy7.fcs comp_pe.cy7.fcs comp_percy.cy55.fcs pbmc_irs005_ill0.fcs pbmc_irs005_il6.fcs	B Conditions - - - - - - - - - - IL10 IL6 LPS	C Doses - - - - - - - - - - -	Doses Nume - - - - - - - - - - - - -	E Doses Units - - - - - - - - - - - - -	F Timepoints - - - - - - - - - -	G Individuals - - - - - - - - - - -	H Plate Shape - - - - - - - - - - - -	I Sample Type - - - - - - - - - - -	J Plate Row - - - - - - - - - - - - - - -	K Plate Columr - - - - - - - - - - - -	L Plate Comp Comp Comp Comp Comp PBMC_LRS005 PBMC_LRS005 PBMC_LRS005	M FCS File Category Compensation Control Compensation Control Compensation control Compensation control Compensation Control Experiment Files Experiment Files
1 2 3 4 5 6 7 8 9 10 11	A FCS Filename comp_alexa 488.fcs comp_pacblu.fcs comp_perg_ory7.fcs comp_percp_cy55.fcs pbmc_lrs005_i10.fcs pbmc_lrs005_jls.fcs pbmc_lrs005_jns.fcs pbmc_lrs005_jns.fcs	B Conditions - - - - - IL10 IL6 LPS Unstim	C Doses - - - - - - - - - - - - -	Doses Nume	E Doses Units - - - - - - - - - - - -	F Timepoints - - - - - - - - - - - - -	G Individuals - - - - - - - - - - -	H Plate Shape - - - - - - - - - - - -	l Sample Type - - - - - - - - - - - - - -	Plate Row - - - - - - - - - - - - -	K Plate Columr - - - - - - - - - - - - -	L Plate Comp Comp Comp Comp Comp PBMC_LRS005 PBMC_LRS005 PBMC_LRS005	M FCS File Category Compensation Control Compensation Control Compensation Control Compensation Control Experiment Files Experiment Files Experiment Files

- ③ Cytobank に戻り、②で dimension を追加した場合は、同名の dimension を Sample tags
 > Add experiment dimensions で作成しておきます。
- ④ Actions>Upload>Upload sample tags で②で保存したファイルをアップロードして読み込み ます。



2-9-5チャネルとパネル設定: Assign panels

2-9-5-1 チャネルとパネルとは

チャネルとはデータ取得時の機器チャネルです。機器の種類やメーカーによって定義は異なりますが、FCS ファイルには2つの情報が記載されています。1つは detector name で Channel Name または short channel name と呼ばれます。もう1つは marker/reagent name で Long channel name とも呼ばれます。

Channel Name	Marker/Reagent
SSC-A	SSC-A
Ax488-A	pStat3
PE-A	CD33
FL3-A	empty
PerCP-Cy55-A	CD20
PE-Cy7-A	CD3

Detector name/Channel Name/Short channel name は機器が自動的につけるもので、通常 は変更できません。Marker/Reagent name/Long channel name は通常測定時に任意に設定 できるもので、オペレーターが実験や研究対象によって自由に設定しています。

この Channel Name と Marker/Reagent のセットをパネルとしています。 Cytobank では上図のようにセットの表で示し、 FCS ファイルをアップロードした時、 そのファイルに記載されている 2 つのチャネル情報を読み取り、同一のパネルを持つファイルをパネルごとにまとめます。 FCS ファイルに記載されているチャネルの数や 2 つの名前が同一でない場合、パネル/チャネルの不一致となります。

- 2-9-5-2パネル/チャネルの確認と編集
- ① メニューバー> Sample tags> Assign panels を選択します。

PBMC Experim	ment (Fluorescence)			
Actions	🌱 Sample tags 🗸 🛛 ۸- Data QC	🗘 Gates	ភ្នំ Advanced analyses	L Illustrations
? Experim	Assign panels (1)	Assign files to panels and adjust	etails	
	III Plate Annotator	t (Fluorescence)		
	Edit Condition tags (4)	: (Fluorescence)		

② Experiment 内のパネルが表示されます。Experiment 内のパネルが自動作成されており、チャネル数と Name (= Detector name/Channel Name/Short channel name)、Reagent (= Marker/Reagent name/Long channel name)が表として表示されます。また、同一のパネルを持つファイルは各パネルボックス下部にまとめられています。

③ ファイル名左のボックスにチェックを入れて選択し、Move ボタンをクリックすると New panel (新規のパネルを作成し移動)や既存の別のパネルに移動を選択することができます。既存の別のパネルに移動したい場合は、移動先のパネルに表示される Move to…ボタンでも可能です。

Panel 1 🥖		\oplus	Panel 2 🥖	
18 channels		٩	18 channels	٩
Name	Reagent		Name	Reagent
FSC-A	FSC-A		FSC-A	FSC-A
FSC-W	FSC-W		FSC-W	FSC-W
SSC-A	SSC-A		SSC-A	SSC-A
PE-A	CD33		PE-A	CD33
PE-Cy7-A	CD3		PE-Cy7-A	CD3
Time	Time		Time	Time
Ax488-A	pStat3		Ax488-A	pStat3
PE-TR-A	PE-TR-A		PE-TR-A	PE-TR-A
PerCP-Cy55-A	CD20		PerCP-Cy55-A	CD20
Ax647-A	pp38		Ax647-A	pp38
Ax700-A	Ax700-A		Ax700-A	Ax700-A
Ax750-A	Ax750-A		Ax750-A	Ax750-A
PacBlu-A	CD4		PacBlu-A	CD4
Qdot525-A	Qdot525-A		Qdot525-A	Qdot525-A
PacOrange-A	CD8		PacOrange-A	CD8
Qdot605-A	Qdot605-A		Qdot605-A	Qdot605-A
Qdot655-A	Qdot655-A		Qdot655-A	Qdot655-A
Qdo Move 1 s	elected file to		Qdot705-A	Qdot705-A
Over 1 s	nel		Move	to Panel 2 🔶
I M Comper	sation Files		Unstim1 - pbr	mc_lrs005_unstim1.fcs
I Panel 2				
🗹 Unstim2 - pbn	nc_lrs005_unstim2	.fcs		

④ デフォルトでは含まれる全てのパネルが visible となっており、解析に使用できるようになっています。
 特定のパネルの visible のチェックを外すとそのパネルに属するファイルは解析に使用できなくなります。

😰 Assign par	nels 🛈	Config	gure control files				😰 Assign par	els 🛈 🛛 🛛) Config	ture control files				
Panel 1			Panel 2		Compensation files		Panel 1			Panel 2		U Visible	Compensation files	
18 channels		9,	18 channels	Visible panels	🗆 6 files	٩	18 channels		9,	18 channels		9,	🔲 6 files	9,
Name	Reagent		Name	Real analysis in the	Alexa 488 - comp_alexa 48	88	Name	Reagent		Name	Reagent		Alexa 488 - comp_alexa 488	3
FSC-A FSC-W SSC-A Time Ax488-A PE-A PE-TR-A PE-CP-Cy55-A PE-CY7-A Ax647-A Ax700-A Ax750-A	FSC-A FSC-W SSC-A Time pStat3 CD33 PE-TR-A CD20 CD3 pp38 Ax700-A Ax750-A		FSC:A FSC:W SSC:A Time Ax488:A PE:-A PE:-R-A PE:-Cy7:-A Ax647:-A Ax700-A Ax750-A	FSC 4 - 000-mmm FSC W SSC A - 1 Time pStat3 cd33 PE-TR-A CD20 CD3 pp38 Ax700-A Ax700-A	Alexa 647-comp.alexa 64 PE-comp.ge PE_Cy7-comp.ge.cy7 PE_Cy7-comp.ge.cy7 PacBu-comp.gatchu PerCP_Cy55-comp.percp	47 p_cy55	FSC-A FSC-W SSC-A Time Ax488-A PE-A PE-A PE-TR-A PE-CP-Cy55-A PE-Cy7-A Ax647-A Ax700-A	FSC-A FSC-W SSC-A Time pStat3 CD33 PE-TR-A CD20 CD3 pp38 Ax700-A Ax750-A		FSC:A FSC:W SSC:A Time Av488-A PE:-A PE:-R-A PE:-Cy7-A Av467-A Av700-A Av700-A	FSC-A FSC-W SSC-A Time pStat3 cd33 PE-TR-A CD20 CD3 pp38 Ax700-A Ax750-A		Alexa 647 - comp. alexa 647 PE - comp.pe (sy7 PE(sy7 - comp.pe, cy7 PacBlu - comp.pacblu PerCP_Cy55 - comp.percp.	r cy55
□ 4 files		Q	One file	Q			□ 4 files		Q	One file		Q,		
 IL10 - pbmc_li IL6 - pbmc_lrs LPS - pbmc_lr Unstim2 - pbr 	rs005_il10 :005_il6 s005_lps mc_lrs005_uns	tim2	Unstim1 - pbn	nc_Irs005_unstim1			 IL10 - pbmc_lit IL6 - pbmc_lrs LPS - pbmc_lrt Unstim2 - pbm 	rs005_il10 005_il6 s005_lps nc_lrs005_unstin	n2	Unstim1 - pbr	nc_Irs005_ur	stim1		

- ⑤ パネルの Reagent 欄は Marker/Reagent name/Long channel name を読み取った情報 が表示されているため、オーバーカーソルで表示される鉛筆マークをクリックすることで名前を任意に変 更することができます。
- ⑥ パネルの削除はできません(属する FCS ファイルがなくなった場合自動でパネルもなくなります)。

2-9-5-3パネル/チャネルの不一致の解消方法

Cytobankの解析では基本的に同一のパネルであることを前提としています(Advanced Analysis では設定で使用するチャネルが同一であれば実行できるものも例外としてあります。)。下図のようなメッセージが表示されたら、パネル/チャネルの不一致があります。

There are currently panels that contain conflicting channels. These conflicts must be resolved $^{\times}$ before illustrations may be drawn.

不一致のケースとその解消は以下の4つとなります。Assign panels 機能で行います。

- 測定時に設定した Reagent 名の入力ミスや大文字小文字、スペースの有無、半角全角の違い による不一致の場合(例"CD38" vs "cd38")
 - ・ メッセージをクリックすると Assign panels に移動し、解消提案が表示されますので、参考にしてパネルを同一にします。
- ある Channel name について異なるマーカーを使用したファイルが含まれているため Reagent が 不一致の場合(例"pSTAT3"と"pp38")
 - ・ パネルを分けます(または別のパネルを新規作成してそこへ片方を移動)。
 - 両方を含む Regent 名(例"pSTAT3&pp38")のパネルを作成してまとめます。どちらの Reagent かはファイル名や下流の解析で判断します。
- チャネル数がファイルによって不一致の場合
 - ・ パネルを分ける(または別のパネルを新規作成してそこへ片方を移動)。
 - ・ 機器から FCS ファイルを出力し直します。
 - ・ <u>サポートまでお問い合わせ</u>ください。
- Reagent は同一だが、Channel name が不一致の場合(例"FL1" vs "FL-1")
 機器ソフトウエアのバージョンアップ等によって出力される Channel name 名 (Detector name/Channel Name/Short channel name) が変わってしまった場合など
 - ・ パネルを分ける(または別のパネルを新規作成してそこへ片方を移動)。
 - ・ 機器から FCS ファイルを出力し直します。
 - サポートまでお問い合わせください。

2-9-6ファイル属性の管理

2-9-6-1ファイル属性とは

Cytobank では FCS ファイルを下記の3つの属性に分類します。

- 1. Experiment Files: Experiment のすべての機能で表示され、使用することができます。
- Compensation Controls: <u>Compensation</u> 画面以外では表示されず、解析画面でも選 択対象になりません。また、<u>Assign panels</u> 画面にも表示されません。

Other Controls : FMO や isotype control、その他コントロールファイルのための分類です。
 <u>Gating Editor</u>、<u>illustration Editor</u>、<u>Sample tags</u>では表示され、使用できます。一方、
 Advanced Analysis や Export statistics では表示されず使用もできません。

FCS ファイルをアップロードする時、ファイル名に beads や comp、FMO などが含まれていると自動的に 分類されます。その他のファイルは Experiment Files に分類されます。

2-9-6-2 Experiment にある FCS ファイルが解析画面で表示されない

コントロールファイルとご認識してしまうような文字列がファイル名にある場合や、機器から出力された FCS ファイルでないファイルをアップロードした時、または、コンペンセーションコントロールファイルだがファイル名に com を含まずコントロールファイルとして認識されない時には、手動で属性を変更することができます。

① メニューバー> Sample tags> Assign panels を開きます。

3	Cytobank	Experiments	Projects			
	PBMC Experim	nent (Fluorescence) Sample tags -	-A⊷ Data QC	🛱 Gates	Advanced analyses	L. Illustrations
	Experim	Assign panels (3)		Assign files to panels and adjust	etalls	
		III Plate Annotator		t (Fluorescence)		
5		🗲 Edit Condition ta	gs (4)	:(Fluorescence)		
2			nt dimensions			
2						
22	Notes					

- ② Assign panels 画面上部の Configure control files をクリックします。
- ③ Configure control files ウィンドウで、各ファイルがどの属性かをラジオボタンで設定します。属性 名のボタンを押すと全てのファイルをその属性にすることができます。
- ④ Apply changes ボタンを押します。

Assign panels 🛈	Configure contro	Ifiles				
Configure control files	Q Filter displ	ayed				
Files mark	ed as Other Controls are exch	uded from advanced analy.	ies.			
	Filename	Tube / sample name	Experimental	Other Control	Compensation Control	
	comp_alexa 488.fcs	Alexa 488	0	0	۲	
	comp_alexa 647.fcs	Alexa 647	0	0	۲	
	pbmc_lrs005_il10.fcs	IL10	۲	0	0	
	pbmc_lrs005_il6.fcs	IL6	۲	0	0	
	pbmc_lrs005_lps.fcs	LPS	۲	0	0	
	comp_pe.fcs	PE	0	0	۲	
	comp_pe_cy7.fcs	PE_Cy7	0	0	۲	
	comp_pacblu.fcs	PacBlu	0	0	۲	
	comp_percp_cy55.fcs	PerCP_Cy55	0	0	۲	
p	bmc_lrs005_unstim1.fcs	Unstim1	۲	0	0	
р	bmc_lrs005_unstim2.fcs	Unstim2	۲	0	0	

※コンペンセーションコントロールファイルを含む<u>データ QC: PeacoQC</u>の実行後は、オリジナルのコンペン セーションコントロールファイルは Other Control になり、QC 後のファイルのみが Compensation Control になります。これはこの後自動コンペンセーションを行う場合に混乱を避けるためです。

2-1 0 Figure 作成 (Illustrations; Event level plot)

Cytobank ではひとつのデータからのグラフを作成して並べるのではなく、Experiment 内のデータセットを サンプルタグを含む項目(dimension)でレイアウト表示する仕様になっています。

Illustration Editor では図や統計表を作成することができます。操作による変更は、デフォルトではすべて自動保存されるようになっています。illustrationを別名で保存することで Experiment 内に複数の illustration を作成し保存することができます。

メニューバーの Illustrations で画面を表示します。

Cytobank	Experiments	Projects				Help 🚺 Cyto	bankSupport
PBMC Experime	ent (Fluorescence)						Sharing
Actions	Y Sample tags	1 Gates	க் Advanced analyses	L Illustrations	Scales	A Compensation	Private

2-10-1 新規 illustrator の作成

- ① X=1-N-> Illustrations> New illustration をクリックします。
- ② Illustration の名前を入力します。
- ③ Illustration Editor 画面に切り替わります。
- ④ Plots、Layout で illustration を作成し、Save や Export で保存設定や出力します。

PBMC (F	Fluorescence) ns 🏾 🌱 Sam	ple tags	🗘 Gates	🚓 Advan	ced analyses	L Illustration	ns	🐨 Scales	A Compensati	on i	Shar Pri	ing ivate
() II	ustration Edi	tor — Ne	w illustratio	n //								
	🏄 Plots 🚽	₽ *।	Layout 👻	💾 Save 👻	Expo	rt -		l	Templates -	8		8

2-10-2 Plots 設定

- ① Illustration Editor で Plots をクリックします。
- ② Plot type でプロットの種類を選択します。ここでは Event level plot について説明します。

2 - 1 0 - 2 - 1 Dot Plot : ドットプロット



- Overlaid:オーバーレイ表示のオン/オフします。
- Concatenate: <u>ファイル統合プロット</u>の有無/設定をします。
- Color by: プロットのカラー表示を以下から選択して設定します。
 - ・ Black=黒色単色
 - Density = 細胞密度でカラー表示
 - Z-axis channel=Z-axis に設定したチャネル強度のカラー表示
 - ・ Overlaid dimension=オーバーレイカラー表示(自動的にオーバーレイもオンになります)
- Palette:カラー表示のパターンを設定します。
- Show gates : Gates ツールでのゲーティング表示をオン/オフします。 Overlaid や
 <u>Concatenate</u> で <u>Tailored par file ゲーティング</u>をしている場合は、最初のファイルのゲート位置 が表示されます(統計値等は Tailored の位置で計算されます)。
- Size:プロットサイズを選択します。
- <u>More plot settings</u>: プロットの詳細設定ができます。

2 - 1 0 - 2 - 2 Histogram : ヒストグラム



以下の項目を設定できます。

- Overlaid:オーバーレイ表示のオン/オフします。
- Concatenate: <u>ファイル統合プロット</u>の有無/設定をします。
- Fill color: プロットのカラー表示を以下から選択して設定します。
 - ・ None:塗りつぶしなし
 - ・ Fixed: 左のカラーウィンドウで塗りつぶし色を指定します。



※線の色を指定したい場合は More plot settings の Outline color で可能です。

 (Overlaid 使用時のみ)Statistics:各ヒストグラムの統計値で色付けします。Fill opacity%:透明度、Statistic:色付けに使用する統計(統計によってその下の項目が変 わります)、Scale range:最大値・最小値の範囲(global=全体、Local=Table ごと、 Custom=最小値・最大値を任意に設定)を設定できます。



 (Overlaid 使用時のみ) Overlaid dimension:凡例ごとに色を変えます。Fill opacity%:透明度を設定できます。プロットの凡例カラーボックスをクリックして好きな色に変更 することもできます。



- (Overlaid 不使用時のみ)Show gates: ゲートを表示します。<u>Concatenate</u>で
 <u>Tailored par file ゲーティング</u>をしている場合は、最初のファイルのゲート位置が表示されます
 (統計値等は Tailored の位置で計算されます)。
- Size:プロットサイズを選択します。
- <u>More plot settings</u>: プロットの詳細設定ができます。

2-10-2-3Contour:等高線



- Concatenate: <u>ファイル統合プロット</u>の有無/設定をします。
- Color by: プロットのカラー表示を以下から選択して設定します。
 - ・ None:黒色の等高線
 - ・ Density:細胞密度についてカラー等高線で表示します。Palette でカラーパターン を選択できます。
 - ・ Z-axis channel: Z-axis に設定したチャネル強度のカラー表示
- Show gates : Gates ツールでのゲーティング表示をオン/オフします。 Overlaid や
 <u>Concatenate</u> で <u>Tailored par file ゲーティング</u>をしている場合は、最初のファイルのゲート位置 が表示されます(統計値等は Tailored の位置で計算されます)。
- Size:プロットサイズを選択します。
- <u>More plot settings</u>: プロットの詳細設定ができます。等高線のスムージングや%、低密度ドット 表示の設定もできます。



2 - 1 0 - 2 - 4 More plot settings

左側でプロットの詳細な設定ができます。上部は選択した設定で細かく異なりますため、お気軽に<u>サポートまでお問い合わせ</u>ください。共通する General セクションより下の項目について説明します。

utliers start at 10	•	Preview plot	
eneral			
lot size Small		▶ -	
now gates			
lack background			
xis markings		- <u>2000</u>	
xis tickmarks			
cale numbers			
how axis labels 🛛 🔍 🗌	()	- FSC	
Axis labels			
atistics		FSC-A 👻	
tatistic Median			
how tables of statistics 🛛 🗸 🔍		File / sample IL10 - pbmc_lrs005_il10 -	
ighlight control cells	~	Population Ungated -	

変更は Apply changes ボタンで反映されます。

- General:以下の設定をすることができます。
 - ・ Plot size : プロットのサイズ
 - ・ Show gates : ゲート表示
 - ・ Black background : プロットの背景を黒色表示
- Axis markings:以下の表示非表示を選択できます。
 - ・ Axis tick marks : 軸目盛り
 - ・ Scale numbers : 軸の数値
 - ・ Show axis labels : 軸ラベル (X/Y)
 - ・ Axis labels: 開くと Row/Column 毎に軸ラベルを表示するかを設定できます。
- Statistics : 以下を設定できます。
 - ・ Statistic:使用する統計値を選択できます。
 - Event count : 細胞数
 - Percentile: パーセンタイル値
 - Percent of population : %ポピュレーション
 - Median : 中央値
 - Mean: 平均値
 - Geometric mean:幾何平均
 - Standard deviation:標準偏差
 - Variance:分散
 - Coefficient of variation: 変動係数

Minimum:最小値 Maximum:最大値 Pearson correlation (X vs Y) :相関係数 Percent in gate:ゲート内% (Gate を選択します) Percentile distance(5to95) :パーセンタイル 5 と 95 の間の値 Channel range:チャネル範囲

- ・ Show tables of statistics : 画面下部の統計表の表示非表示
- ・ Highlight control cells:統計表のコントロールセルの色の有無

2-10-3 Layout 設定

- ① Illustration Editor の Layout をクリックします。
- Rows と Columns に割り当てるカテゴリー(例: FCS ファイル、populations など)をクリックして選択します。サンプルタグの dimension も割り当てることができます。
- ③ カテゴリーの右の数字 of 数字をクリックし、レイアウト表示する項目を選択します。



- ④ ③の右にオーバーカーソルで現れる網掛けをドラッグ&ドロップすると割り当てを上下に移動する ことができます。
- ⑤ (必要な場合)作成した Rows, Columns をカテゴリーでさらに縦または横に並べたい場合は、Add dimension をクリックして Table を作成します。
- ⑥ Tableの右の両矢印をクリックして並べ方を縦・横と変更できます。Concatenateを設定している場合は、矢印マークを統合マークに変更することでファイル統合プロット表示にすることができます。
- ⑦ Table も Rows, Columns と同様の変更を行うことができます。また、網掛けの右のXボタンで、削除することもできます。
- ⑧ Layout Preview ボタンを押すと並び方のレイアウトが表示されます。

※Columns, Rows, Tables のいずれかに Channels を設定した場合、X,Y,Z と選択できるメニューが表示され、各チャネルを X 軸、Y 軸、Z 軸(色強度)に設定できます。

	Colum	ns ↔	Table 1	\leftrightarrow					
	🕍 Ch	iannels	Ŷ	Populations ~					
	D pp3i D pSta D CD3	B it3 33 3	CD3- CD3- CD3- CD3- CD3- CD3- CD3- CD3-	CD33+:DC etc CD33+:Monocyte H:Tcell	Add	d colun	nn dimension		
	4	of 18 selected 🗈	3	of 8 selected 🖻					
FCS Files	Previ	ew Q						← Table 1: P	opulations
Constimation Constitution Const	Previ	ew Q	Population: Cl	D3-CD33+:DC et	tc		р	← Table 1: P opulation: CD3	opulations - -CD33+:Mono
Unstim1 LPS IL0 L205 Selected d	Previe	ew Q	Population: CI Columns: X C	D3-CD33+:DC et	tc		P	← Table 1: P opulation: CD3 Columns: X C	opulations - -CD33+:Mono hannels
Z FCS Files Unstim1 LPS IL10 IL0 2 of 5 selected ₢	Previo	ew Q X Channel: pp38	Population: CI Columns: X C XChannel: pStat3	D3-CD33+:DC en Channels X Channel: CD33	X Channel: CD3		P X Channet: pp38	← Table 1: P opulation: CD3 Columns: X C X Channet: pStat3	opulations - -CD33+:Mono hannels X Channel: CD33
D FCS Files Unstim1 UPS IIL10 UB 2 of 5 selected @ ③ Add row dimension	- Previe	ew Q XChannel: pp38 File: Unstim1	Population: CI Columns: X C X Channel: pStat3 File: Unstim1	D3-CD33+:DC et Channels X Channel: CD33 File: Unstim1	X Channel: CD3 File: Unstim1	es	P. X Channel: pp38 File: Unstim1	← Table 1: P opulation: CD3 Columns: X C X Channet: pStat3 File: Unstim1	opulations - -CD33+:Mono hannels X Channel: CD33 File: Unstim1
COS Files	- Previe	X Channel: pp38 File: Unstim1 X Channel: pp38	Population: CI Columns: X C X Channel: pStat3 File: Unstim1 X Channel: pStat3	D3-CD33+:DC et channels X Channel: CD33 File: Unstim1 X Channel: CD33	X Channel: CD3 File: Unstim1 X Channel: CD3	ows: Files	P X Channel: pp38 File: Unstim1 X Channel: pp38	← Table 1: P opulation: CD3 Columns: X C X Channet pStat3 File: Unstim1 X Channet pStat3	opulations - -CD33+:Mono hannels X Channel: CD33 File: Unstim1 X Channel: CD33

2-10-4 ファイル統合プロットの表示: Virtual concatenation

視覚的に理解しやすい figure を示すために、いくつかの FCS ファイルを統合したプロットを表示する必要 があることがあります。例えば健常者と疾患者、治療前と後、寛解者と応答なし者、などのグループ間のポ ピュレーション分布の違いを強調したい場合などです。

2-10-4-1バーチャルファイル統合(Virtual concatenation)とは

複数の FCS ファイルを1つの FCS ファイルに統合する方法は R や Kaluza など他の解析ソフトで提供されていますが、Cytobank ではファイル自体を統合することなくソフトウエア上でバーチャルに統合しプロット 表示をすることができます。実際に統合ファイルを作成するのでないため、色々な観点のグループ間比較を するための統合プロットを簡易に作成することができます。

2-10-4-2統合するためのグループの設定

Sample tags 機能で FCS ファイルにグループを割り当てます。グループの分類(dimension)はいく つも作成できるので、治療前後、治療種類、治療後のタイムポイント、検体背景(性別や年代、リスク スコア等)など様々な観点でのグループ分けができます。

2-10-4-3基本操作のガイドビデオについて

Cytobank サーバーの日本語サポート Experiment の子 <u>Experiment : v10.3 日本語ガイドビデオ</u>を開き、Experiment summary ページの Attachments セクションからダウンロードできます。

≡ A	ctions	🌱 Sample tags	-∕∿- Data QC	🗘 Gates	क्तै Advanced a	nalyses	L Illustrations	10 ¹ Scales	A Compen	isatio
?	Experime	nt summary	Clone	🥒 Edit	details					
c	🤊 Linked	experiments								
					_	⑦日本語 Cytobank Japa	サポート support for nese			
					€ v (10.3日本副 virtual con	語ガイドビデオ ncatenation)			
	Total	size of all linked: 5	9.9 MB							
1	Illustra This experi	ntions	New any illustrations yet.							
	🕼 Crea	te an illustration								
6	🖉 Attach	ments								
	File Na	ne			Download	Date	Uploaded By	Size	md5sum	
	Cytoba No desc	nk_v10.3ガイドビラ ription	デオ_MAPSS-LS-20	02308-43.mp4		Aug 27	🙎 Yuko Nakane	41.3 MB	75ae3e2	ŵ

2-10-4-4基本的な操作方法

バーチャルファイル統合(VC)は illustration 機能内にあります。 Dot plot、 histogram、 contour で使用することができます。

- ① Illustration Editor を開きます。
- ② Plots 設定を開きます。
- ③ Dot plot、histogram、contour のいずれかを選択します。
- ④ Overlaid がオフになっていることを確認します。Concatenate と Overlaid は併用できません。
- ⑤ Concatenate で以下を選択します。
 - All multi-file positions: Layout に割り当てられた dimension や sample tags で複数ファイルがあるものをすべて統合します。
 - Rows または Columns: Layout の Rows または Columns に割り当てられているものだけ を統合します。

Actions	🌱 Sample tags	-∿- Data QC	🏠 Gates	ភ្លំ Advanc	ed analyses	L Illustrations
🕕 Illustra	tion Editor — C	ytobank Suppor	rt illustration			
<u>iá</u>	Plots 🛑 🛒	Layout - [Save 🚽	Export		
(Plot type	🚈 Dot 🗲	ď			
	Overlaid				CI	045
	Concatenate	All multi-file positions	s		600	
	Color by	Density				Lympn
	Palette	En Fuego				
	Show gates	 ✓ ○ 				
	Gate label	Gate Name		1_patient1_BCR-XL	angt]	
	Size	/ledium	ı_pati			CD45+
	🗹 More	e plot settings			-[
					C	:D45
					ft i	
		PBI	MC8_30min_pati	ient2_BCR-XL		
						CD45+
					=	
					C	:D45

- Layout 設定を開きます。
- ⑦ Rows, Columns, Tables に各々dimension や sample tags を割り当てます。
- ⑧ 統合または表示対象とするか対象外とするかを x of y をクリックで選択することができます。



⑨ どのファイルが統合されたプロットかを確認するにはプロットオーバーカーソルで表示されるウィンドウ上部
 に示されます。



⑩ VCを解除/設定したい場合は、レイアウトの統合(Concatenated)マークを縦矢印または横矢 印にします。

₽ Layout	💾 Save 👻	D Evport	
		It Export	
(i)			ď
Rows	1 🗹 FCS Files	8 of 16	
Columns	Orientation of R	lows	CD45
Table 1	Vertical		CD3+Lymph
Tuble 1	Concatenated	•	
	Rows Columns Table 1	Rows Image: Columns Columns Orientation of R Table 1 Image: Concatenated	Rows Image: Columns Drientation of Rows Columns Orientation of Rows Table 1 Image: Concatenated

⑪ (Plots 設定の Concatenate が All multi-file positions の場合) プロットのオーバーカーソルで 表示されるウィンドウの Concatenate を設定することでも統合プロットにすることができます。

နို Advanced analyses	L Illustrations	There are multiple (8) files assigned to this layout position: PBMC8_30min_patient1_Reference, PBMC8_30min_patient2_Reference, PBMC8_30min_patient3_Reference, PBMC8_30min_patient5_Reference, PBMC8_30min_patient5_Reference, PBMC8_30min_patient5_Reference,
	Ungated	PBMC8_30min_patient8_Reference
		Layout position Row – Condition: Reference Column – Population: Ungated X axis: pStat5 (pStat5(Nd150) * Y axis: Cell_length *
		Plot settings I More plot settings Plot type
		Overlaid Concatenate None Color by Density Palette En Fuego Show gates Size Medium

※ヒストグラムや等高線プロットでも同様に設定できます。

※マニュアルゲーティング解析だけでなく、次元削減マップでも同様に設定できます。

2-10-4-5簡単なチェック方法

表示されているプロットの統合を確認するには、プロット下の Statistics で Displaying を Raw、 event count に設定すると細胞数が表示されるので簡単に確認できます。

) Hide		
	Displaying 🛈 Raw Ever	nt count 🗸 💽
	F	CS Files
	single	33553
	CD3+	23745

2-10-4-6次元削減マップのファイル統合プロット表示

不均一性の高いデータセットの場合、次元削減マップの分布がファイルで偏りデータセット全体を把握し難いことがままあります。そのような時にファイル統合プロット(VC)にすると、全体と各ファイルや各グループの分布を確認することができます。

※VCの他に、Rを使用して FCS ファイルを1つにまとめた統合ファイルを作成し、それを Cytobank へ 再アップロードする方法のご案内が必要な場合はサポートまでお気軽にお問い合わせください。

- ① 次元削減(tSNE または UMAP)解析結果 Experiment を開きます。
- ② グループごとに統合したい場合は Sample tags でグループの tags を作成しておきます。
- メニューバーillustrations > New illustration をクリックし、新規 illustration を作成します。
- ④ DR マップを表示します。
- ⑤ Plots 設定で Concatenate を統合したいレイアウト(Rows または Columns)に設定するか、Layout 設定で統合した項目の矢印マークを統合マーク(Concatenated)に設定します。



 ⑥ または、Plots 設定で Concatenate を All multi-file positions に設定すると、レイアウトで 割り当てた dimension や sample tags で複数ファイルのものを統合して表示してくれます。



⑦ 統合プロット上にマウスオーバーするとウィンドウが現れ、どのファイルの細胞が統合されたプロットかを確認することができます。



2-10-5 Save 設定

Illustration EditorのSave 設定では、以下を設定できます。

	Plots -	,∰* Layout ⊸	💾 Save 👻	Export	
			🗹 Auto-sav	/e enabled	
			🗞 Clone ne	ew copy	
		CD3-CD33+:I	📰 Save set	tings template	(j
	pp38	pStat3	🗊 Reset		
ſ			💼 Delete		

- Auto-save enabled : illustration の自動上書き保存のオン/オフ。オフにすると Save ボ タンが表示され、押した時にだけ上書き保存されるようになります。
- ② Clone new copy: 現在の illustration の複製を作成します。
- ③ Save settings template : Plots 設定をテンプレートとして名前をつけて保存します。保存 したテンプレートは右側の Templates ボタンから呼び出すことができます。
- ④ Reset: illustration を初期状態に戻します。
- ⑤ Delete: illustration を削除します。

2-10-6 Export 設定

Illustration EditorのExport設定では以下を設定し、Exportボタンを押すことでFigureとしてエクスポートできます。



- Illustration の名前部分をクリックしてエクスポートファイル名の変更、右のプルダウンメニューで ファイル形式を SVG、PDF、PNG から選べます。
- ② Export options を開くと以下の項目をエクスポートファイルに含めるかを設定できます。
 - Save a copy as experiment attachment: エクスポートファイルを experiment
 の attachment セクションに自動添付保存します(デフォルトオン)。
 - Include illustration name header : illustration 名をヘッダーにします(デフォルトオン)。

- Include experiment details : experiment の details 欄に記載した内容を含め ます(デフォルトオン)。
- Include tables of statistics : illustrations 画面下部の統計表を一緒にエクスポ ートします(デフォルトオフ)。
- ・ Include QR code link: URLのQRコードを含めます(デフォルトオフ)。

2-10-7 Templates ボタン

Illustrations のテンプレートを呼び出して適用することができます。

	ີ ມີ Scales	🖄 Compensat	ion 🚦	Sharing I <mark>Private</mark>
	E	🗄 Templates 👻	-	8
Personal viSNE Dura IM Basic heatmap Public Pairwise Plots Dot Overlays Channel-colored Do Cytobank Classic D Cytobank Classic H Traditional Histogra	ot ensity Dot istogram O ım Overlay:	verlays		
				× 🖓

- Personal:ご自身で保存したテンプレートやアクセス共有されたテンプレートがリスト表示 されます。削除したい場合は右のゴミ箱ボタンを押します。また、他のユーザーと共有したい 場合は、その隣にある共有ボタンを押し、ユーザーネームを入力、選択します。
- ・ Public:公開されている experiment のテンプレートを使用することができます。

2-10-8 印刷ボタン

印刷ビューをブラウザの別タブで開くことができます。

ີ Scales	⚠ Compensation	Sh	aring Private
5 5 5	🗄 Templates 👻 🌔		8

2-10-9 ロックボタン

デフォルトではロックがかかっており、Experimentを他のユーザーとフルアクセス共有している場合、他の ユーザーが illustration を編集できないようになっています。 他のユーザーと編集を共有したい場合は、ロックボタンをクリックして解錠します。各ユーザーからの変更は 上書きになります。



2-10-10 プロット上でのレイアウト変更操作

illustrations 上のプロットやその周辺(軸ラベルや Columns, Rows, Tables の項目名)をクリック やドラッグ&ドロップすることで、使用する項目を変更したり、順番を変更したりすることができます。

2-10-11 Statistic table : プロット下の統計表

プロット下部にプロットと同じレイアウトの統計表が表示され、定量的な検討をすることができます。表の表示/非表示は Plots 設定の more plot settings で設定できます。

表上部の Displaying セクションでは統計表と上部プロットで扱う統計値を以下から選んだり、統計対象を変更したりすることができます。

- Raw:生の値
- Log10 ratio:Log10比
- Fold:フォールド
- Log2 ratio:Log2比
- Arcsinh ratio: Arcsinh 変換後の比
- Transformed ratio:スケール設定で設定されている変換(リニア/Log/Arcsinh)後の比
- Difference: 差分

※統計グラフで複数のファイルの統計値を一つのセルの値で表示する場合は、Raw 以外は選択できな くなります。



CSV ボタンを押すと、CSV ファイルとしてダウンロードし、エクセル等で数値	፤を扱うことができます。
---	--------------

A1		* 1 2	$\sim J_X$	Condit	ion														
4	- A	в	с	D	E	F	G	н	1.1	 к	L.	м	N	0	P	Q	R	s	
1	Condition	X channel	Population	FCS Filenan	Median														
2	Unstim	CD33	Ungated	Unstim1	89.68313														
3	Unstim	CD33	8 cells (Sca	Unstim1	46.24519														
4	Unstim	pStat3	Ungated	Unstim1	38.89934														
5	Unstim	pStat3	B cells (Sca	Unstim1	28.63679														
6	1.6	CD33	Ungated	11.6	89.65573														
7	11.6	CD33	8 cells (Sca	11.6	44.96155														
8	11.6	pStat3	Ungated	IL6	219.0314														
9	11.6	pStat3	B cells (Sca	IL6	32.46771														
10																			
11																			
12																			

2-11統計グラフと統計的有意差検定(Illustrations; summary plot)

統計グラフとしては、Heatmap、Box plot、Violin plot、Bar chart、Line chartとSummary dot があります。



Sample tags 機能と組み合わせることで多様な作図ができます。Sample tags 機能では実験上の 個体番号や条件、サンプルの種類などの変数を"dimension"として扱い、各 FCS ファイルにアノテーショ ンとして割り当てることができます。統計プロットや検定をする場合はこの Sample tags との組み合わせ が特に重要で、例えばサンプルグループの dimension を作成し、健常者と疾患患者や治療前と後、治 療効果のありとなし、など各グループの tags を FCS ファイルにアノテーションすれば、グループ間の違いを 示すためのボックスプロットやバイオリンプロットを表示したり、有意差検定の結果を p 値やシンボル(*) でプロット内に表示したりすることもできます。

2-11-1統計グラフ設定の基本

統計チャートは、イベントレベルプロットとは対照的に、ひとつのボックスやバイオリン、バー等にサンプルグル ープの情報をまとめて表示するようデザインされています。多くの場合、ボックスやバイオリン、バーごとにある 実験変数を反映する表示を目指すことでしょう。したがって、有意義な統計チャートを作成するには、 Sample tags でファイルに実験変数やそのカテゴリーをアノテーションしておくことが必須です。 サンプルグループでまとめた表示をしたくない場合は、FCS ファイルを Layout 設定で選択し、Sample tags で作成したものは選択から除外します。 例として下図に、サンプルファイルごとの統計チャート(上)と Conditions でまとめられたグループでの統 計チャート(下)を示します。



Plot 設定の Plot type でボックスプロット、バイオリンプロット、ラインチャートまたはサマリードットを選択す ると、Layout 設定の Rows と Columns は、各々x-axis groups と Subgroups に表示が変わりま す。Subgroups では、ボックスやバー、バイオリンにサンプルをどうまとめるか(グループ間比較のグループ 分け)を設定し、x-axis groups ではその比較を何を対象に X 軸に並べるかを設定します。

subgroups に Conditions を設定し、その Sample tags である unstimulated と stimulated に ついてボックスプロットを表示する設定になっています。 X-axis groups には Channels を設定し検討し たいチャネルである I F N gとT N F a を選択して表示しています。(Conditions, Channels の順 序は変更可能)。



97

最大 50 個の統計チャートを Table として並べて表示することができます。 50 以上のチャートを使用して レイアウトを構成しようとすると、「The maximum number of summary charts supported is 50. Please adjust your layout. サポートされているサマリーチャートは最大 50 です。 レイアウトを変 更してください」というエラーメッセージが出ます。

2-1 1-2Heatmap : ヒートマップ

📴 Plots 🗟	,⊞* Layout →	💾 Save 🚽
(i) Plot ty	pe 🔛 Heatmap	ď
Statis	tic Percent in gate	
Gá	te 🏠 Cells	
Scale ran	ge Global	
Pale	tte Yellow	
1: >; •	Viewthrough plot	
ľ	More plot settings	

2-11-2-1基本のヒートマップ

Plots > Plot type を Heatmap にします。

- Statistic : 使用する統計値を選択できます。
 - Event count:細胞数
 - ・ Percentile:パーセンタイル値
 - Percent of population: population A /population B ※BはAの上流ポ ピュレーションという前提。計算は厳密には(A⊆B)/Bで算出される。
 - ・ Median:中央値
 - ・ Mean:平均値
 - Geometric mean:幾何平均
 - Standard deviation : 標準偏差
 - Variance:分散
 - Coefficient of variation:変動係数
 - Minimum:最小值
 - ・ Maximum:最大値
 - Pearson correlation (X vs Y): 相関係数
 - Percent in gate:指定したゲート内イベントが何%か
 - ・ Percentile distance (5to95): パーセンタイル 5と95の間の値

- ・ Channel range: チャネル範囲
- Scale range:ヒートマップカラー範囲の設定を選択できます。
 - ・ Global: すべてのヒートマップ表での最小値と最大値で表示



Local: 各ヒートマップ表ごとの最小値・最大値で表示



・ Custom:最小値・最大値を任意に指定して表示



- Palette:ヒートマップカラーを選択できます。
- Viewthrough plot:ヒートマップをオーバーカーソルしたときにポップアップするプロットの設定
- <u>More plot settings</u>: 詳細設定

ヒートマップのカラー表示に使用する値は、生値の他にプロット下部の<u>Statistic table</u>の Displaying セクションでも変更できます。使用する統計値等を変更することで色強度表示の範囲も大きく影響を受 けるため、より適切な設定をお使いください。



2-11-2-2統計値ヒートマップ

Plots 設定で Collapse tables をオンにするか Layout 設定の Table で Collapsed を選択すると、 Layout で Collapsed に割り当てた dimension の sample tags に従ってサンプルをまとめ、 median か mean の統計値ヒートマップ表示することができます。



- Median か mean かは Plots 設定の More plot settings の Aggregate で選択できます。

_



 プロット上で Collapsed 項目をクリックし、ポップアップメニューの Header settings の Color で背 景色を変更できます。



Table1、Table2…と複数をCollapsed にすると複数層の統計値ヒートマップ表示もできます。



2-11-3Box Plot:箱ひげ図



Plots > Plot type を Box にします。

- Show dots: オンにするとグループ内の各サンプルファイルの値をドットで表示します。
- Dot size: ドットのサイズを大きくまたは小さくします。
- Show lines: 対応ありデータの場合、同一サンプルをドットで結びます。プロットの下に、 サンプルがどう対応しているかに関する説明が表示されます。
- Gray shading: ドットはカラー表示のまま、ボックスをグレーで表示します。
- Jitter: ドットの並び方を変更します。0 は全てのドットを垂直方向に、1 はドットを最大 限離して表示します。
- Y axis scales: Y軸の最小・最大スケールを選択します。Local = Table ごと、
 Global=表示データセット全体、または Custom (ユーザーが値を入力) が選択可能
 です。
- Width: ボックスの幅を設定します。幅は、固定、自動拡大、カスタムの3つから選べます。
- Statistic: Y軸に使用する統計値を選択します。
 - Event count : 細胞数
 - ・ Percentile:パーセンタイル値
 - Percent of population: population A /population B ※BはAの
 上流ポピュレーションという前提。計算は厳密には(A⊆B)/Bで算出される。
 - ・ Median:中央値
 - ・ Mean:平均値
 - Geometric mean:幾何平均
 - Standard deviation : 標準偏差

- Variance:分散
- Coefficient of variation: 変動係数
- ・ Minimum : 最小値
- ・ Maximum:最大値
- Pearson correlation (X vs Y):相関係数
- Percent in gate:指定したゲート内イベントが何%か
- ・ Percentile distance (5to95): パーセンタイル 5と95の間の値
- Significance test: <u>有意差検定</u>の方法を選択します。



CD4+T 細胞に対する刺激の効果を IFNg 及び TNFa の median 値で示すボックスプロット例。ドットは各サンプルファイルの値を示す(青は未刺激、オレンジは刺激)。

2-1 1-4 Violin plot:バイオリンプロット

Plots		×
Plot type	↓ [†] • Violin	
Show dots	\checkmark	
Dot size	ó 0	
Jitter	0.3	
Show lines	\bigcirc	
Show boxes		
Gray shading	\bigcirc	
Y axis scale	Local	
Width	Fixed	
Statistic	Event count	
Generation Significance test	None	

Plots > Plot type を Violin にします。

- Show dots: オンにすると各サンプルファイルの値ドットで示します。
- Dot size: ドットのサイズを大きくまたは小さくします。
- Show lines: 対応ありデータの場合、同一サンプルをドットで結びます。プロットの下 に、サンプルがどう対応しているかに関する説明が表示されます。
- Gray shading: ドットはカラー表示のまま、バイオリンをグレーで表示します。
- Jitter: ドットの並び方を変更します。0は全てのドットを垂直方向に、1はドットを最大 限離して表示します。
- Show boxes : ボックスを表示します。
- Y axis scales: Y軸の最小・最大スケールを選択します。Local = Table ごと、
 Global=表示データセット全体、または Custom (ユーザーが値を入力) が選択可能
 です。
- Width: ボックスの幅を設定します。幅は、固定、自動拡大、カスタムの3つから選べます。
- Statistic: Y軸に使用する統計値を選択します。
 - Event count : 細胞数
 - ・ Percentile:パーセンタイル値
 - Percent of population: population A /population B ※BはAの 上流ポピュレーションという前提。計算は厳密には(A⊆B)/Bで算出される。
 - ・ Median:中央値
 - ・ Mean:平均値
 - Geometric mean:幾何平均
 - Standard deviation : 標準偏差
 - ・ Variance:分散
 - Coefficient of variation: 変動係数
 - ・ Minimum:最小値
 - ・ Maximum:最大値
 - Pearson correlation (X vs Y):相関係数
 - Percent in gate:指定したゲート内イベントが何%か
 - ・ Percentile distance (5to95): パーセンタイル 5と95の間の値

- Significance test: 有意差検定の方法を選択します。



CD4+T 細胞が刺激後に IFNg 及び TNFs の増加を示すバイオリンプロット例。ドットは各サンプルファイルの値を示す(青は未刺激、オレンジは刺激)。

2-11-5Bar chart : 棒グラフ

(i) Plots			×
Plot type	Jul Bar		
Stacked bar	\bigcirc		
Cut off	None	\odot	
Y axis scale	Local		
Width	Fixed		
Error bars	Standard deviation		
Summary method	Median		
Statistic	Event count		
(i) Significance test	None		

Plots > Plot type を Bar にします。

- Stacked bar: このオプションをオンにするとスタックバー(積み上げ)表示になります。
- Stacked method: absolute=積み上げ、relative=100%積み上げを選択できま す。
- Cut off: 入力した値に点線を表示します。

- Y axis scales: Y軸の最小・最大スケールを選択します。Local = Table ごと、
 Global=表示データセット全体、または Custom (ユーザーが値を入力) が選択可能
 です。
- Width: ボックスの幅を設定します。幅は、固定、自動拡大、カスタムの3つから選べます。
- Error bars: エラーバーを、None=表示しない、Standard deviation=標準偏差、
 Standard error=標準誤差、95th confidence interval=95%信頼区間から 選べます。
- Summary method: Statistic で選択したパラメータについて、バーの表示を複数ファ イルの平均値にするか中央値にするかを選べます。
- Statistic: Y軸に使用する統計値を選択します。
 - Event count:細胞数
 - ・ Percentile:パーセンタイル値
 - Percent of population: population A /population B ※BはAの 上流ポピュレーションという前提。計算は厳密には(A⊆B)/Bで算出される。
 - ・ Median:中央値
 - ・ Mean:平均値
 - Geometric mean:幾何平均
 - Standard deviation:標準偏差
 - Variance:分散
 - Coefficient of variation: 変動係数
 - ・ Minimum:最小値
 - Maximum:最大值
 - Pearson correlation (X vs Y):相関係数
 - Percent in gate:指定したゲート内イベントが何%か
 - ・ Percentile distance (5to95): パーセンタイル 5と95の間の値
 - Significance test : <u>有意差検定</u>の方法を選択します。



CD4+T 細胞に対する刺激の効果を、IFNgとTNFaの値で示す棒グラフ例。オレンジは刺激、青は未 刺激コントロール。

2-11-6Line chart & Stain index:線グラフとステインインデックス

(i) Plots		×
Plot type	✓ Line	
Stain index	\bigcirc	
Show dots	\checkmark	
Dot size	6 0	
Y axis scale	Local	
Width	Fixed	
Error bars	Standard deviation	
Summary method	Median	
Statistic	Event count	
(i) Significance test	None	

Plots > Plot type を Line にします。

- Stain index:ステインインデックスのオン/オフ。ステインインデックスの作成方法は、後述のステインインデックスの項をご覧ください。
- Show dots: オフにするとドットが消え線のみの表示になります。
- Dot size: ドットのサイズを大きくまたは小さくします。
- Y axis scales: Y軸の最小・最大スケールを選択します。Local = Table ごと、
 Global=表示データセット全体、または Custom (ユーザーが値を入力) が選択可 能です。
- Width: ボックスの幅を設定します。幅は、固定、自動拡大、カスタムの3つから選べます。

- Summary method : Statistic で選択したパラメータについて、ライン表示を複数フ ァイルの平均値にするか中央値にするかを選べます。
- Statistic: Y軸に使用する統計値を選択します。
 - Event count:細胞数
 - ・ Percentile:パーセンタイル値
 - Percent of population : population A /population B ※Bは A の上流ポピュレーションという前提。計算は厳密には (A⊆B)/B で算 出される。
 - ・ Median:中央値
 - ・ Mean:平均値
 - Geometric mean:幾何平均
 - Standard deviation : 標準偏差
 - Variance:分散
 - Coefficient of variation: 変動係数
 - ・ Minimum:最小値
 - ・ Maximum:最大値
 - Pearson correlation (X vs Y):相関係数
 - Percent in gate:指定したゲート内イベントが何%か
 - ・ Percentile distance (5to95): パーセンタイル 5と95の間の値
- Significance test : <u>有意差検定</u>の方法を選択します。



刺激の有無による CD4 + T 細胞での IFNg (オレンジ) と TNFa (緑)の 発現強度中央値のグループ 内複数ファイル中央値の違いを示すラインチャート例(左:エラーバーは標準偏差で表示)。レイアウト 設定は右に示すとおり。

- ステインインデックス

ステインインデックスは使用している機器において様々な蛍光色素の相対輝度を測定します。この値は蛍 光色素の輝度の比較や順位付けに使用でき、質の高いサイトメトリーデータの取得には欠くことのできな い、抗体タイトレーションによって最適濃度を決定するのに役立ちます。抗体の使用量が少なすぎると標
識が不完全になってしまいますが、使用量が多すぎると蛍光バックグラウンドの増加や、親和性の低いサイトへの抗体の結合(擬陽性)の原因となってしまう上に実験コストが高くなります。

ステインインデックスは、ポジティブポピュレーションとネガティブポピュレーションとの分離を、ネガティブポピュレ ーションの標準偏差の 2 倍で割った値と定義されます。



図:ステインインデックス:ポジティブポピュレーション(緑)とネガティブポピュレーション(黒)との 分離をネガティブポピュレーションの標準偏差の2倍で割った値と定義されます。

Cytobank ではステインインデックスチャートを使用して、タイトレーションから最適な抗体濃度を決定する ことができます。バックグラウンドノイズを最小限に抑えつつ、明るいシグナルを得るための最適な濃度や希 釈率はステインインデックスチャートを基に決定されます。ステインインデックスチャートを作成する手順は下 記の通りです。

- タイトレーションのサンプルデータを new experiment にアップロードし、Gating Editor で対象の 抗体のチャネルでポジティブとネガティブのゲートをかけます。マーカー発現は抗体濃度によって異なる ことがあるため、必ず全てのファイルを確認してゲートを調整します。プロット左上の Apply gates ボ タンをクリックしてゲートを illustrations に適用します。
- ジ サンプルファイルに Doses と Reagents の dimension での sample tag を付けます。Sample tags>Add experiment dimensions をクリックし、次のウィンドウで Tag files with Doses と Tag files with Reagents をクリックして各々設定します。現在サポートされているのはひとつの



illustration につき1 試薬です。同一の Experiment で複数の試薬を検討する場合は、試薬ご とに新規の illustration を作成し、ステインインデックスチャートを作成ください。

			e		Dose Actions
nter a comma separated list of Doses b	o add:				Dases
Dose 1, Dose 2, Dose 3, etc.		Add Doses			2X
					12
Doses 2X 1X 1.2	14 18 116 132 0				24
All Dotes					118
g files with Doses by dragging them to the	box for the desired tag. You can also filter the	les displayed and move them in bulk using "Move	effies to		1:32
Untagged	2X Tagged Files 🛞	TX Tagged Files 🛛 🛞	12 Tagged Files 🛞	1:4 Tagged Files 🛞	unstained
Fiter	05-Well-A3.4cs	01-Well-A2/cs	01-Well-A3.hts	01-Well-A4.fcs	Import Doses from another experiment
(MAR 10 10					Duplicate a File
					01-We8-A1.5cs -
					Duplicate
					D Reset Doses
	18 Terred Files	116 Tarred Files	122 Tarred files	0 Tarrend Files	
	0530645557	(Manual Article	01/00LATex	OT AMPLAN for	
	11 110 1000	(a) memory	(41 mile 14 14)	011101000	
Reagents			0		Researt Actions
Reagents	rom the following list to	create tage (names can	Accion	nanale) —	Reagent Actions
Reagents ect Reagent name(s) fr	rom the following list to	create tags (names can	edited using Assign	panels) —	Reagent Actions Reagents
Reagents ect Reagent name(s) fr lect Reagent name(s)	rom the following list to	create tags (names can	be edited using Assign	panels) —	Reagent Actions Reagents CD4-PacBlue V405-450BP-A
Reagents ect Reagent name(s) fr lect Reagent name(s)	rom the following list to	create tags (names can hts	De edited using Assign	panels) —	Reagent Actions Reagents CD4-PacBlue V405-450BP-A
Reagents ect Reagent name(s) fr liect Reagent name(s)	rom the following list to Add Reager	create tags (names can Its	be edited using Assign	panels) —	Reagent Actions Reagents CD4-PacBlue V405-450BP-A Duplicate a File
teagents ct Reagent name(s) fr lect Reagent name(s) covents	Add Reager	create tags (names can Its	De edited using Assign	panels) —	Reagent Actions Reagents CD4-PacBlue V405-450BP-A Duplicate a File 01-Well-A1.fcs
teagents ct Reagent name(s) fr lect Reagent name(s) eagents CD4-PacB	Add Reager	create tags (names can Its	De edited using Assign	panels) —	Reagent Actions Reagents CD4-PacBlue V405-450BP-A Duplicate a File 01-Wel-A1.fcs Dumlisete
teagents ct Reagent name(s) fr lect Reagent name(s) eagents CD4-PacB All Reagents	Add Reager	create tags (names can Its	be edited using Assign	panels) —	Reagent Actions Reagents CD4-PacBlue V405-450BP-A Duplicate a File 01-Well-A1.fcs Duplicate
eagents ct Reagent name(s) fr lect Reagent name(s) eagents CD4-PacB All Reagents les with Reagents by drag	rom the following list to Add Reager like V405-4508P-A aging them to the box for the	ecreate tags (names can tts	be edited using Assign the files displayed and move	panels) — : them in bulk using 'Move files to	Reagent Actions Reagents CD4-PacBlue V405-450BP-A Duplicate a File 01-Well-A1.5cs Duplicate
eagents ct Reagent name(s) fr lect Reagent name(s) eagents CD4-PacB All Reagents les with Reagents by drag	Add Reager	desired tag. You can also filter	be edited using Assign the files displayed and move	panels) — = them in bulk using 'Move files to	Reagent Actions Reagents CD4-9x68lue V405-4508P-A Duplicate a File U1-Wel-A1.5c Upplicate
Reagents cct Reagent name(s) fr lect Reagent name(s) ccdt Reagent name(s) CD4-PacB All Reagents Mies with Reagents by drag Untagged	Add Reager Add Reager Hue V405-4508P-A aging them to the box for the CD 4-Pa	desired tag. You can also fifter cellue V405-4508P-A ®	be edited using Assign the files displayed and move	panels) — them in bulk using "Move files to	Reagent Actions Reagents CD4-PacBlue V405-4508P-A Duplicate a File 01-Well-A1.5cs Duplicate
Reagents cct Reagent name(s) fr lect Reagent name(s) reagents CD4-PacB All Reagents Untagged Untagged Filter	Add Reager	desired tags (names can desired tag. You can also fifter CBLue V405-4508P-A Files	be edited using Assign the files displayed and move	panels) — them in bulk using 'Move files to	Reagent Actions Reagents CO-Pardiau V405-4508P-A Drughiceta # File 01-Well-A1.fcs Duplicate Duplicate
Reagents cct Reagent name(s) fit lect Reagent name(s) lect Reagent name(s) cD4-PacB cagents CD4-PacB cagents lites with Reagents by drag Untagged Timer Move files to	Add Reager	desired tag. You can also fifter CBLue V405-450BP-A ® Files	be edited using Assign the files displayed and move	panels) — = them in bulk using 'Move files to	Reagent Actions Reagents CD4-PacBlue V4054508P-A Duplicate a File 01-Well-A1.5cs Duplicate Duplicate
teagents ct Reagent name(s) fr iect Reagent name(s) cD4-PacB cD4-PacB all Reagents ies with Reagents by drag Jatagged itter Move files to	Add Reager Add Reager Add Reager	create tags (names can ita desired tag. You can also fifter cBlue V405-4508P-A Files	be edited using Assign the files displayed and move	panels) —	Reagent Actions Reagents CO+ParaBlue V405-4508P-A Duplicate a File Di-Veel-A1.5cs Duplicate Duplicate
teagents ct Reagent name(s) fit lect Reagent name(s) fit lect Reagent name(s) cD4-PacB cD4-Pa	Add Reager Add Reager	create tags (names can ts desired tag. You can also filter citible V405-1500P-A @ Files HA1fcs HA2fcs	be edited using Assign the files displayed and move	panels) —	Reagent Actions Reagents CD4-PacBlue V405-4508P-A Duplicate File Duplicate Duplicate
Reagents ct Reagent name(s) fi fiect Reagent name(s) fi fiect Reagent name(s) CD4-PacB All Reagents fier with Reagents by drag Untage Untage Untage	Add Reager Add Reager Add Reager	create tags (names can its desired tag. You can also fitter cibbe V405-4500 P-A Fittes II-A1fcs II-A2fcs 	be edited using Assign the files displayed and move	panels) —	Reagent Actions Reagents CO4-Pacillus V405-4508P-A Duplicate a File 01-Well-A1.5cs Duplicate Duplicate
Reagents ct Reagent name(s) fi dect Reagent name(s) dect Reagent name(s) CD4-PacB All Reagents Uses with Reagents Uses with Reagents Wave files to	Add Reager Add Reager Add Reager	create tags (names can tag) desired tag You can also fitter cible V405 can also fitter fittes IFA1fcs IFA3fcs IFA3fcs	be edited using Assign the free slipplayed and move	panels) —	Reagent Actions Reagents CD4-PadBur V405-4508P-A Douglicates File 01-Well-A1.5cs Duplicate Duplicate
Reagents Ct Reagent name(s) fit lead Reagent name(s) lead Reagent name(s) CD4-PacB AII Reagents View with Reagents by drag Jintagged View Tee Tee to	Add Reager Add Reager Add Reager Add Reager	create tags (names can its desired tag. You can also filter Cellule V406-4500P-A RA155 RA155 RA155 RA155	be edited using Assign the fries displayed and move	panels) — them in bulk using 'Move files to	Reagent Actions Reagents CO4-Pacillar V405-4508P-A Duplicate a File 01-Well-A1.5cs Duplicate Duplicate
Reagents tct Reagent name(s) fr eagents CD4-PacB All Reagents Untagged Untagged Wove files to	Add Reager Add Reager Add Reager to V405-4508P.A ging them to the box for the CD4-P CD4	create tags (names can tag) desired tag You can also fitter cititue V405-4500P-A @ HA1/cs HA1/cs HA1/cs HA1/cs HA1/cs	be edited using Assign the files displayed and move	panels) —	Reagent Actions Reagents CO+ParaBlue V405-450BP-A Duplicate a File Di-Veel-A1.5cs Duplicate Duplicate
Reagents text Reagent name(s) fn text Reagent name(s) text Reagent name(s) text Reagents text Reagents text Reagents text with Reagents text with Reagents text reagent text R	rom the following list to Add Reager the V405 4508P.A ging them to the loss for the Tagging them to the loss for the O1-WE O1-WE	create tags (names can test as to be as a second of the se	be edited using Assign the files displayed and move	panels) —	Reagert Actions Reagerts C04-PastlacV405-4508P-A Duplicate a File 01-Well-A1.fcs Duplicate Duplicate

Tag files with Dosesの画面(上)とTag files with Reagentsの画面(下)

- ③ Sample tags 機能で、希釈情報 Doses をファイルにタグ付けします。抗体を 2X(推奨の抗体 濃度の 2 倍)から 1:32 に希釈した例を下に示します。
- ④ ファイルをタグ付けしたら画面上部の水色のナビゲーションバーから Illustrations をクリックし、新規 illustration を作成します。Illustration Editor で以下のように Plots と Layout を設定しま す。Plot type として Line チャートを選択し、Stain index をオンにします。X-axis group に Doses、Table 1 に Reagents を設定します。希釈順序(並び順)は Doses の右側(x of y (x と y は数字))をクリックして変更できます。

🖲 Illustration Editor — Stain Index 🥒	() Illustration Editor -	– Stain Index 🧷			
🕍 Plots 🆽 Layout	🕌 Plots 🚽	⊞* Layout ⊸	💾 Save 🚽	Export	,
③ Plot type ▲ Line		(i) X-axis groups	🖉 Doses	8 of 8	ď
Stain index 🔽					
Show dots		Table 1 1	Reagents	1 of 1	
Dot size 6				10.1	
Y axis scale Local			Add dimension	n	
Width Fixed			☑ [®] Layout Preview		
l					

ステインインデックスチャートの Plots 設定(左)と Layout 設定(右)

⑤ ステインインデックスチャートは、Illustration Editor で直接保存やコピーができます。Reagent で 指定した試薬の各 Dose のステインインデックス値は、チャートグラフ下の Stain Index Statistics に表示されます。



2-1 1-7 Summary dots : サマリードット

(i) Plots		×
Plot type	Summary dot	
Dot size	6	0
Swarm	\bigcirc	
Jitter	0.3	\odot
Mean/median line	None	
Y axis scale	Local	
Width	Fixed	
Statistic	Event count	
(i) Significance test	None	

Plots > Plot type を Summary dot にします。

以下の項目を設定できます。

- Dot size: ドットのサイズを大きくまたは小さくします。
- Swarm:オンでなるべく一列に並べます。
- Jitter: ドットの並び方を変更します。0は全てのドットを垂直方向に、1はドットを最大 限離して表示します。
- Mean/median line:平均または中央値を示すラインを表示します。

- Y axis scales: Y 軸の最小・最大スケールを選択します。 Local = Table ごと、 Global = 表示データセット全体、または Custom(ユーザーが値を入力)が選択可能 です。
- Width: ボックスの幅を設定します。幅は、固定、自動拡大、カスタムの3つから選べま す。
- Summary method: Statistic で選択したパラメータについて、ライン表示を複数ファ イルの平均値にするか中央値にするかを選べます。
- Statistic: Y軸に使用する統計値を選択します。
 - Event count:細胞数
 - ・ Percentile:パーセンタイル値
 - Percent of population : population A /population B ※BはAの 上流ポピュレーションという前提。計算は厳密には (A⊆B)/B で算出される。
 - ・ Median:中央値
 - ・ Mean:平均値
 - Geometric mean:幾何平均
 - Standard deviation:標準偏差
 - Variance:分散
 - Coefficient of variation: 変動係数
 - ・ Minimum:最小値
 - ・ Maximum:最大値
 - Pearson correlation (X vs Y):相関係数
 - Percent in gate:指定したゲート内イベントが何%か
 - ・ Percentile distance (5to95): パーセンタイル 5と95の間の値
- Significance test : <u>有意差検定</u>の方法を選択します。



CD4+T 細胞に対する刺激の効果を、IFNgとTNFa の値で示すサマリードット例。緑は刺激、オレンジは未刺激コントロール。

2-11-8統計的有意差検定機能(共通)

統計チャートで統計検定を実施するには、 Plots 設定> Significance test で使用しているデータセットに適した検定方法を選択します。



Cytobankの推定的統計ツールでは、生物学のデータ解析で広く使われているいくつかの統計検定が実装されています。マニュアルゲートの結果についてでも、FlowSOM などの結果についてでも、Cytobank上でそのまま有意差検定ができます。

下がそのリストです。

None				
	# IV	# Groups	Parametric	Paired
Student's t-test	1	2	~	
Mann-Whitney U test	1	2		
Paired student's t-test	1	2	~	~
Wilcoxon signed-rank test	1	2		~
Kruskal-Wallis H test	1	> 2		
One-way analysis of variance	1	> 2	~	
Two-way analysis of variance	2	> 1	~	

2-1 1-8-1

2-11-8-2有意差検定

Cytobank に実装されている有意差検定方法

- Student's t-test : Student's t-test はひとつの独立変数の2つのグループの平均において違いがあるかを判定するのに用いられるパラメトリック検定です。データは正規分布に従うと仮定されています。

- Mann-Whitney U test : Mann-Whitney U test はひとつの独立変数の2つのグループの分 布が等いかそうでないかを判定するのに用いられるノンパラメトリック検定です。データが正規分布に従っているなどの分布の仮定や制限はありません。
- Paired student's t-test : Paired student's t-test は、同一個体での治療前と後など対応 ありデータでの2グループ間に有意差があるかを判定するのに用いられるパラメトリック検定です。 student's t-test と同様 paired student's t-test もデータは正規分布に従うという仮定のも と、2つのグループの平均が等いかそうでないかを判定します。
- Wilcoxon signed-rank test : Wilcoxon signed-ranks test は、ひとつの独立変数の対応 ありの 2 グループについて、その分布が等いかそうでないかを判定するのに用いられるノンパラメトリック 検定です。Mann-Whitney U test と同様 Wilcoxon signed-rank test も 2 つのグループが 同じ分布を持つかどうかを判定します。
- Kruskal-Wallis H test : Kruskal-Wallis test は正規分布でないと仮定されるデータでのノン パラメトリック検定です。ひとつの独立変数について2以上のグループを比較し、すべてのグループの分 布が同じかそうでないかを判定するのに用いられます。
- One-way analysis of variance (One-way ANOVA): One-way ANOVA はひとつの独 立変数の2以上のグループについてのパラメトリック検定で、サンプルグループ間の平均に違いがある かを判定するのに用いられます。
- Two-way analysis of variance (Two-way ANOVA): The two-way analysis of variance analysis はパラメトリック検定で、2以上のグループでの2つの独立変数について同時 に分析することができます。Cytobankでは2つの独立変数の交互作用は考慮しません。

2-11-8-3多重検定と補正

同時に複数の統計検定を行う際には p 値を補正または調整することをおすすめします。この補正は多重 検定問題として知られている、検定を繰り返すことによって偽陽率が増える問題を回避することに役立ち ます。サイトメトリーデータの場合、複数の細胞集団からなるサンプルの免疫システムについて考えます。サ ンプルを比較する際には、各細胞集団について繰り返し統計検定される多重検定が通例ですので、その ような時に有用です。Cytobank では Layout settings で設定される dimension は Table ごとに 処理されます。多重検定補正も各 Table 内で dimension に対して施されます。P-値補正は Table 間では実施されませんのでお気をつけください。

 2グループ検定法の場合(Student's t test, Mann-Whitney U test, Paired student's t test, and Wilcoxon signed-rank test)は、BonferroniとFDR 補正法が Multiple test methods メニューで選択できます。Benjamini & Hochberg procedure は偽検出率 (False discovery rate)を調整できるよう設計された多重検定補正法です。Bonferroni method と比較するとFDR-controlling method は Type1 エラーに対しての厳しさが緩くなっ ています。

- ・ Bonferroni correction : Bonferroni correction は多重検定問題への対処として一般 的な調整方法です。一緒に実行された検定数に基づいて補正を適用します。
- FDR correction Benjamini & Hochberg procedure, a false discovery rate (FDR) correction method
- Kruskal-Wallis H test または Two-way ANOVA test の場合、p-値の調整と同じくすべてのグループのペアワイズ比較を実施するための2グループ検定法と補正法の組み合わせが選択できます。2グループ以上の時には、他のグループに対して有意に異なる特定のグループを特定します。Noneを設定した場合には、結果は独立変数間の有意性を示す全体的な p-値のみになります。
- Kruskal-Wallis H test の場合:
 - Mann-Whitney U test + Bonferroni : Mann-Whitney U test によるすべてのペアワ イズ比較と Bonferroni correction による補正が実施されます。
 - Mann-Whitney U test + FDR : Mann-Whitney U test によるすべてのペアワイズ比 較と FDR correction による補正が実施されます。
- Two-way ANOVA test の場合:
 - Student's t test + Bonferroni : Student's t test によるすべてのペアワイズ比較と Bonferroni correction による補正が実施されます。
 - Student's t test + FDR : Student's t test によるすべてのペアワイズ比較と FDR によ る補正が実施されます。
- one-way ANOVA が用いられる時は、Tukey's HSD test では各グループの独立変数についてのペアワイズ比較が可能です。
 - Tukey's HSD (honestly significant difference) test : Tukey's HSD test は Tukey's range test や Tukey's test という名称でも知られており、ワンステップで多重検 定問題を補正しペアワイズ比較を実行します。どのグループの独立変数が他と有意にことなるか を特定するためにすべてのペアのグループ平均を比較します。

2-11-8-4 有意差検定の実行方法

推定的統計ツールは illustration Editor に実装されています。

メニューバーの Illustrations> New illustrations で新規の illustration を作成し、Plots にて Plot type を Box か Violin, Bar, Line, Summary dot に設定します。Significance test のドロップダ ウンリストを開くと各検定法についての独立変数数(#IV)、グループ数(#Groups)、パラメトリック かノンパラメトリックか(parametric)、対応ありかなしか(Paired)が一覧で表示されます(下 図)。ここからデータセットに適した検定法を選択します。

Actions	🌱 Sample tags	🗘 Gates	ដាំ Advanced	analyses	🛦 Illu	ustrations	
🛈 Illust	ration Editor —	Statistical Infer	ence 🧷				
l	🕌 Plots 🚽 🛛 🖡	🕈 Layout 🚽	💾 Save 🚽	🗊 Ex	port -		
	(i) Plot type	∮ ∔∯ Box	ď				
	Show dots	\bigcirc					
	Y axis scale	Local					
	Width	Fixed					
	Statistic	Percent of popula	ti				
	Population A	T regs			3		
	Population B	CD45+ cells					
	Significance test	None	•	+ cells	2.5		
			None	e			
				# IV	# Groups	Parametric	Paired
		-	Student's t-tes	t 1	2	~	
		Ma	nn-Whitney U test	t 1	2		
		Pair	ed student's t-tes	t 1	2	~	~
		Wilcoxor	signed-ranks test	t 1	2		~
		Kri	uskal-Wallis H test	t 1	> 2		
		One-way a	nalysis of variance	9 1	> 2		
		Iwo-way a	nalysis of variance	e 2	> 1	~	

2-11-8-5有意差検定を含む統計ツールの設定手順

- ① Illustration Editor の Plots でどの統計値と検定を用いるかを選択します。
 - i. Plot type のプルダウンメニューで Summary plot に属するプロットを選択します。選択したプロットに関連した設定項目に表示が変わります。
 - ii. Statistic ドロップダウンメニューで用いたい統計値を選択します。各サンプルのプロットにはこの 統計値が用いられます。また有意差検定にも使用します。
 - iii. 仮定に基づいて Significance test ドロップダウンメニューから用いる検定を選択します。対応 ありの検定(Paired Student's t test または Wilcoxon signed-rank test)を選択した 場合、Cytobank は sample tags の情報をもとに自動的に対応の情報を抽出して使用し ます。
 - iv. 各プロット内で複数の 2 グループ検定を実行するか、Kruskal-Wallis H、ANOVA (One-way または Two-way) を選択した場合には、適切な多重検定法を選択します。
- ② Layout 設定でプロットのレイアウトを設定します。
 - ・ 比較のために Sample tags でグループ分けを定義した dimension を独立変数(I.V.)として割り当てます。デフォルトでは Layout の上から 2 つ目 Subgroups に設定することで適用

されます。※独立変数(I.V.)が Layout の1つ目に設定される場合もありますため、確認ください。

- ii. 2つ以上の Sample tags を表示対象に選択し、I.V.の比較グループに設定します。
 - Layout で2グループ検定法を選択した場合 (Student's t-test, Mann-Whitney U test, Paired student's t-test, または Wilcoxon signed-rank test)、
 Layout で I.V.に設定した dimension が2つのグループ (Tags) が選択されている ことをご確認ください。グループ数が2以外の場合、Layout 内にエラーメッセージが表示 されます。
 - two-way ANOVA を選択した場合は2つの独立変数(I.V. #1とI.V. #2: Layout の一番上と2番目の項目)に各々dimensionを割り当てます。

🚯 Illustration Editor — Statistical Inference 🧪		 Illustration Editor 	— Statistical Infe	rence 🥒				
Step #1	🕌 Plots 🚽	∰* Layout ⊸	💾 Save 👻	潘 Plots 👻	₽ Layout →	💾 Save 👻	Export -	
	i Plot t	ype	ď		i			ď
	Show d	lots		Step #2	X-axis groups	Blood types	1 of 4	
	Y axis so	cale Local			Subgroups	Conditions	2 of 2	
	Wi	dth Fixed			Table 1	Add dimension	1	
	Stati	stic Percent of popula	ati			-		
	Populatio	on A T regs	v			R Lavout Preview		
	Populatio	on B CD45+ cells	-			E Edjournement		
	Significance t	test None				ų		
			to					
			95 VE					
			2 State					

2-11-8-6結果の表示方法

有意差検定の結果は2つの表示形式があります。

① プロット上に P 値かアスタリスクシンボルで表示

プロット内または横に P 値かアスタリスクシンボルで表示します。表示される値またはアスタリスクは 選択された検定方法に依存します。P 値の範囲とシンボル(アスタリスクの数)の関係は以下の 通りです。

P-value	Symbol
0.01 <p<0.05< td=""><td>*</td></p<0.05<>	*
0.001 <p<0.01< td=""><td>**</td></p<0.01<>	**
0.0001 <p<0.001< td=""><td>***</td></p<0.001<>	***
<0.0001	***

- 2グループ検定(Student's t test、Mann-Whitney U test、Paired student's t test、Wilcoxon signed-rank test):
 - Show p-values で表示の項目を選択した場合、比較グループに跨るラインとその上
 中央に数値またはシンボルが表示されます。すべての比較を表示するか、有意差のあるものだけとして p <0.05 だけ表示するかを選択できます。
 - Multiple testing で補正方法を選択した場合は、補正後の P 値またはシンボルが 表示されます。



図: Student's t-test 検定結果を箱ひげ図上に表示. 補正なし(左)とBonferroni 補正(右)

- One-way ANOVA、kruskal-Wallis H test:
 - Show p-values で表示の項目を選択し、比較したいグループを X-axis group に 設定した場合、グループ間に有意差のあるものの上部に、グループを跨る線と P 値また はシンボルが表示されます。 X-axis group 以外に設定した場合、すべてのグループ 間に有意差があるものの上部に線と P 値またはシンボルが表示されます。
 - Multiple testing を選択した場合は、プロット横に表が表示されます。この表は有意
 差(p<0.05)のあったペアワイズとその検定結果をシンボルで表示します。



図:棒グラフ上に表示された one-way ANOVA 結果. Tukey HSD なし (左) と Tukey HSD 適用後 (右)

- Two-way ANOVA :
 - 有意差のある独立変数名がその検定結果のシンボルとともにプロット横の表に表示されます。
 - Multiple testing の設定はプロット表示には反映されず、その下の Statistic Inference table にのみ反映されます。



※データが欠損していたり、変化がない場合、Show p-values 設定が P values(all)または Symbols (all) となっている場合は NaN が表示されます。

② Statistical inference 表での検定結果の表示

有意差検定を行った場合に illustration のプロットの下に表示される Statistical Inference というセクションで表になっています。

- 2グループ検定: Student's t-test, Mann-Whitney U test, Paired student's t test, Wilcoxon-signed rank test では、表の各行がひとつの比較にまとめられています。
 - Mean A と Mean B: A と B に割り振られている独立変数グループについて算出された平均値(Plots 設定集計方法で Mean が選択されている場合)
 - Median A と Median B: A と B に割り振られている独立変数グループについて算 出された中央値(Plots 設定集計方法で Median が選択されている場合)
 - SD A と SD B: 独立変数グループ A と B の 各 々の標準 偏差
 - ・ nAとnB(nは数値):独立変数グループAとBの各々のデータポイント数
 - p-value:補正なしのP値
 - Adjusted p-value: 多重検定や補正後のp値. Kruskal-Wallis H, Oneway or Two-way ANOVA test で multiple testing を実行した場合は補正後のP値だけが表示されます。
 - ・ Significance p-value:補正なし P 値に対応したシンボル
 - ・ Adjusted significance p-value:補正後のP値に対応したシンボル
- Kruskal-Wallis H, One-way ANOVA, Two-way ANOVA: これらの検定方法を 選択し、multiple testing を使用した場合は2つの結果の表が表示されます。プロット下 に表示される Statistical inference の表は上述の2グループ検定と同様です。上部 (プロットの横)の表では、検定方法名が記され、有意差のあった独立変数グループのすべ

ての組が表示されます。

- Table 名: Layout 設定の Table を使用している場合、その名前が表示されます。
- X-axis 名: Layout 設定で x-axis を1I.V.に割り当てている場合、その名前が表示されます。 Two-way ANOVA では x-axis が I.V.に割り当てられます。
- ・ 独立変数名:グループ比較の独立変数名
- ・ P-value: 各組み合わせ、行の P 値
- ・ Significance p-value: P 値に対応したシンボル

例えば検定方法を One-way ANOVA で、multiple testing を Tukey HSD にする と、上(プロット横)の結果表は one-way ANOVA での結果、下の statistical inference 表は Tukey HSD の結果が表示されます。

2-11-9統計グラフのエクスポート

統計チャートの作成が完了したら、Cytobank experiment 内に illustration として保存(デフォルト では自動保存になっています)するだけでなく、illustration editorの図として PDF、PNG、SVG 形式 でエクスポートできます。プロット下部にある統計検定の表と統計表は各々csv ボタンにて csv 形式でエ クスポートすることも可能です。

2-12共有・プロジェクト・権限

Cytobank のデータや Experiment は他のユーザーと安全に共有することができます。Experiment を まとめたプロジェクトには、共有メンバーに権限付けもできます。

※共有できるユーザーは同一サーバーのユーザーに限ります。Premium Cytobank ユーザー同士共有 できます。Enterprise Cytobank ユーザーは同じ Enterprise サーバー内のみの共有となります。異な る Cytobank サーバーヘデータ移行されたい場合はご相談ください。

2-12-1ユーザーコネクション(Premium cytobank のみ)

アカウント間のアクセス許可を双方で確認し、user personal identification information (PII) privacy に準拠するため、Experiment や Project の共有をする前にユーザーコネクションを確立します。

- 2-12-1-1-1ユーザーコネクションにユーザーを追加
 - ① My profile ボタンをクリック> account setting をクリックし、My account ページを開きます。
 - スクロース下の User connections セクションの Request new user connection のボックス にコネクションに入れたいユーザーの Cytobank 登録 E-mail を入力します。

My account	/ Edit			
User connect	ions			
Request ne	w user conne	ction		
user@email.com	🕀 St	ubmit		
Existing co	nnections			
User	Date connected	# shared experiments	# shared projects	
gara	20-April-2022	1	0	20
0	pre-existing	0	0	20

③ リクエストを受けたユーザーはメール通知を受け取るとともに Experiment manager に User connection requests が表示され、Accept か Decline かを選択できます。

Experiment Manag	er	٩	🛛 🗹 Gro	S Group linked		
New experiment	• New experiment					
Q All	957	User	Email	Sent at	Respond	
Mine .	861	Bob Jones	bobjones@gmail.com	Sep 13, 2021	✓ Accept X Decline	
Shared with me Public	88 23	Jimmy Smith	jimmy_the_smith@yahoo.net	4 hours ago	✓ Accept X Decline	

 Accept されると Existing connections リストに追加され、Experiment や Project の共有 が可能となります。

※ユーザーコネクションが確立された時点では共有が可能となっただけで、アカウント内の Experiment はこの時点では共有されていません。

2-12-1-2ユーザーコネクションからのユーザーの削除

- ① そのユーザーと共有しているすべての Experiment で共有を解除します。
- ② My account> Existing connections でそのユーザーとの#shared experimentsと #shared projects が0となっているのを確認します。
- ③ その右の×マークをクリックします。

2-12-2データセキュリティ・デフォルトのアクセス権限について

Cytobank ヘアップロードされたデータは基本的にはアップロードしたアカウントでのみ閲覧・操作可能です。他のアカウントとの共有には以下の操作によるいずれかの許可が必要となります。

2-12-3Experimentの共有

① 共有したい Experiment を開きメニューバー右の Sharing Private をクリックします。

TEST Actions	🌱 Sample tags	🗘 Gates	ភីង Advanced analyses	L Illustrations	The Scale	es 🛆 Compensation	Sharing Private
(2)	Experim	ent Sha	ring の Share	with a User 7	ボックスに共有者の Cvt	obank ユーザ	-名ま

② Experiment Sharing of Share with a Oser かうえに共有者の Cytobank ユーリーを たは登録 Email アドレスを入力します。

③ すでに共有したことのあるユーザーやユーザーコネクションに含まれているユーザーの場合、候補者リ ストが表示されるので、選択します。



ユーザーコネクションのないユーザーの場合は、No users found の下に表示される Connect with a user ボタンをクリックし、ユーザーコネクションに追加します。

share with a user				
	Q, sf)		
	No users found			
Tŀ	Connect with a user	7 		

④ 共有者はフルアクセスとなります。アカウント容量は所有者に帰属し、共有者の容量は消費しません。

※共有者が再解析する場合は、illustrations や Advanced analysis はアカウントごとに分かれて 作成されますが、gating や Scale 等は上書きになってしまうため、Clone を作成して別にすることもお すすめします。

2-12-4アクセス権限を付加した共有ができるプロジェクトの作成と管理

① 画面上部の Projects から New を選択します。



- ② Name にプロジェクト名を入力します。
- ③ 必要であれば Description を記入します。
- ④ Sharing level でフルアクセス以外のプロジェクトの共有者(メンバー)の illustrations への制 限を選択します。
 - View Only:保存された illustrations の閲覧のみ
 - Illustrations: フルアクセス以外のユーザーも illustrations を作成可能
- ⑤ Cloning level で他のユーザーが Clone を作成するときの制限を選択します。
 - Disable:メンバーの Clone の作成と生データのダウンロードを許可しません。
 - Clone Files Only: 生データ以外 (解析結果など)のダウンロードや Clone 作成は許可しません。 生データのみの Clone とダウンロードは可。
 - Full Clone: 生データ以外も含めた Clone の作成を許可します。

⑥ Create Project ボタンを押します。

New project	
Name	Project demo
Description	demo
Sharing level	View only View of View
Cloning level	Disabled
	Create Project

- ⑦ 詳細設定が表示されます。
 - Edit Project Members で付与する権限別にユーザーを追加し、プロジェクトのメンバーに することができます。各々のボックスに Cytobank 登録 Email アドレスを入力し、候補リスト から選択することで追加できます。各メンバーの右の×マークをクリックするとメンバーから外す こともできます。候補が表示されない場合、ユーザーコネクションにユーザーを追加ください。
 - Manager: プロジェクトメンバーやプロジェクト Experiment の追加や削除を管理できます。また、共有や Clone 権限の変更もできます。
 - Leaders: Manager によって設定された共有・Clone 制限下でプロジェクトの
 Experiment にアクセスできます。アクセス権限は Members と同じです(組織体制に対応するために名称を分けた)。
 - Members: Manager によって設定された共有・Clone 制限下でプロジェクトの
 Experiment にアクセスできます。アクセス権限は Leaders と同じです(組織体制に対応 するために名称が分けられている)。

⑧ Edit Experiments でプロジェクトに含める Experiment を選択することができます。リスト上部のボックスでリストにフィルターをかけることができます。左のリストがアクセス可能な Experiment、右のリストがプロジェクトに含まれている Experiment になります。リスト下の Add Experiment(s)/Remove Experiment(s)ボタンで出し入れができます。

ytobank	Experiments	Projects				Help
Edit project	t <i>'Project dei</i>	mo'				
Edit Details						
	Name	Project demo				
	Description	demo				
					.::	
	Sharing level	View only	~			
	Cloning level	Disabled	~			
Update p	project					
Edit Project	Members					
Managers	an add and ramo	a project members leaders	managars	and experiments. Managers can a	ter sharing and cloping p	ermissions for membe
and leaders.	an add and remov	re project members, redders, r	munugers, c	and experiments. Managers can a	ter sharing and cioning p	annissions for membe
🖸 Yuko	Nakane					
Add Mar	nager:					
[a						
Leaders	access project ex	maximants according to the s	haring and	cloning permissions set by Manag		
Add Lea	der:	periments according to the si	nunng unu	cioning permissions set by munug	615.	
Q						
Members						
Members co	an access project e	experiments according to the	sharing an	d cloning permissions set by Mand	igers.	
Add Mer	mber:					
[4						
Invite a ne	wuser					
Add experim	ents currently un	assigned to projects:				
Filter displa	yed:	lassigned to projects.				
[12571] 2	- 0210210Cutof P	PMC/TIL Paul Fisher - for f	fler w	Remove experiments from the	s project:	
[13572] 20	0210210Cytof P 0210210Cytof P	BMC/TIL Paul Fisher - fcs f	files w	Filter displayed:		
[13754] A	PI demo					^
[13832] TI	EST					
[13833] TI	EST2					
			\sim			
						~

※例) プロジェクトAは5つの Experiment からなっており、各 Experiment に解析担当がいる(解析担当者5名+PI)。各解析担当者は、担当する Experiment にはフルアクセスできるが、担当以外の Experiment については生データのみの Clone 作成かダウンロードしかできないようにしたい。この場合は、2-10-3 でプロジェクトAを作成し、以下の設定にします。 -Cloning level を Clone Files Only に設定

-Edit Project Members で、各解析担当者をLeaders(または Members)に割り当て -Edit Experiments で 5 つの Experiment を選択

その後、2-10-2の方法で各 Experiment を各々の解析担当者とだけフルアクセス共有します。

3 Advanced analyses (機械学習解析)

3-1機械学習による解析

サイトメトリー技術の発達により測定から得られるデータの複雑性と規模が拡大する中で、バイアスを減ら しマニュアルゲーティングでは気づけなかったサブセットへの気づきを得ることができる手法として、機械学習 による解析が用いられています(Keyes, T.J et. al. "A Cancer Biologist's Primer on Machine Learning Applications in High-Dimensional Cytometry". Cytometry, 97;782-799)。 Cytobank では教師なし学習である次元削減アルゴリズムを4種類(<u>tSNE-CUDA、UMAP、opt-</u> <u>SNE、viSNE</u>)とクラスタリングアルゴリズム2種類(<u>FlowSOM、SPADE</u>)、教師あり学習 <u>CITRUS</u> を実装しています。

また、従来の手法であるマニュアルゲーティングは結果のばらつきの主な要因のひとつであるとされてきました(Maecker et al. 2005)。解析の信頼性と再現性を高めるために、データ解析の集中化やゲーティングストラテジーの標準化、オペレーション手順の標準化などが提案されていますが、時間のかかるものとなってしまっています(Macchia et al. 2020)。サイトメトリーデータの機械学習による解析は非常に有用ですが、仮説や知見に基づく独自のゲーティングストラテジー解析を機械学習で自動化する方法は欠けている現状があります(Hu et al. 2021)。Cytobankでは、少数のサンプルでの独自のゲーティングストラテジーから教師あり機械学習によってモデルを構築し、プログラミングなしで、機械学習による自動ゲーティングによって、解析者間のポピュレーション結果のばらつきを軽減するとともに、解析ワークフローを合理化することができます

これら機械学習解析はメニューバーの Advanced analyses から実行することができます。

Actions	🌱 Sample tags	A+ Data QC	🗘 Gates	🖧 Advanced analyses 🔹	L Illustrations
? Experim	ent summary	Clone	n Ed	Dimensionality reduction	
-	Name: Cloned from:	0- Demo Gating illu 0- Demo Gating illu	ustrations (Cloi Istrations	Clustering	
_	Project: Labels:	None Demo		CITRUS	
Notes				Automatic gating ④ ਨੇ New analysis	

3-2自動ゲーティング

3-2-1 自動ゲーティングとは

Cytobankの自動ゲーティング機能は、ユーザー独自の細胞集団の同定やゲーティング時の境界線を用いることを可能にします。この機能は2つのステップから構成されています。ひとつめのステップは、マニュアルゲートされたデータを用いてモデルをトレーニングします。モデルはユーザーの与えたゲート情報から細胞をどのように細胞集団に割り当てるかを学習します。ふたつめはinference(推論)ステップと呼ばれ、モデルを新規データに適用して、マニュアルゲートの調整作業なしに各細胞イベントを細胞集団に分類(割り当て)します。



最低 10 ファイルのマニュアルゲーティングされたデータがモデルのトレーニングには必要です。必要な場合、 ファイルや親ポピュレーションによって微調整した tailored gate 機能も使用してください。適切なゲーティ ングストラテジーが確立したら、Gating Editor の Apply gates ボタンを押します。 作成されたモデルは、他の Experiment に適用することが可能なだけでなく、そのモデルを含む Experiment を共有することで他のユーザーとも共有することができます。

良好に動作するモデルをトレーニングするには、解析データに含まれているばらつきがトレーニングデータにも 含まれていることが重要です。ばらつきとは、試薬のロットや機器の設定などのテクニカルなものとサンプル毎 のマーカー発現レベルの違いなどの生物学的なものとを含みます。また、このばらつきがモデルトレーニング の各段階であるトレーニング、バリデーション、テストの全てのセットに反映されるトレーニングデータセットで モデルの学習が行われる必要があります。

3-2-2トレーニングデータの準備

- 実行可能な範囲でなるべく多くのサンプルをトレーニングデータに使用します(最低 10 ファイル)。
- トレーニングデータにはできる限り適切なゲーティングをします。
- トレーニングデータには予想されるばらつき、つまり試薬や機器設定、サンプルの生物学的要因などが可能な限り反映されるようにします。

3-2-3モデルトレーニングの設定・実行

- メニューバー> Advanced Analysis > Automatic gating > Train new model を選択
 > Create をクリックします。
- ② Analysis name に名前を入力します。Create ボタンで設定画面に移動します。



③ Population でモデル構築の対象とするポピュレーションを選択します。選択されたポピュレーション の上流のポピュレーションは、そのポピュレーションを分類するのに必要な段階として自動的に対象 となります。選択されたポピュレーションを同定するのに使用した全ての階層のチャネルがモデルのト レーニング時に考慮されます。デフォルトでは、各ポピュレーションの重要度は全て同じとされていま す。モデルトレーニングの設定では、ポピュレーションの重み(KPI weighting)の値を増加させ ることもできます。weight 値を増加させた特定のポピュレーションは、全体の KPI を算出する際 にそのポピュレーションの結果への影響を大きくすることとなります。Magnificationを増加させて 複数のモデルをトレーニングして比較する時に、最も重要なポピュレーションの weight を増加させ ると、どのモデルが最も求めるパフォーマンスかを知るのに役に立ちます。



④ FCS Files で、トレーニングに使用するファイルを選択します。
 必要なファイル数は最低 10 ファイルですが、実行可能な範囲でなるべく多くのサンプルをトレーニングデータに使用することをおすすめします。
 ファイルには予想されるばらつき、試薬や機器設定、サンプルの生物学的要因などが可能な限り反映されるようにします。

Cyto	Autopring theme Autopring theme RACione IM Basic Standardization St	E Select FCS files	×
	ictions 🌱 Sample tags		
< 2	Autogate training analysis	Q, All None Inverse 12	
	Analysis name: Qianju	Site 2 No wash 1]Tube1.fcs_file_internal_comp_Ungated.txt - [Site 2 No wash 1]Tube1.fcs_file_internal_comp_Ungated Site 2 No wash 1]Tube1.fcs_file_internal_comp_Ungated	
	Populations	1)Tube2.fcs_file_internal_comp_Ungated	
	CD3-CD19-	[Site 2 No wash 1]Tube3.fcs_file_internal_comp_Ungated.txt - [Site 2 No wash 1]Tube3.fcs_file_internal_comp_Ungated	
	CD4+	[Site 2 No wash 1]Tube4.fcs_file_internal_comp_Ungated.txt - [Site 2 No wash 1]Tube4.fcs_file_internal_comp_Ungated	
	CD8+	[Site 2 No wash 1]Tube5.fcs_file_internal_comp_Ungated.txt - [Site 2 No wash	
	CD14+	1]lube5.tcs_tile_internal_comp_Ungated	
	CD19+ CD45+	[Site 2 No wash 2] lube1.fcs_hle_internal_comp_Ungated.txt - [Site 2 No wash 2]Tube1.fcs_file_internal_comp_Ungated	
	CD45+ CD3+	[Site 2 No wash 2]Tube2.fcs_file_internal_comp_Ungated.txt - [Site 2 No wash 2]Tube2.fcs_file_internal_comp_Ungated	
	CD56+CD16+NK cells CD56bright NK cells	[Site 2 No wash 2]Tube3.fcs_file_internal_comp_Ungated.txt - [Site 2 No wash 2]Tube3.fcs_file_internal_comp_Ungated	
	Cells	[Site 2 No wash 2]Tube4.fcs_file_internal_comp_Ungated.txt - [Site 2 No wash 2]Tube4.fcs_file_internal_comp_Ungated	
		[Site 2 No wash 2]Tube5.fcs_file_internal_comp_Ungated.txt - [Site 2 No wash 2]Tube5.fcs_file_internal_comp_Ungated	
		[Site 2 No wash 3]Tube1.fcs_file_internal_comp_Ungated.txt - [Site 2 No wash 3]Tube1.fcs_file_internal_comp_Ungated	
	14 of 14 selects	Site 2 No wash 3]Tube2.fcs_file_internal_comp_Ungated.txt - [Site 2 No wash	
			Close
	Event compline		ciose

⑤ Event Sampling で、データのサンプリング方法を選択します。
 Equal: 各ファイルの細胞数の違いは無視して同数の細胞をサンプリングします。
 Proportional: 各ファイルの細胞数の違いの比を反映してサンプリングする細胞数を決めます。
 Use all events:すべての細胞を用いる。

Event sampling	
 Equal Proportional 	Desired total events: 500000
O Use all events	Actual events to use: 3,093,243 (no sampling necessary)

自動ゲーティングモデルのトレーニングで使用するダウンサンプリングは、Smart downsampling という手法で、ランダムピックアップと異なり、トレーニングセットサンプル全ての中のポピュレーションの 多様性に最大限最適化されています。マニュアルゲーティングで同定された各ポピュレーションが全 て漏れなく確実にトレーニングにサンプリングされるようになっています。これにより希少な細胞のポピ ュレーションを見逃すリスクを回避できます。



Optional settings:

6 Random Seed

機械学習モデルのトレーニングには乱数が多く組み込まれており、トレーニングデータのサンプリン グやモデルトレーニングのパラメータなどが含まれます。トレーニング結果の再現性を確認したい 場合は、乱数シードの設定をランダムの変わりに任意のシードにすることで、複数のトレーニング でも同一のサンプルとパラメータ選定で実行することができます。しかし、一般的にはランダムで 実行することを推奨します。 ⑦ Optimal clusters

トレーニングデータセットがいくつのグループになる背景を持っているかの見積もりを設定します。 (例:4機器からのファイルの場合4グループ)

(8) Training magnifications

magnification を増加させると、異なるパラメータを持つ複数のモデルを同一のトレーニングデ ータで実行することができ、最も KPI の高かったモデルが結果として返されます。 Magnification を1より大きい値にして特定のポピュレーションの重み(weight)を増加さ せると、そのポピュレーションにとってより適したパフォーマンスをするモデルが選択されるようにする ことができます。注意点としては、magnification は計算時間がその値に比例して長くなること です。

Random see	ed 🚺	
(Random)		
Optimal clu	sters 🕕	
3		
Training ma	gnification 🕕	
1		

(9) Compensation & Data scaling

自動ゲーティングモデルはコンペンセーション済、スケール調整済のデータのパターンを学習します。 他のフローサイトメトリー解析と同じく、適切なコンペンセーションとスケール調整は重要です。コンペ ンセーションおよびスケール調整のクオリティは、モデルのトレーニングに使用するデータセットとモデル を適用して自動ゲーティングする新規データセットとで同等に処理されている必要があります。自動 ゲーティングを行う前には必ずコンペンセーションとスケール調整が適切に処理されているかをご確 認ください。

Test view

 ① Create an experiment with blind test files and their inferred populations のボッ クスにチェックを入れると、モデルトレーニングのブラインドテストで用いたファイルとマニュアルゲーティン グ、自動ゲーティング結果を含めた新規 Experiment として作成します。モデルトレーニングのテ スト結果の評価・確認をすることができます。

> **Test set review** (i) Create experiment with blind test files and their inferred populations

3-2-4新規データへのモデルの適用方法

トレーニングしたモデルを用いて新規の FCS ファイルを自動ゲーティングするには、新規 FCS ファイルのある Experiment を開きます。モデルトレーニングを行った Experiment と同じでも別でもどちらでも問題ありません。必ず、スケールやコンペンセーション、パネルとチャネル名設定が適切に前処理されているかを確認してください。

- メニューバーの Advanced Analysis > Automatic gating: Select Run inference using model > Create をクリックします。
- ② Analysis name に名前を入力します。Create ボタンを押すと、設定画面に移動します。



③ FCS Files ボックスで対象の FCS ファイルを選択します。この後選択するモデルを適用して自動 ゲーティングしたいファイルを全て選択します。



 ④ FCS ファイルを選択すると、Trained Autogating Model セクションで自動ゲーティングモデルを 選択することが可能になるのでモデルを選択します。



- ⑤ Clone the gates from the current experiment にチェックを入れると Experiment 内の マニュアルゲーティングを自動ゲーティング適用後の結果 Experiment に引き継ぐことができます。 実行後の自動ゲーティング解析の結果をマニュアルゲーティングと比較して評価するのに便利で す。
- 画面上部緑色の Run Autogate inference analysis をクリックして実行を開始します。実行が完了し結果の Experiment が作成されると登録 E-mail に通知が来ます。



 ・通知メールのリンクをクリックするか、親 Experiment の Experiment summary ベージの
 Linked Experiment などから結果 Experiment にアクセスできます。

3-2-5自動ゲーティング(モデル作成/適用)の結果の評価

トレーニングモデルが作成されると Analysis run info が表示され、作成されたモデルで行ったブラインド テストの結果が表示されます。また、Create an experiment with blind test files and their inferred populations にチェックを入れるとブラインドテストで用いたデータファイルとトレーニングの正解 であるマニュアルゲーティング、自動ゲーティング結果を含んだ Experiment が作成されるので、 illustrations でグラフ化して評価することができます。

作成したモデルを使用して新規ファイルに自動ゲーティングを適用した場合、Clone the gates from the current experiment にチェックを入れると結果 Experiment にマニュアルゲーティングの情報も引き継ぐことができます。

- 3-2-5-1作成した自動ゲーティングモデルのパフォーマンスの評価
- Analysis run info での評価

モデルのトレーニングが完了すると、トレーニング設定ページに Analysis run info が表示され、実 行時間やモデルの KPI を確認することができます。



v.4.100

モデルの KPI の決定には、テストセットの全てのイベントとモデルトレーニングでの全てのポピュレーション において、各イベントが有しているユーザーのマニュアルゲートによるポピュレーション属性と、そのイベント について自動ゲーティングが予測したポピュレーション属性が一致するかどうかを Cytobank が採点しま す。

テストデータセットでのポピュレーションごとの情報は、以下の項目について一覧で表示されます。

Population	Actual	Predicted	True positives	Precision	Recall	F1 score
Singlets	231600	231608.0	231539.0	99.97	99.97	99.97
CD45+	207150	207152.0	207053.0	99.95	99.95	99.95
Cells	207345	207356.0	207254.0	99.95	99.96	99.95
Lymph	74682	74688.0	74617.0	99.9	99.91	99.91
CD4+	34655	34656.0	34622.0	99.9	99.9	99.9
CD45+ CD3+	58349	58356.0	58294.0	99.89	99.91	99.9
CD14+	21049	21048.0	21022.0	99.88	99.87	99.87
CD8+	19901	19907.0	19869.0	99.81	99.84	99.82
CD19+	7680	7679.0	7662.0	99.78	99.77	99.77
CD3-CD19-	8444	8452.0	8422.0	99.65	99.74	99.69
CD56+ CD16+ NK cells	6420	6435.0	6404.0	99.52	99.75	99.63
CD56bright NK cells	552	554.0	541.0	97.65	98.01	97.83

Actual: マニュアルゲーティングでのイベント数

Predicted: 自動ゲーティングでのイベント数

True positives: マニュアルゲーティングと自動ゲーティングと両方ともがこのポピュレーションであるとしたイベント数。真陽性。

Precision: 自動ゲーティングがこのポピュレーションと予測した全イベントのうち真陽性だったものの割 合(%)。値が大きいほど偽陽性または第一種過誤が少ないことを示します。

Recall: マニュアルゲーティングでこのポピュレーションとされたイベントのうち真陽性だったものの割合 (%)。値が大きいほど偽陰性または第二種過誤が少ないことを示します。

F1 score: Precision (適合率)とRecall (再現率)の調和平均で、次の式で計算されます。



 $2x \frac{(Precision \ x \ Recall)}{(Precision \ + \ Recall)}$

経験的に、F1 スコアが 90%以上の場合は very good、80 から 90%は good、50 から 80%は OK、50%以下は not good を示すとされています。

一覧表の他に、ポピュレーション毎の予測スコアについてテストセットファイルの分布をボックスプロットで確認することも可能です。



- Create experiment with blind test files and their inferred populations チェックボック スにチェックをいれて実行していた場合での作成された Experiment での評価
- 作成された Experiment へは、自動ゲーティングの setup ページの View created experiment からアクセスすることができます。

x

```
▲ Autogate training analysis setup Copy settings Analysis complete! ✓ View created experiment
② 自動作成された Experiment には、ブラインドテストセットに割り当てられた FCS ファイルと自動ゲ
ーティングポピュレーションごとにそのポピュレーションのためのチャネルパラメータが「auto_gate_ポピュ
レーション名」で追加されています。Gating Editor の軸パラメータを開くと確認することができます。
```

channel	FSC-A	Selecte
channel		
⊷ Lo	ViaKrome FVD 808 [FL21-A] FSC-Width	^ G
	TIME auto_gate_Singlets auto_gate_LIVE auto_gate_Cells auto_gate_Leukocytes auto_gate_Lymph auto_gate_Granulocytes auto_gate_Mono auto_gate_Bcells auto_gate_CD19- lymph	A A A A A A A A A A A A A A A A A A A
	<	``` ,

③ この追加チャネルでクラスターゲートとして同定され、「ポピュレーション名_auto_cluster_gate」という名前のゲートが自動的に作成されています。



④ Population manager では自動ゲーティングで作成されたポピュレーションには「ポピュレーション名 [auto]」と名前がつけられています。

Population Manager	Populatio	in Tree	Population	Sunburst													
Activate Population	Delete Popula	tion Cop	r Population	Add Populat	on (Q, lym	ph		Clear	Q. Filler Gal	83	Clear						
↓Population \Gate→	Singlets	Cells	CD45+	Lymph	CD45+ CD3+	CD4+	CD8+	NK marker (UR)	NK marker (UL)	NK marker (LL)	NK marker (LR)	CD19+	CD3- CD19-	CD14+	Singlets_auto_cluster_gate	Cells_auto_cluster_gate	CD
Lymph		2															
Lymph [auto]																2	
4																	

- ⑤ <u>illustrations 機能の統計グラフ</u>を用いて自動ゲーティングモデルのパフォーマンスを検証するには、 マニュアルゲーティングのポピュレーションと自動ゲーティングのポピュレーションを event count で比較 します。
- ⑥ Illustrations メニューを開いて自動作成された Autogating boxplot…という名前の illustration を開きます。



- ⑦ 特定のポピュレーションについてのブラインドテストのパフォーマンスを検証するには、Autogating boxplot の illustration editor メニューの <u>Save >Clone new copy でコピーを作成した</u>後そ の他のポピュレーションの表示選択を解除するか、新規 illustration を作成します。
- ⑧ 特定のひとつのポピュレーションについてマニュアルゲーティングと自動ゲーティングの結果をボックスプロットで比較するには、Plots 設定画面の Show dots と Show lines をオンにします。プロットの統計値は csv ファイルとしてエクスポートすることができ、マニュアルゲーティングと自動ゲーティング結果の相関性を他の専用ソフトウエアで詳しく検討することもできます。

) Plots	×	() Layout	×	
Plot type		X-axis groups		
Show dots		Subgroups ① Populations 20	f 29 🔛 ×	
Dot size 6		Table 1 Add dimension		
Show lines				
Gray shading				
Yaxis scale Local				
Width Auto-expand				
Statistic Event count				
Significance test None				
		Ef Layout Preview		
a				
		8000		
		6000		
		count		
		0 tu 4000		
		2		

- ⑨ イベント(細胞)レベルで結果を図示するにはオーバーレイドットプロットを使用します。
- Plots 設定で Plot type を Dot に、Overlaid をオンにして、Layout 設定の Overlaid に

 Populations が割り当てられるようにします。Populations では表示したい自動ゲーティングポピュレーションのみにチェックを入れて選択します。
- マニュアルゲーティングで使用した X、Y 軸を設定して Show gates をオンにすると、自動ゲーティン グで作成した各ポピュレーションのオーバーレイプロットと、マニュアルゲーティングでどの位置に境界線 を引いたかを同時に確認することができます。<u>Tailored par file ゲーティング</u>をしている場合は、最 初のファイルのゲート位置が表示されます(統計値等は Tailored の位置で計算されます)。
- ② プロットの解像度を変更したい場合は Plots 設定の Size や Dot size で調整できます。



- 3-2-5-2モデルを適用した自動ゲーティング結果の評価
- ① トレーニングモデルを適用した自動ゲーティングの結果 Experiment には、Set up 画面の View created experiment をクリックしてアクセスします。

Autogate inference analysis setup Copy settings	Analysis complete!	View created experiment	
---	--------------------	-------------------------	--

② 結果 Experiment の FCS ファイルには、自動ゲーティングパラメータがチャネルに追加されており、その名前は auto_gate_ポピュレーション名(トレーニング時にマニュアルゲートで与えられたポピュレーション名)となっています。Gating Editor ではこのチャネルによる自動ゲートが作成されており、そのゲート名はポピュレーション名_auto_cluster_gate となっています。



- この自動ゲートと一緒にポピュレーションも作成され、ポピュレーション名[auto]と名前がつけられます。 Population manager や Illustration editor、 Population treeでこれまでのポピュレーションと同様に扱うことができます。
- ④ マニュアルゲーティングと比較したい場合は、Population tree と Sunburst を用いると便利です。



- ⑤ また illustration Editor では統計比較も可能です。自動作成されている Autogating box plot of event counts という名前の illustration を開きます。
- ⑥ Layouts 設定の subgroup に populations を割り当て、1 つの自動ゲーティングポピュレーションと対応するマニュアルゲーティングポピュレーションの計2 つを選択します。
- ⑦ Plots 設定で、Show dots と Show lines をオンにします。

 Plots 	×	() Layout	
Plot type		X-axis groups	
Show dots		Subgroups ① Populations 2 o	f 29 🔛 ×
Dot size 6		Table 1 Add dimension	
Show lines			
Gray shading			
Yaxis scale Local			
Width Auto-expand			

- ⑧ 統計的有意差検定機能で自動ゲーティングとマニュアルゲーティングの相違を確認します。
- ⑨ イベント(細胞)レベルで結果を図示するにはオーバーレイドットプロットを使用します。
- Plots 設定で Plot type を Dot に、Overlaid をオンにして、Layout 設定の Overlaid に

 Populations が割り当てられるようにします。Populations では表示したい自動ゲーティングポピュレーションのみにチェックを入れて選択し、マニュアルゲーティングで使用した X、Y 軸を設定します。



- Show gates をオンにすると、自動ゲーティングで作成した各ポピュレーションのオーバーレイプロット と、マニュアルゲーティングでどの位置に境界線を引いたかを同時に確認することができます。
 <u>Tailored par file ゲーティング</u>をしていた場合は、最初のファイルのゲート位置が表示されます(統計値等は Tailored の位置で計算されます)。
- ② プロットの解像度を変更したい場合は Plots 設定の Size や Dot size で調整できます。

3-2-6自動ゲーティング後の解析例

- 3-2-6-1統計グラフを用いたグループ間比較
- ① Sample tags 機能で、各 FCS ファイルに比較したいグループ(群)の情報(例えば、タイムポイントや刺激や治療の有無や前後、疾患や変異の有無、洗浄の有無など)をアノテーションします。

Illustration Editor で新規の illustration を作成し、Plots 設定で Box plot などの統計プロットを選択後、Layout 設定の Subgroups に比較したい sample tags の dimension を設定します。



- Plots 設定の <u>Significance test</u> で使用したい統計検定方法を選択し、グループ間の有意差を 検討します。
- 3-2-6-2次元削減マップでのバックゲーティング 自動ゲーティングで作成されたポピュレーションを次元削減マップ上にオーバーレイ表示することで、デー タセットの全容をクリアに包括的に確認・把握することができます。また、使用したモデルが適用した新規 データでも問題なくワークしているかを確認するのにも役立ちます。
- 自動ゲーティングの結果 Experiment で、メニューバー> Advanced Analyses > New
 <u>dimensionality reduction</u>を選択します。どの次元削減アルゴリズムを選択するかは、ガイドビデ オやアプリケーションノートを参照ください。
- ② 次元削減アルゴリズムが完了したら、結果 Experiment が自動作成されます。
- 次元削減アルゴリズム結果 Experiment で、メニューバー> illustrations > new illustration を選択します。
- ④ <u>Illustration Editor</u>では、<u>Plots 設定</u>で Plot type: Dot plot、Overlaid をオンにします。
- Layout 設定で Overlaid に Populations を割り当て、自動ゲーティングで作成されたポピュレーションをすべて表示対象とします。



3-2-7トレーニング上限と計算時間の目安、トラブルシューティング

3-2-7-1トレーニング上限と計算時間の目安

Cytobank 簡易マニュアル

他の機械学習アルゴリズム同様、サイズの大きすぎるデータをインプットするとサーバーでの計算処理が Failed となります。モデルトレーニングでは、自動ゲーティングに使用するチャネルパラメータ数、細胞イベン ト数、対象ポピュレーション数が大きく影響します。これらを少なくすると、Failed になる頻度を効果的に 減少させることができます。

v.4.100

モデルのトレーニング時の目安として、249 ファイル、計 3700 万細胞のデータセットを用いた検証結果を 下図に示します。縦軸がモデルトレーニングに使用するチャネル数、横軸がインプットする細胞イベント数 で、緑色が問題なく完了できる範囲、黄色が Failed となる可能性のある範囲、赤色がメモリオーバーで Failed になる範囲です。



黄色の領域の設定がされている場合は、設定画面の Event sampling セクションに Warning メッセージが表示されます。この時点でイベント数や対象ポピュレーションを削減することも可能です。

Populations	FCS	lles		
Cho	ose Solu	orted files	Total events	Choose
CD5+ Ly (Ly) CD19-		CLS-MF10C-005 Dx-B 001-AV16220 B-M 00008 903 2017-11-16 1136 101	56,676	56,676
CD19- CD5-		CLS-MF10C-006 Dx-B 001-AV16220 B-M 00008 907 2017-11-16 1146 105	57,493	57,493
CD19+ B cells		CLS-MF10C-007 Dx-B 001-AV16220 L-N 000089 15 2017-11-16 1214 113	42,876	42,876
CD19+ CD10+		CLS-MF10C-008 Dx-B 001-AV16220 B-M 00008 911 2017-11-16 1156 109	52,524	52,524
CD45+ cells		CLS-MF10C-128 Dx-B 001-AV16220 W-B 00009 669	52,443	52,443
Gra Kappa+		CLS-MF10C-146 Dx-B 001-AV16220 W-B 00010 240	53,945	53,945
Lambda+		CLS-MF10C-148 Dx-B 001-AV16220 W-B 00010	48,279	48,279
10 01 10 selected		334 of 455 mes selected — Sampling 22,07	0,017 total eve	ints .
Event sampling				
O Equal	Total ev	ents in selected files: 22,070,017		
Use all events		Warning – input data size approaching memo	ry limits. If you	ur 🛛

赤色の領域の設定がされている場合、設定画面の Event sampling セクションに、メッセージが表示されます。設定したポピュレーション数での上限細胞イベント数と使用するチャネルパラメータ数が表示されま

す。メッセージに従い、上限細胞イベント数以下になるように変更するか、細胞イベント数を優先したい場合は使用するチャネルパラメータ数がメッセージに記されている数以下になるように対象ポピュレーションを 変更することをおすすめします。

\land Autogate training analysis set	IP Copy settings	Status: New	Run Autogate trai
Gra Kappa+	MLL-MF10C-012 Dx-B 012-BA38564 W-B 0000 0236	62,301	62,301
Lambda+	MLL-MF10C-013 Dx-B 013-BA38564 W-B 0000	47,382	47,382
18 of 18 selected	453 of 453 files selected – Sampling 29,	189,151 total events	ŝ
Event sampling C Equal Proportional Use all events	otal events in selected files: 29,189,151	elected populations i	is
	22,100,000 . The selected populations are defined using Reduce sampled events, or select less populations to red gating channels to 8 or less.	11 gating channels. luce the number of	

モデルのトレーニングに必要とする時間は対象ポピュレーションの数と細胞イベント数に依存します。一般 的なサイズのデータセットであれば、高速で計算処理されます。各細胞イベント数の時のポピュレーション 数に応じた目安時間を下図に示します。



この他に Advanced options の Training magnification を1以上に設定すると、その数値の倍数 分の計算時間が必要となります。

(Rand	om)			
dvar	ced options			
Opti	mal clusters 🛈			
3				
□Cr	eate experiment with l	lind test files an	d their inferred popul	ations
Trair	ing magnification (
1				

トレーニングしたモデルを適用するステップ (Autogating inference run) では、データサイズの上限はありません。

3-2-7-2トラブルシューティング

- モデルトレーニングの設定画面でマニュアルゲートしたポピュレーションが表示されない場合

Cytobank のゲーティング機能(Gating Editor)では、Apply gates ボタンを押すことでそれまで の操作でされたゲートとポピュレーションの情報を Experiment 全体つまり Gating Editor 以外の機 能にも適用し、Experiment gate としてバージョン保存します。自動ゲーティング機能は Gating Editor 外の機能のため、Apply gates ボタンを押した Experiment gate のバージョンが反映され ます。

自動ゲーティングのモデルトレーニングの設定画面でポピュレーションが表示されない場合、 Experiment gate が更新されていない可能性が高いため、Gating Editor に移動して Apply gates ボタンを押してください。

- モデルのトレーニングで、あるひとつのポピュレーションの F1 スコアが0の場合
 自動ゲーティングアルゴリズムではモデルトレーニング時いくつかの工程でランダム性を用います。その際たまにモデルが、あるポピュレーションを認識できないことがあります。このような場合は Copy settings をクリックし Random seed がランダムになっているのを確認して、トレーニングを再実行してください。
- モデルトレーニング結果のプロットを見ると、各ポピュレーションのスコアがばらばらの場合
 マニュアルゲーティングでも難しいようなポピュレーションはどうしてもばらつきが大きくなってしまいます。このような場合はトレーニングデータの一貫性を向上させるため、与えているマニュアルゲート位置を確認して修正をご検討ください。
- モデルのトレーニングでひとつまたはいくつかのポピュレーションだけが F1 スコアが悪かった場合
 発現パターンが変調するポピュレーションはマニュアルゲーティングが難しく、これが F1 スコアの低成績を しばしば招きます。トレーニングデータに与えたマニュアルゲート位置の確認と修正をご検討ください。
 また、前述のように自動ゲーティングアルゴリズムではモデルトレーニング複数の工程でランダム性を使用 するため、使用した乱数の組み合わせによってモデルのパフォーマンスが低い結果となることが稀にありま す。この場合には Copy settings で、使用する seed をランダムにしてトレーニングを再実行してください。
 あるいは影響を受けるポピュレーションの重みを増やして、それらに寄ったモデルトレーニングを実行さ せることもできます。
- トレーニングモデルの適用を実行した時にいくつかのポピュレーションについての自動ゲーティングが失敗する場合

トレーニングによってモデルはデータの情報と構造を学習します。この構造とは、チャネルパラメータのスケ ール設定やコンペンセーション設定によって大きく影響されます。コンペンセーションマトリックスとスケール 設定は完全に同一のものを使用する必要はありません。むしろトレーニングに使用するデータとモデルを 適用するデータのスケールとコンペンセーションは処理後の見え方(データ構造)が整合するように前 処理してください。同一のパネルが使われ、一貫性のある機器設定で測定されたデータの場合、既存 の Experiment のスケール設定をインポートすれば適切です。

- スケールとコンペンセーションの設定が適切であることを確認したのに、モデルでの自動ゲーティングが 完璧でない場合

クアドラント(四分割)ゲートでトレーニングされたモデルを使用する際、細胞イベントが複数のクアドラントに割り当てられることがあります。これは Cytobank がクアドラントゲートの4つを独立した四角形ゲ

ートとして扱うからです。このため自動ゲーティングでは、境界線上のイベントなどは複数のゲートポピュレ ーションに割り当てることが予想されます。
3-3次元削減アルゴリズム(tSNE-CUDA、UMAP、opt-SNE、viSNE)

機械学習の教師なし学習に属するアルゴリズムで、多次元情報が有するデータの背景にあるパターンや 分布を低次元(2次元)上に反映して可視化する手法です。マーカー発現パターンにおいて類似性の 高い細胞イベントが2次元プロット上で近い位置に配置されるように可視化されるため、これまでのマニュ アルゲーティングとは異なるアプローチでのサブポピュレーション解析や比較が可能となります。 Cytobank では tSNE-CUDA、UMAP、opt-SNE、viSNEの4種類が実装されています。

3-3-1軸スケールの調整(前処理)

次元削減アルゴリズムでは演算中にゼロ(0)やマイナス値が発生した場合にエラーとなるため、Log ス ケールでのインプットができない仕様となっています。ほとんどの FCM 機器からの fcs ファイルでは、

Cytobank アップロード時に自動でスケールを Arcsinh に変換するようになっていますが、解析前にご確認ください。

一般的にマニュアルゲーティングなどの解析をするのに適したスケールに設定することで、良好に実行できま す。または、以下の方法でスケールを調整します。

- ① メニューバー右の Scales をクリックします。
- ② Experiment Scales 画面で蛍光チャネルのパラメータの Scale が Arcsinh になっているかをご確 認ください。



- ③ チャネル名の目玉マークをクリックして左にヒストグラム表示します。
- ④ 必要な場合、Arg. (= 0 近傍のリニア表示幅のパラメータ)や min. 、max.の数値を変更し、
 ネガティブの集団がシングルのピークとなり 10^0 付近にくるようにします。

目安値: フローサイトメトリーArg.150以上、CytoFLEX Arg.3000、マスサイトメトリーArg.5 ⑤ すべての蛍光チャネル(解析に使用する)について調整をします。

3-3-2ゲーティングによる解析対象の絞り込み(前処理)

死細胞やダブレットが含まれていると解析に影響するため、解析を始める前に<u>ゲーティング</u>で除いておきま す。とトでは見分けにくい機器測定時の異常イベントを<u>データ QC</u>で自動検出・除去しておくと更に良いで す。また、特定の細胞集団内を解析対象としたい場合は、ゲーティングしてポピュレーションを作成しておき ます。Gating Editor でゲーティング完了後、<u>Apply gates ボタンを押してゲートポピュレーションの情報</u> を Experiment 全体に反映するのを忘れずに行ってください。

3-3-3設定と実行

- ① メニューバー > Advanced analyses > Dimensionality Reduction > New analysis を 選択します。
- ② 実行したい次元削減アルゴリズムを選択し名前を入力します。
- 3-3-3-1 Basic settings (共通基本設定)
- Population でインプットする対象の細胞集団を指定し、FCS Files でファイルを選択します。 (viSNE は population を複数選択することができ、選択した各 population についてファイルを 選択します。)
- ② Event sampling でサンプリング方法を選択します。Desired total events はインプットする細胞数を表示します。各アルゴリズムでの上限イベント数を下に示します。

Num Events	<u>tSNE</u> -CUDA Max Num Channels	UMAP Max Num Channels	opt-SNE Max Num Channels	<u>ViSNE</u> Max Num Channels
9-10M	25	15		
8-9M	50	25		
7-8M	75	35		
6-7M	100	40		
5-6M	100	50		
4-5M	200	70		
3-4M	200	100		
2-3M	200	100	30	
1.5-2M	200	200	60	Enterprise (*)
1-1.5M	200	200	100	Up to 1.3M Premium (*)
250K-1M	200	200	200	*
Up to 250K	250-800	250-800	250-800	*

- Proportional:各ファイルの細胞数/全ファイルの総細胞数を反映してサンプリングする細胞数 を決めます。各ファイルの割合を Desired total Events に反映して各ファイルからのサンプリ ング数を定めます。

selected		
Population - FCS File	Event Count	Proportional Sampling
CD3+ T cells		
pbmc_lrs005_il10.fcs - IL10	170788 39.4%	8658 39.4%
pbmc_lrs005_unstim1.fcs - Unstim1	177718 40.1%	9010 40.1%
CD20+ B cells		
pbmc_lrs005_il10.fcs - IL10	42323 9.8%	2146 9.8%
pbmc_lrs005_unstim1.fcs - Unstim1	43120 9.9%	2186 9.9%
intact cells		
CD3+CD4+ T cells		
CD3+CD4- T cells		
CD33- lymphocytes		
CD33+ monocytes		
Total Events Selected	: 433949	22000 to sample
t Sampling 🕕		
Proportional Equal		

 Equal:各ファイルから同数の細胞をサンプリングします。Desired per file Events (t SNE-DUDA, UMAP, opt-SNE) または Desired total Events (viSNE)を元に各ファ イルから同数をサンプリングします。細胞数の少ないファイルがある場合はその数に合わせて同 数を他のファイルからサンプリングします。

Population ①	FCS files				(3
Choose					Choo	ose
singlets	Selected files		Total events	singlets	To sample	
	Tube1		39.125	36.457	25.000	
CD14+CD16lo	Tube2		28.084	25.477	25.000	
CD4+	Tube3		32.379	30.419	25.000	
CD45+CD3+	Tube4		47.851	43.631	25.000	
CD8+	Tube5		27.531	26.526	25.000	
monocytes	Tube6		30.619	28.985	25.000	
Ungated	Tube7		81.245	76.947	25.000	
	Tube8		42.073	39.878	25.000	
		Total:	328.907	308.320	200.000	
9 present		R of R files selected - Samplin	vg 200 000 tota	levents		
opresent		o or o mes selected — sampin	15 200.000 tota	il evento		_
Event sampling 🚯						
Equal Proportional	esired events per file: 2500	0 × 8 files = 200.000 des	ired events			
Use all events	ctual events to sample: 25.0	000 × 8 files = 200.000				

- All events: すべての細胞を用います。
- ③ Channels でサンプルの細胞集団サブセットやグループ分けに用いるパラメータを選択します。
 - 一般的に除外すべきパラメータ:
 - FSC, SSC
 - ・ 核 (DNA) 染色
 - 生死判定

- ・ ビーズ
- Time
- ・ ハウスキーピング遺伝子
- · 外因性制御遺伝子
- ・ (プレゲートをしている場合)プレゲートとその上流のパラメータ
- ・ (他のアルゴリズムと組み合わせている場合)結果として付与されたパラメータ
- すべてのファイルに共通する Channel のみ選択できます。ファイルのパネルが1種類でない場合、パネル内で解析に用いるすべてのファイルで共通している Channel のみ選択できます。
- ④ 必要な場合、Advanced settings を変更します。Advanced settings については次のセクションに記載します。
- ⑤ (viSNE 以外で Advanced settings の下) Normalizing scales が必要な場合はチェックを 入れます。ここでのノーマライズは各チャネルの平均を0、標準偏差1で規格化したものとなります。 フローサイトメトリーデータでチャネル間のダイナミックレンジが異なる場合に有効なことがあります。
- 6 Data scaling と compensation は事前に確認・調整してあることをお薦めします。
- ⑦ Run Analysis ボタンを押して実行します。
- ⑧ 画面上部にステータスバーが表示されたらサーバーでの計算が始まっています。端末では Cytobankを終了しても大丈夫です。
- ⑨ 完了すると通知のメールが登録のメールアドレスに届きますので結果の URL をクリックします。



- 3-3-3-2 Advanced settings (アルゴリズム別)
 - tSNE-CUDA

tSNEを GPU で高速処理するため、最も高速に次元削減することができます。Cytobank で は以下の各種パラメータ設定(Advanced settings)を自身で調整することも、自動的に 調整させることもできます。

•

Itera	tions 🛈	
750		
Ζ Αι	utomatic 🛈	
Perp	lexity 🛈	
30)
Lear	ning rate	e (i)
166	66	
Ζ Αι	utomatic	
Early	/ exagge	ration 🛈
12		

 Iterations は tSNE アルゴリズムのグラジエント演算の繰り返し計算数です。収束すると 計算を終了する opt-SNE などとは異なり、設定した iteration だけ計算を実行します。 Automatic ボックスのチェックを入れると、対象イベント数を 1500 で除算した数(最小 750) に自動調整されます。iteration 数が小さすぎると細胞集団やサブポピュレーショ ンが tSNE-CUDA マップ上に十分反映されません。収束していない tSNE-CUDA マップ の特徴は下図のようなフィラメント状や引き伸ばしたような状態で見られることがあります。 このような場合は、Automatic ボックスのチェックを外し、iteration 数に任意の数を入 力して、iteration 数を増加します。



Perplexityは、高次元での各イベントの近接点との類似性の計算の際に、近接する何 点について考慮するかを設定します。デフォルトは 30 で、ほとんどのシングルセルデータセッ



トに適していますが、大規模なデータセットの場合は検討する必要がある場合があります。 概ね 5 ~50 での検討をお薦めします。 Learning rate は、機械学習の最適化においてどのくらい値を動かすかの設定をします。Automatic ボックスにチェックを入れると、対象イベント数を early exaggeration (後述)で除算した数(最小 200)に自動調整します。基本的にはデフォルトの自動 調整をお使いいただくのをお薦めしますが、手動調整も可能です。Learning rate は小 さすぎても大きすぎても細胞集団が適切に分離したマップになりません。また、iteration とも密接に関係しているため、learning rate を下げる時は iterations を増やして検討 することをお薦めします。



- Early exaggeration は、グラジエント演算の初期段階において、類似性の計算に系 数を乗算することで、イベントの動きを増加させ、グローバル構造的な主要な分布パターン を先に反映させるとともにクラスター(ランド)の分離を向上させるために設定します。
 Early exaggeration のフェーズの後は、局所構造の反映のみを実行するようになりま す。初期段階の計算において、ギャップが増加傾向を示す場合は early exaggeration か learning rate が大き過ぎる可能性があります。
- UMAP

tSNEと同様に、高次元情報の各イベントの類似性(サイトメトリーデータの場合はマーカー発 現パターン)の分布構造を、低次元(2次元)上に反映して表示(可視化)することがで きます。経験的に、UMAPの方がtSNEよりもグローバル構造が反映されたマッピングが可能 です。3つのパラメータを設定することができます。

Num neighbors は最も重要なパラメータで、初期の高次元グラフを構築する時に用い る最近傍のイベントの数を設定します。 UMAP の局所構造とグローバル構造のバランスに 寄与します。 値が小さいと局所構造をより維持したマップを作成します。 値が大きいとグロ ーバル構造をより維持するマップを作成しますが、 詳細構造が失われる可能性もありま す。 デフォルトは 15 で、 範囲は 2 ~100 です。



Minimum distance は、UMAP マップに表示するイベント間の距離を設定できます。
 値が小さいと密度が高いように表示し、値が大きいとイベント間距離を緩めて表示できます。
 デフォルトは 0.01 で、範囲は 0 ~0.99 です。



Collapse outliers は、外れ値のイベントをまとめることができます。Z スコアが3より大きい場合は外れ値と見なされ、チェックが入っている場合には自動的に最大値または最小値にまとめられます。ほとんどのデータが狭い範囲に絞られているような時には、チェックを外して検討すると良い場合があります。

- opt-SNE

tSNE アルゴリズムのグラジエント演算時、計算の収束を自動的に判断して終了します。大規 模データセットでの解析の際に、結果の品質は維持しながらパラメータの最適化をするための試 行錯誤にかかる時間を節約することができる可能性があります。

Advanced settings 🚯
Perplexity 🕕
30
Learning rate i
16666
Automatic 🛈
Early exaggeration
12
Seed
(Random)
Max iterations 🔅
1000

・ Perplexity は、高次元での各イベントの近接点との類似性の計算の際に、近接する何 点について考慮するかを設定します。デフォルトは 30 で、ほとんどのシングルセルデータセッ



トに適していますが、大規模なデータセットの場合は検討する必要がある場合があります。 概ね5~50での検討をお薦めします。

 Learning rate は、機械学習の最適化においてどのくらい値を動かすかの設定をします。Automatic ボックスにチェックを入れると、対象イベント数を early exaggeration (後述)で除算した数(最小 200)に自動調整します。基本的にはデフォルトの自動 調整をお使いいただくのをお薦めしますが、手動調整も可能です。Learning rate は小 さすぎても大きすぎても細胞集団が適切に分離したマップになりません。また、iteration とも密接に関係しているため、learning rate を下げる時は iterations を増やして検討 することをお薦めします。



- Early exaggeration は、グラジエント演算の初期段階において、類似性の計算に系数を乗算することで、イベントの動きを増加させ、グローバル構造的な主要な分布パターンを先に反映させるとともにクラスター(ランド)の分離を向上させるために設定します。
 Early exaggeration のフェーズの後は、局所構造の反映のみを実行するようになります。初期段階の計算において、ギャップが増加傾向を示す場合は early
 exaggeration か learning rate が大き過ぎる可能性があります。
- Seed は tSNE アルゴリズムの 2 次元乱数マップを作成する際に用いる乱数シードを設定することができます。 つまり、同一データ・同一設定・同一 seed で複数回アルゴリズムを実行すると、まったく同一の結果のマップを得ることができるので、ワークフローの検証やアルゴリズムの再現性の確認に用いることが可能です。 実行前に指定しない場合は、 Cytobank がランダムに seed を設定して実行します。 計算が完了して結果が表示されると、 どの seed が使用されたのかを結果 Experiment の setting 画面で確認することができます。 Seed は tSNE アルゴリズムの客観性の基盤となるパラメータのため、以下 2点にご注意ください。

- 1. 特定の seed をいつも(または複数の Experiment にまたがって)使用すること。tSNE アルゴリズムのランダム性に人為的に制限をかけることになります。
- Seed を入力して実行する場合は、無作為の数値を設定するようにします。
 その時のお財布の中の金額やレシートの数値などが時々使われます。
- Max iterations は、tSNE アルゴリズムのグラジエント演算が収束しなかった場合に終 了させる繰り返し計算回数を設定します。opt-SNE では、Max iterations よりも少な い計算回数でも、収束に至ると最適化されたと判断して自動的に計算を終了します。
- viSNE

tSNE に Barnes-Hut を実装しており、Cytobank の次元削減アルゴリズムではもっとも早く に搭載されています。Premium Cytobank のトライアルユーザーでは、Advanced settings はデフォルトで変更ができない仕様となっています。

Advanced Settings (i)	۰
Seed	(Random)
# Iterations	1000
Perplexity	30
Theta	0.5

- Seed は tSNE アルゴリズムの 2 次元乱数マップを作成する際に用いる乱数シードを設 定することができます。 つまり、同一データ・同一設定・同一 seed で複数回アルゴリズム を実行すると、まったく同一の結果のマップを得ることができるので、ワークフローの検証やア ルゴリズムの再現性の確認に用いることが可能です。実行前に指定しない場合は、 Cytobank がランダムに seed を設定して実行します。計算が完了して結果が表示され ると、 どの seed が使用されたのかを結果 Experiment の setting 画面で確認すること ができます。 Seed は tSNE アルゴリズムの客観性の基盤となるパラメータのため、以下 2 点にご注意ください。
 - 1. 特定の seed をいつも(または複数の Experiment にまたがって)使用する こと。tSNE アルゴリズムのランダム性に人為的に制限をかけることになります。
 - 2. Seed を入力して実行する場合は、無作為の数値を設定するようにします。その時のお財布の中の金額やレシートの数値などが時々使われます。
- # iterations (number of iterations)は、tSNE アルゴリズムのグラジエント演算の 繰り返し計算回数で、デフォルトでは 1000 に設定されています。 データの種類、サンプ

ル種類、パネルや細胞数によってはデフォルトでは計算が十分でない場合があります。フロ ーサイトメトリーデータでは細胞数が10万細胞あたり#iterations1000を検討の目安 となります。マスサイトメトリーの場合はもう少し少ない回数が目安となりますが、パネル

(チャネル数)も影響することにご留意ください。Iterations が小さ過ぎると、下図のよう にボール状やフィラメント状、引き伸ばされたようなマップが作成されます。また、tSNE1, tSNE2 の各軸の範囲が 30 以下の場合も iteration 不足の目安となります。計算の 収束は tSNE アルゴリズムの乱数マップの初期値にも依存するため、十分な iterations も同様であることにご留意ください。



Perplexity は、高次元での各イベントの近接点との類似性の計算の際に、近接する何 点について考慮するかを設定します。デフォルトは 30 で、イベント数 100 以上であれば ほとんどのデータセットで良好な結果となります。設定できる値の上限はイベント数に依存 し、(解析対象イベント数-1)/3よりも大きい値は入力できません。一般的に Perplexity を大きくすると viSNE マップの空間的な分離が向上することが期待されます が、データセット内に分離できる細胞集団が十分ない場合は、perplexity を大きくしても ランドが形成されることはありません。 •



Theta は、Barnes-Hutの影響度の大きさを設定し、基本的にはデフォルトで変更しないことをお薦めします。値が小さいほど tSNE オリジナルに近く計算に時間がかかり、値が大きいほど Barnes-Hutの影響が強くなり計算は早く終了します。



3-3-4複数のアルゴリズム結果の比較表示

次元削減アルゴリズムが完了すると、その結果は元の Experiment に紐づいた result Experiment として Cytobank に存在します。解析前の Experiment(親 Experiment)はメニューバーの帯が水 色ですが、 結果の Experiment は紫色です。

同一の親から複数のアルゴリズムを実行し、結果 Experiment が並列に子となっている場合、どの結果 が最も良さそうかを並べて比較することができます。

① 結果 Experiment の Settings ボタンを押します。

U Cytobank UI	MAP analysis					-	UMAP	Sharing
Actions	🌱 Sample tags	🗘 Gates	ភ្លំ Advanced analyses	L Illustrations	Scales	A Compensation	• Settings	Private

- ② Analysis run info が開きます。
- ③ プロット上部の Compare results with other analysis をクリックします。または、右上 黄緑色の View results> Compare with other analyses をクリックします。



- ④ Compare dimensionality reduction analysis results ウィンドウが開きます。
- Select analysis to compare セクションで表示したい結果 Experiment にチェックを入 れます。
- ⑥ 左側メニューで、比較表示するファイルやマーカー、More settingsを開くと Zoom やカラー パレット、表示ドットの設定変更や背景色を変更できます。



⑦ (必要な場合) Analysis run info の下部は実行した設定を確認することができます。

Population	FCS files			U	channels
Singlets	Selected files	Total Events	Singlets	To Sample	110-CD3
	Marrow1 01 Basal1	351,619	189,316	50,000	115-CD45
Unselected population:	nselected population: Marrow1_06_BCR 472,082			50,000	139-CD45RA
CD123+ DCs	Marrow1_12_IL7	507,989	275,521	50,000	142-CD19
CD19+	Та	otal: 1,331,690	732,509	150,000	144-CD11b
CD3+T cells					145-CD4
CD33+ Monocytes CD4+ T cells					146-CD8
CD45+					147-CD20
CD8+T cells					148-CD34
Memory CD4+ T cells					158-CD33
Naive CD4+ T cells					160-CD123
Naive CD8+ T cells					167-CD38
Ungated					170-CD90
14 present	3 of 3 files selected — Sam	pling 150,000 total e	vents		13 of 41 selected
Advanced settings Num neighbors ① 15 Minimum distance ③ 0.01 Collapse outliers: No ④					
Transformations Data scaling Data will be scaled accordin	ng to the scale settings of the experiment. Edit scales 🗹				
No U					
Compensation					
The experiment compensation will be used: File-Internal Compensation Edit					

3-3-5結果のマップ上でのゲーティング

結果の Experiment では、解析に用いたイベントについて元の Experiment の情報をすべて有したまま、チャネルパラメータに次元削減パラメータ(tSNE1とtSNE2/UMAP1とUMAP2)が追加されています。



Gating Editor の X・Y 軸パラメータを tSNE1・2/UMAP1・2 にすることで、tSNE・UMAP マップとして 表示し、ゲーティングすることが可能です。

- ① メニューバーの Gating をクリックします。
- ② Active Population でプロットに表示したい集団を選択します。
- ③ Plot type を Dot-Color by channel にします。
- ④ X 軸を tSNE1/UMAP1 に設定します。
- ⑤ Y 軸を tSNE2/UMAP2 に設定します。
- ⑥ Coloring Channel をカラースケールの色で表示したいパラメータに切り替えます。
- ⑦ パラメータを細胞集団を同定するマーカーに切り替えると、プロット上のどの細胞がそのマーカーを強く発現しているかを色で確認することができます。
- ⑧ どのランド(プロット上の島)がどの細胞集団であるかが判断できたら、ゲーティングツールで細胞 集団を囲い、名前を付けます。
- ③ マップ上でゲーティングした集団も、マニュアルゲーティングの集団と同様ゲートリストに追加されます。
- ⑩ Apply gates ボタンを押します。
- Illustration に移動しますので、マップ上でゲーティングしたポピュレーションを用いて、サンプル間 比較やヒートマップ表示などを行うことができます。



3-3-6結果の良い/悪いの見方

機械学習処理である次元削減アルゴリズムでは、インプットするデータが適切な前処理をされていなかったり、設定が適切でなかったりすると良くない結果となります。

良くない例としては、たとえば viSNE では下図左のように奥行がありランドに分かれていないプロットが出 力されます。



このような結果が得られた場合は以下をご確認ください。

- 設定した細胞集団は適切か(デブリやダブレットを含んでいないか)。
- 使用したパラメータのスケール設定は適切か。
- コンペンセーションは適切か。

上記に問題ない場合は、演算不足の可能性が考えられますため、viSNE 設定の Advanced setting を開いて、Iterations をデフォルトの 1000 から 3000 または 5000 に増やしてお試しください。



他のアルゴリズムについては、デフォルト設定でほとんどの典型的なデータセットに良好なようになっていますが、Advanced settings もご参照ください。

3-3-7結果の illustrations での表示と比較

特定のマーカーの発現強度のサンプル間での違いを、次元削減アルゴリズムのマップで比較したい場合、 illustrations でプロットを並べることで分かりやすい Figure にすることができます。 ここでは例として、viSNE 結果の刺激なしと LPS 刺激後のサンプルとの間で、CD33+の細胞集団にお いて pp38 リン酸化強度が異なっている figure の illustrations 設定を展開します。 背景:刺激なしサンプルと LPS 刺激ありサンプルについて、細胞集団を同定するマーカー CD3,4,8,20,33 を Channels に設定し、viSNE を実行しました。viSNE 完了のメールを受け取り、 そのリンクを開きました。

3-3-7-1 Channel colored map でサンプル間の発現変動を比較

① Gating を開きます。



- ② Coloring Channel を切り替えて各マーカーの色強度から、どのランドがどの細胞集団かを判断し(または対象の集団のみを)ゲーティングします。
- ③ Apply gates を押します。
- ④ メニューバー> illustration > New illustration で illustration Editor に移動します。
- ⑤ Plots と Layout を開き、下図のように設定します。
- ⑥ Layout の各項目は数字 of 数字部分(上図赤枠)をクリックして以下に設定します。
- ⑦ FCS files では、刺激なし(Unstim1)と LPS 刺激あり(LPS)のサンプルファイルを選択します。
- ⑧ Channels では、CD33とpp38のみを選択します。
- ⑨ Populations では、singlets(プロットに表示したい細胞集団)を選択します。
- 10 プロットが自動で変更されます。
- ① CD33+のランド(②のゲーティングでは viSNE CD33++) において、刺激なしのサンプルでは pp38 は上がっていないが、LPS 刺激ありのサンプルでは pp38 が上がっていることが示されてい ます(赤矢印)。

(i) Plots	×	(i) Layout			×
Plot type	Dot	Rows 1	FCS Files	2 of 8	
Overlaid		Columns 👄	Channels Z	2 of 24	
Color by	Z-axis channel	Table 1 ↔	☆ Populations	1 of 9	
Zaxis	Varying by Columns				
Palette	Rainbow				
Plotting method	Binned Pixels				
Show gates					
Size	Medium				
🗹 Mo	ore plot settings		🗹 Layout Preview		
ation					
	Unstim1	Single cells 2: p38 4564 -1302 -359 -54 -153 -153 -153 -154 -153 -154 -155 -154 -154 -155 -154 -155 -154 -155 -1	1463 -590 -222 -44 -97 1463 -590 -222 -44 -44 -97		

② Plots 上で Show gates を ON にするとプロットに②のゲートを表示することもできます。

表示するラベルは More plot settings をクリックして Gate label 設定で、表示なし(None)、イベント数(Event count)、ゲート名(Gate Name)、パーセント(Percentages)を選択することができます。



Overlaid や <u>Concatenate</u> で <u>Tailored par file ゲーティング</u>をしている場合は、最初のファイルのゲート位置が表示されます(統計値等は Tailored の位置で計算されます)。

3-3-7-2各細胞集団のヒートマップ表示で比較

- ① illustrations > New illustration \tilde{c} illustration Editor へ移動します。
- ② Plots と Layout を下図のように設定します。

(i) Plots	×	(i) Layout		×
Plot type 🔛 Heatmap		Rows 1	☆ Populations	4 of 13
Statistic Median		Columns \leftrightarrow	🕌 Channels 🛛	6 of 24
Scale range Global		Table 1 ↔	FCS Files	2 of 8
Palette Yellow			Add dimension	
 Viewthrough plot 				
☑ More plot settings			🗹 Layout Preview	

- ③ Layout の各項目は数字 of 数字部分(上図赤枠)をクリックして以下に設定します。
- ④ Populations では作成したゲートからヒートマップに使用したいポピュレーションを選択します。ここでは viSNE マップ上で作成したゲートとします。
- 5 Channels では細胞表面マーカーと機能性マーカーをすべて選択します。

 ⑥ FCS files では刺激なしサンプル(Unstim1)と LPS 刺激ありサンプル(LPS)のファイルを選 択します。



- ⑦ ヒートマップが自動変更され、表示されます。
- 8 LPS 刺激サンプルで、viSNE マップ上で作成した CD33+のポピュレーションでのリン酸化機能性 マーカーpp38の発現が増加していることが示されています。
- ⑨ ヒートマップ上にオーバーカーソルをするとポップアップが表れ、セル内イベントのプロットが表示されます。
 Viewthrough plot 部分でポップアッププロットの設定を変更することもできます。
- ・ ヒートマップに用いる統計値は Plots の Statistics で変更するか、ヒートマップ下部の Statistics 欄 Displaying 部分で変更可能です。
- 3-3-7-3マニュアルゲーティングを結果のマップ上に表示
 - ① Gating Editor でマニュアルゲートをしておきます。
 - ② illustrations > New illustration で illustration Editor へ移動します。
 - ③ Plots と Layout を下図のように設定します。

④ Layout 設定の Populations で①のマニュアルゲートポピュレーションを選択します。



- ⑤ プロットの X 軸 Y 軸が tSNE1 & tSNE2 でない場合はプロットをクリックし、ポップアップの Setting で X axis と Y axis を変更します。
- ⑥ 凡例の名前をドラッグ&ドロップで上下を変更することができます。凡例の一番上にある項目がオ ーバーレイの最下層に配置され、並び順に配されています。

3-3-7-4ファイル統合プロットでマップの全体像を表示する 次元削減マップのファイル統合プロット表示をご参考ください。

3-4 FlowSOM

3-4-1 FlowSOM とは

処理が速く質の高いクラスタリングができる Self-Organizing Maps (SOM s)アルゴリズムを用いて各 マーカーが全細胞でどのように振舞っているかを可視化する手法です。設定されたパラメータをもとに細胞 をクラスター化した後、クラスターの SOM を作成し、これを Minimum Spanning Tree(MST)にしま す。MST 上のクラスターには設定された数のグループに自動的に分けられメタクラスターとして割り当てられ ます。Cytobank では、FlowSOM 結果として、ポピュレーション細胞数とマーカー強度についてをパイチャ ート、スターチャートやチャネルカラープロットなどの形式の SOMs と MTSs で出力するほか、新規 Experiment として作成し、元のチャネルにクラスターとメタクラスターの ID を付加します。

3-4-2tSNE/UMAP マップ上に FlowSOM メタクラスターを表示する場合

FlowSOM で作成されたメタクラスターを tSNE/UMAP マップ上に表示したい場合は、先に tSNE/UMAP を施しておきます。流れとしては、データのインポート> コンペンセーション、スケール調整、 データ QC、クリーンアップ> (プレゲート)> tSNE/UMAP>FlowSOM となります。この時、 FlowSOM は tSNE/UMAP の結果 Experiment(帯が紫色)から始めてください。

3-4-3 FlowSOM 設定·実行

- ① メニューバー> Advanced Analyses> Clustering > FlowSOM > New Analysis を選択 します。
- ② New FlowSOM analysis name;に解析名を入力します。
- ③ Population で、対象の細胞集団(ゲートポピュレーション)を選択します。
- ④ FCS Files で、ファイルを選択します。
- ⑤ Clustering channels でクラスタリングに用いるチャネルを選択します。

G FIOWSOFF Analysis Set	cup Copy Settings	Status: New	Run FlowSOM Ana	lysis
Analysis Name: FlowSOM de	FCS Files			ViSNE 解析名の編集
Ungated Unselected population: (2024) (2024) (2024): 8 cell 解析対象集 団(ゲート 名)を選択	Unselected files 11.10 11.6 LPS Unstim1 Unstim2	Total Eve 8,315 8,272 8,407 8,402 8,409 ファイルを選択	nts Ungated 8,315 8,222 8,407 8,402 8,409	- First select some files クラスタリング チャネルを選 択
17 present		0 of 5 files selected		

- ⑥ Event Sampling で、データのサンプリング方法を選択します。
 - Equal: 各ファイルの細胞数の違いは無視して同数の細胞をサンプリングします。
 - Proportional: 各ファイルの細胞数の違いの比を反映してサンプリングする細胞数を決め ます。
 - Use all events:すべての細胞を用いる(上限 400 万細胞)。

Algorithm settings

⑦ SOM creation で、用いる SOM を選択します。

- Create a new SOM:新規に SOM を作成します。
- Use an existing SOM from another run:既存の SOM を用います。選択すると使用 する SOM の Experiment の FlowSOM を指定できるウィンドウが表示されます。
- ⑧ Clustering method で、クラスタリング方法を選択します。Hierarchical Consensus を推 奨します。
- ⑨ Number metaclusters で、メタクラスターの数を設定します。
- ⑩ Number clusters で、クラスターの数を設定します。
- ⑪ Iterations で、機械学習の繰り返し数を設定します。10を推奨します。
- Seed で、乱数シードの設定をします。Random を推奨します。

Transformations

- ③ Data scaling で、スケールノーマライズの有無を設定します。
- ⑭ Compensation で、適用されているコンペンセーションを確認できます。

🚍 Actions 🛛 🔟 Illustrations 🛛 🌱 Sample Tags 🛛 📅 Advanced Analyses 🔹 🟠 Gating	
FlowSOM Analysis Setup Copy Settings	5
Algorithm settings	
SOM Creation ③	用いる SOM の選択(新規/既存)
Clustering method ()	クラスタリング方法:Hierarchical Consensus 推奨
k-Means Number metaclusters () 10	メタクラスター数
Number clusters () 100 •	クラスター数
Iterations ① 10	機械学習の繰り返し数:10 推奨
Seed () (Random)	乱数シードの設定 : Random 推奨
Transformations	
Data Scaling () FlowSOM scales data according to the scale settings of the experiment. Edit scales (Normalize Scales	スケール設定
Compensation ① File-Internal Compensation ①	コンペンセーション設定(表示のみ)

PDF output settings

ID Populations for pie charts で、PDF 出力するパイチャートに表示する細胞集団を選択します。

- ① Clustering sizing で、クラスターの表示サイズを選択します。
 - Relative: クラスターの細胞数に応じて相対的に変化させます。
 - Fixed:固定
 - Relative and Fixed:両方
- 18 Output file type で、出力するファイル形式を選択します。PNG の方が解像度は高いです。
- 19 Toggle metacluster background in plots で、メタクラスター表示するプロットを選択しま

	Run HowSow Analysis
PDF output settings	
	バイチャート表示する集団設定
Choose Choose	
CD20+ PacOrange-A	
CD3-CD20+:B cell Qdot525-A	
CD3-CD33+:DC etc Qdot605-A Odot655-A	MTS で表示するチャネル設定
CD3-CD33+:Monocyte Qdot705-A	THU CANNY OF THURK
CD3+:Tcell Time	
(CD3+:Coll) (CD3+:Tcell) (SNE1	
16 of 16 selected 6 of 20 selected	J
Cluster sizing (i)	
Relative Max relative cluster size (pixels)	クラフターの表示サイズ
© Fixed 15	
Relative and Fixed	
Relative and Fixed	
Relative and Fixed Output file type ③	出力ファイル形式
Relative and Fixed Output file type ③ PDFs PDFs (another interpreted file also)	出力ファイル形式
Relative and Fixed Output file type ③ PDFs PNGs (greatly increases file size) Reth type:	出力ファイル形式
Relative and Fixed Output file type ① PDFs PNGs (greatly increases file size) Both types	出力ファイル形式
Relative and Fixed Output file type ① PDFs PNGs (greatly increases file size) Both types Toggle metacluster background in plots ①	出力ファイル形式
 Relative and Fixed Output file type () PDFs PNGs (greatly increases file size) Both types Toggle metacluster background in plots () On channel-colored MSTs 	出力ファイル形式 メタクラスタリングするプロットの選択
 Relative and Fixed Output file type ① PDFs PNGs (greatly increases file size) Both types Toggle metacluster background in plots ① On channel-colored MSTs On legend plots 	出力ファイル形式 メタクラスタリングするプロットの選択

- 20 画面上部緑色の Run FlowSOM Analysis ボタンを押して実行します。
- 22 画面上部にステータスバーが表示されたら Cytobank を終了しても大丈夫です。
- ② 計算が完了すると通知のメールが登録のメールアドレスに届きますので URL をクリックします。

	9 1 2 0 5 5 5
	[Premium Cytobank] Your FlowSOM analysis is complete ▷ ₩₩₩₩₩₩
+	admin@cytobank.org 13:48 (43 分前) ☆ ヘ To 自分 マ
	ズA 英語 ◆ > 日本語 ◆ メッセージを翻訳 ^{次の言語で無効にする:}
	Hi Yuko,
	The FlowSOM analysis FlowSOM demo for experiment <u>vISNE_157552_vISN</u> result.(Clone) is now complete
	<u>Click here to view your FlowSOM analysis</u> . The PDF output can be found atta to the analysis settings page or on the Experiment Summary page of the pare experiment.
	- The Cytobank Team

3-4-4 FlowSOM 結果の出力ファイル

- ① 通知メールの URL を開くと FlowSOM 結果画面が開きます。
- 画面上部の Download the run info and plots をクリックし、FlowSOM 結果フォルダをダ ウンロードします。

Analysis complete!	View created experiment	B Download run info and plots

- ③ ダウンロードしたフォルダの中には、FlowSOM 設定にもよりますが、以下のファイルが入っていま す。
- FlowSOM run info file:設定や解析の情報が入っています。
- Supporting_files: 結果の解析には用いません。FlowSOM クラスターをもとに新規 FCS ファ イルを作成したり、解析プロセス時間の記録が含まれています。
- Results フォルダ: PDF や PNG 形式の結果ファイルが入っています(詳細は次のセクション)。
 ファイル形式は FlowSOM 設定 Output file format によって異なります。
- 3-4-4-1 Result フォルダ内結果ファイルの見方

FlowSOM 設定にもよりますが、典型的な results フォルダは以下のファイルを含みます。設定で Output format を PNG にした場合は、 PDF でなく同じ名前ルールで PNG ファイルとなります。

Name	Туре
channel_colored_MSTs	File folder
mst_definition	File folder
🛃 legend	Adobe Acrobat Document
💫 metaclustering_comparisons	Adobe Acrobat Document
Population_pie_charts	Adobe Acrobat Document
ଌ star_plots	Adobe Acrobat Document

- Legend.pdf: クラスターとメタクラスターの MST と SOM フォーマットでの ID を見ることができます。
- Channel_colored _MSTs フォルダ: FlowSOM 設定の PDF output settings で選択したチャネルについてすべてのサンプルを統合した MST を表示する
 aggregated_channel_coloerd_MSTs.pdf を含みます。メタクラスターの表示の有無は設定の toggle the meta-cluster background で変更可能です。
- Metaclustering_comparisons.pdf: hierarchical consensus clustering・ hierarchical clustering・k-Meansの3つのクラスタリング方法によるメタクラスターMSTが入 っており、最適化のための比較ができます。設定で指定したものについてはラベルに記載がありま す。

Star_plots.pdf:スタープロットでのMSTとSOMフォーマットによる結果です。スタープロットでは、クラスター内のすべての細胞についての、(FlowSOM設定で指定した)clustering markerの平均強度が示されています。扇形になっている各マーカーの高さが強度を示しており、 扇が大きいものほど発現強度が高いということになります。クラスタリングチャネルとスタープロットで 表示しるチャネルは別々に設定できるため、Seedを同一にして再実行することで、見せたいチャ ネルを強調することができます。



 Population_pie_charts.pdf:マニュアルゲーティングとの比較ができます。パイチャートでは、 FlowSOM が自動的に作成したクラスター内に、マニュアルゲーティングでの各ポピュレーションが 何%含まれているかを示します。MST と SOM フォーマットの両方が出力されます。クラスターの周 りを囲む色で示されているメタクラスター表示の有無は設定の toggle the meta-cluster background で変更可能です。



- mst_definition フォルダ:MST の構築に関連する情報が入っています。

※csv ファイルの出力はなくなりました。FlowSOM 結果 Experiment から <u>Export statistics</u> にて出 カください(テンプレートを使用すると便利です)。

3-4-5 FlowSOM 結果の Cytobank への出力

FlowSOM が完了すると、元の Experiment の情報に FlowSOM cluster ID と metacluster ID (下図) が追加された FlowSOM 結果 experiment が作成され元 Experiment の子 Experiment としてリンクされます。



FlowSOM 結果 Experiment では、metacluster_id を番号ごとに自動ゲーティングし(上図右)、 ポピュレーションとします。また、tSNE などの次元削減アルゴリズムの結果 Experiment から FlowSOM を実行すると、illustrations に Metacluster Box plot、Metacluster Dot overlays、 Metacluster Heatmap の3つの illustration が自動的に作成されます。

3-4-5-1 クラスターの自動ゲーティング: Cluster gates

FlowSOM メタクラスターは自動でポピュレーションとして作成されますが、FlowSOM クラスターや他のア ルゴリズムのクラスターパラメータや、または任意の複数のクラスターを ID 番号で自動的にポピュレーション 化することができます。その他 file number などのナンバリングパラメータでも実行可能です(一度に作 成できる数は 200 までなので、それ以上の場合は複数回に分けて実行します)。

- ① メニューバー> Gates > Gating Editor をクリックします。
- ② Gating 画面上部の Cluster gates をクリックします。

✓ © Gating Editor
 ☑ Apply gates
 ☑ Ø ☆ + +- +- ☆ Import gates
 ☆ Cluster gates
 ♡ Reset all

- ③ Add Cluster Gates ウィンドウが表示されます。
- ④ Select a channel でクラスターチャネル (例 FlowSOM_cluster_id) を選択します。

- ⑤ Gate name prefix: に名前を入力します。
- ⑥ Define specific gates: では必要であれば以下のような設定ができます。
 - 特定の ID のクラスターだけゲートポピュレーションにしたい場合は、ID 番号をコンマ区切りで 入力します(例クラスターID 1と2と3を各々ゲートポピュレーションにしたい→1,2,3)。
 - 連続した番号を指定する場合は最初の番号と最後の番号をコロンで繋ぎます(例: クラス ターID 5 から 10 までを各々ゲートポピュレーションにしたい→5:10
 - クラスターを統合してゲートポピュレーションにしたい場合は()でくくります(例:クラスターID
 6と8と10を統合したい→(6,8,10)、20から40までをすべて統合したい→
 (20:40)、クラスターID1から3までと10から12までを1つに統合したい→
 (1:3,10:12))。
- ⑦ Apply Change ボタンを一回だけ押します。ここでは画面は変わりません。

Select the channel that contains the cluster IDs: Select a channel Gate name prefix: cluster_ Define specific gates (optional): (1) i.e. 1,2,3,5:10, (6,8,10), (20:40)	$\hat{\Omega}$	Add Cluster Gates	×
Gate name prefix: cluster_ Define specific gates (optional): (1) i.e. 1,2,3,5:10, (6,8,10), (20:40)	Selec	t the channel that contains the cluster IDs: ect a channel •	
Cluster_ Define specific gates (optional): (i) i.e. 1,2,3,5:10, (6,8,10), (20:40)	Gate	name prefix:	
Define specific gates (optional): (i) i.e. 1,2,3,5:10, (6,8,10), (20:40)	clust	er_	
i.e. 1,2,3,5:10, (6,8,10), (20:40)	Defin	e specific gates (optional): 🛈	
	i.e. 1	,2,3,5:10, (6,8,10), (20:40)	
		Cancel Apply changes	5

- ③ ブラウザの再読み込みボタンを押します。Gates リストに作成されたクラスターゲートが表示されます。
- ⑨ Gating 画面の Apply gates ボタンを押します。

3-4-5-2 FlowSOM 結果の評価

FlowSOM 結果の評価方法としては、以下の3つが便利です。

- FlowSOM メタクラスターをtSNE マップ上にオーバーレイ表示
 - FlowSOM を実行する前に tSNE 等の次元削減アルゴリズムを実施し、その結果 Experiment (紫色帯)の Advanced analysis から FlowSOM を実行します。 FlowSOM 結果 Experiment を開きます。
 - メニューバー> Illustrations > View all saved をクリックし Metacluster Dot Overlays を開きます。

③ Plots と Layout を下図のように設定します。

Plots		>	×	(i) Layout			
Plot type	Dot			Overlaid	Ĵ	Ŷ Populations	6 of 18
Overlaid	\checkmark			Columns	\leftrightarrow	FCS Files	2 of 5
Color by	Overlaid dimension					Add dimension	
Dot size	2	\$					
Dot shape	Circle						
Dot opacity %	100	\$					
Show gates	\bigcirc						
Size	Medium						
🗹 Mo	ore plot settings					Layout Preview	

- Plot type : Dot
- Overlaid : On
- Color by : Overlaid dimension
- Layout 設定の Overlaid: Populations
- Columns : FCS files
- ④ Layout の Populations で FlowSOM メタクラスターのポピュレーションを設定します。
- ⑤ プロットの X 軸 Y 軸が tSNE1 & tSNE2 でない場合はプロットをクリックし、ポップアップの Setting で X axis と Y axis を変更します。

Illustration	olot settings		Unstim1 - pbmc_lrs005_unstin _viSNE_FlowSOM	n1_Single_cells	Export
	FlowSOM_metacluster1 FlowSOM_metacluster2 FlowSOM_metacluster3 FlowSOM_metacluster4 FlowSOM_metacluster5 FlowSOM_metacluster6	Unstim1	Overlaid: Column – FCS File: X axis: tSNE1 Y axis: tSNE2 Plot settings Plot type Overlaid Color by Dot size Dot shape Dot opacity % Show gates	6 Populations Unstim1 C More p Dot Overlaid dimen 2 Circle 100	lot settings sion (\$
Statistics (1) Hide			Size	Medium	

 Save> Clone new copy で複製を別名作成し、Layout の Populations でマニュアルポ ピュレーションのみを選択すると、viSNE マップ上でのマニュアルゲーティングの分布と FlowSOM メタクラスターを比較することができます(下図左:マニュアルゲーティングのポピュレ ーションを viSNE マップ上に表示、右: FlowSOM によるメタクラスター)。この時、ブラウザの タブを分けることでモニターに並べて表示することもできます。



- クラスタリングチャネルでのメタクラスターのヒートマップ表示
 - メニューバー> Illustrations > Metacluster Heatmap を開くか、ない場合は New を選 択し、名前をつけます(例 FlowSOM Heatmap)。
 - ② Plots と Layout を下図のように設定をします。Populations では FlowSOM で作成された メタクラスターのポピュレーションを、Channels では各メタクラスターのフェノタイプを同定するため Clustering channel を設定します。

(i) Plots		×	(i) Layout				×
Plot type	🔛 Heatmap		Rows	\$	1 Populations	6 of 18	
Statistic	Median		Columns	\leftrightarrow	🕌 Channels 🛛	6 of 22	
Scale range	Global		Table 1	↔	FCS Files	2 of 5	
Palette	Yellow				Add dimension		
⊁ Vie	wthrough plot						
🗹 Mo	More plot settings				🗹 Layout Preview		

- Plot type : Heatmap
- Statistic : Median
- Layout O Rows : Populations
- Columns : Channels
- Table 1 : FCS Files
- ③ 細胞表面マーカーの強度の情報から、各メタクラスターがどの細胞集団であるかや、どのような 特性を有する集団であるかを確認し、同定することができます(下図の例では FlowSOM_metacluster1はCD3+CD4+、FlowSOME_metacluster2は CD3+CD4-など)。メタクラスターを統合したい場合は、<u>Automatic cluster gate 機能</u>で 統合したゲートポピュレーションを作成し、そちらを illustration で使用します。



■クラスタリングチャネルでのメタクラスターのヒストグラムやドットプロット表示 Illustrations で、下図のようにヒストグラムやドットプロットで比較や同定も可能です。 Populations では FlowSOM で作成されたメタクラスターのポピュレーションを、Channels では各 メタクラスターのフェノタイプを同定するため Clustering channel を設定します。 ヒストグラム例



ドットプロット例

	 Plots 	×	(i) Layout		>	¢
	Plot type Dot		Overlaid 1	☆ Populations	6 of 18	
	Overlaid		Columns ↔	🛓 Channels 🛛 X	6 of 22	
	Color by Overlaid dimensi	ion	Table 1	FCS Files	2 of 5	
	Dot size 2	•		Add dimension		
	Dot shape Circle					
	Dot opacity % 30	¢				
	Show gates					
	Size Small					
	More plot settings			Layout Preview		
Illustrat	ion					
musua						
				actim 1		
			01	ISUITI		
		CD33 (v) CD20 (v)	CD3 (v)	CD4 (v)	pStat3	pp38
	FlowSOM_metacluster1					
	FlowSOM_metacluster2	13	E	E E		
	FlowSOM_metacluster4	1			a a	
	FlowSOM_metacluster5	***				
	Tionsoni_inclucion	CD33 (v) CD20 (v)	CD3 (v)	CD4 (v)	pStat3	pp38
					Processory.	
				LPS		
		00000000000000000000000000000000000000			1000 1000	
		CD33 (v) CD20 (v)	CD3 (v)	CD4 (v)	pStat3	pp38
	FlowSOM_metacluster1	1 1		1	<u> </u>	
	FlowSOM_metacluster3					
	FlowSOM_metacluster4					
	FlowSOM_metacluster6		1	E		
	η	CD33 (v) CD20 (v)	CD3 (v)	CD4 (v)	pStat3	pp38

- 3-4-5-3グループ間で有意差のある FlowSOM メタクラスターを検討
- <u>箱ひげ図</u>でグループ間に有意差のあるメタクラスターがあるかを検討する
 FlowSOM 結果 Experiment に自動作成された Metacluster Box plot という名前の
 illustrations を開き、Plots 設定と Layout 設定を <u>sample tags</u>を用いて調整して、各メタクラ
 スターについて群間比較と<u>有意差検定</u>し、グループ間で有意差のあるメタクラスターを同定することが
 できます。



3-4-5-4 <u>ヒートマップ</u>でマーカーの変動を可視化する

FlowSOM 結果 Experiment に自動作成された Metacluster Heatmap という名前の illustrations を開き、Clustering channel に設定しなかった機能性マーカー(サイトカインやリン酸 化、遺伝子など)を表示する Channel に設定することで、サンプル間の変動を可視化することができま す。



3 - 4 - 5 - 5 複数のファイルをグループにまとめて FlowSOM メタクラスターを次元削減マップ上にオーバ ーレイ表示する

FlowSOM メタクラスターを次元削減マップ上にオーバーレイで統合表示する機能は <u>virtual</u> <u>concatenate</u>に搭載されていないため、R を使用して FCS ファイルを1つにまとめた統合ファイルを作成 し、それを Cytobank ヘ再アップロードする必要があります。 <u>サポート</u>までお気軽にお問い合わせください。

3-5SPADE

3-5-1 SPADE とは

SPADE (Spanning-tree Progression Analysis of Density-normalized Events) とは、表 現形の類似した細胞をクラスタリングし、階層的なツリー構造で表現する可視化手法です。

【参考文献】

Extracting a cellular hierarchy from high-dimensional cytometry data with SPADE. Qiu P, *et al.* Nat Biotechnol (2011) 29(10):886-91

3-5-2 SPADE 設定・実行

- ① Experiment メニューバー> Advanced analyses>SPADE > New analysis を選択しま す。
- ② New SPADE analysis name: に名前を入力します。
- ③ 画面が切り替わります。

	Actions	1 Illustrations	🌱 Sample Tags	🔶 SPADE	🗳 visne	🖉 CITRU	IS 🏠 Gating	
	< ? 🔶 SPAI	DE Analysis <i>'New</i>	SPADE'					
Target Number of Nodes : ノード数の指定 目安)9~15 チャネルの解析で 200 ノード	SPADE A Compensation File-Internal Co Target Numbe 200 Downsample Percent 10 Absolute Num	nalysis Controls 2	SPADE Setup Settings - click Choose Population 1 of 7 selected Cli Ungated	ie/Gate/Setup to o	Configure Clustering Ch 18 clustering cha Nothing selected choose your Clus Channels. Unselected Clu Channels. ESC-A	annels Choose Setup Click to ening stering	Panel 1	Setup
Down sampled Events Target : ダウンサンプルの設定 Percent : 各ファイルで総イベント	Actions Rename / Copy Ana	inalysis lysis Settings	CD3-CD3+IMoney CD3-CD3+IMoney (CD3+CD4+IHelper (CD3+ICell) When you are d	yte T cell done setting up	FSC-W SSC-A CD3 Time pStat3 PE-TR-A CD20	rameters, click to	p run the analysis.	
数あたり何%をインプットするか。 Absolute Number:インプット細 胞数(各ファイル)		/				実	行ボタン]
デフォルト 10%(推奨)	Populati 解析対象 ピュレーシ	on : のゲートポ ョンの指定	Clusterin Channels クラスタリン チャネルの3	ig s: グに使用 指定	する	Fold-C Group フォール ールとな	hange s: ドチェンジのコン るファイルの指	小口定

- ④ Target Number of Nodes: クラスタリングによって形成されるノードの数を指定します。目安としては 9~15 パラメータのデータセットの場合 200 ノードを設定します。
- ⑤ Downsampled Events Target:ダウンサンプル強度を指定します。ダウンサンプルではヒスト グラムで著しく高いピークの上部を切るような方法で SPADE へのインプット細胞数を減らすととも に、低いピークを持つ細胞集団を同定しやすくします。Percent または Absolute number のど ちらかに値を入力します。Percent は、各ファイルの全細胞数を 100%としたとき、何%の細胞

数にダウンサンプルして SPADE ヘインプットするかを設定します。Absolute number は、各ファ イルから何細胞にダウンサンプルしてインプットに用いるかを設定します。10%をデフォルトとして推 奨しています。

- ⑥ Population:解析に使用するゲートポピュレーションを指定します。通常は死細胞・ダブレット・ デブリ除去後のゲートポピュレーションを指定します。
- ⑦ Clustering Channels: クラスタリングに使用するチャネルを指定します。細胞集団を規定する チャネル(細胞表面マーカーなど)をパラメータとして選択します。
- ⑧ Fold-Change Groups:コントロールとなるファイルを指定します。
- の Compensation: コンペンセーションの調整が可能ですが、先にコンペンセーション調整は済ませておいて File internal を使用することを推奨します。
- 11 画面下部の緑色部分をクリックして実行します。

3-5-3 SPADE 結果の表示と解析

① 通知メールの URL を開くと SPADE ツリーのある結果画面が開きます。



- ② 画面左の機能ツールボックス SPADE Tree Rending を開き、ツリー表示するファイルやパラメー タを切り替えてどのノードがどの細胞集団であるかを解析します。
 - Sample/File Name: 表示するサンプルファイルの選択


- Parameter: ノードの色強度として表示するパラメータ
- Metric: ノードの色強度に用いる統計値
- Scale Range: ノードの色強度の最大/最小値の決定方法
 - ・ Local→現在表示されているファイルのみでの最大/最小値
 - ・ Global→全てのファイルの最大/最小値
- Scale Symmetry: 色強度を基準値から上下に対称とするか(Symmetry)しないか (Asymmetry)
- Color Scheme: 使用するカラースケールの選択
- Background Color: 背景色の選択
- Node Size: ノードの直径の最大/最小値の設定
- Scale Range: ノードの直径の最大/最小値の決定方法
 - ・ Local→現在表示されているファイルのみでの最大/最小値
 - ・ Global→全てのファイルの最大/最小値
- Add Bubble : 選択したノードをまとめてバブルを作成します。
- Remove Selected Bubble(s): 選択したバブルの解除
- Select All nodes: すべてのノードを一括選択します。
- Rotate Selected Nodes: ノードやバブルを選択しボタンを押して回すと任意の角度に
 回転します。
- Save Tree:結果から追加したバブルやツリーの修正を保存します。
- Reset visualization:メール通知されたときの初期の SPADE ツリーに戻します。

3-5-4バブルの作成

- ① 表現形の似たノード同士を囲って細胞集団などの名前をつけることができます。
- ② バブルとして囲いたいノードを複数選択します。選択すると、ノードの縁が太くなります。Shift ボタ ン+クリックでノードごとに選択解除できます。
- ③ Add Bubble ボタンをクリックします。
- ④ バブルの名前(細胞集団名など)を入力し OK を押します。

バブルを選択して Remove selected bubble ボタンをクリックするとバブルを解除することができます。

3-5-5選択したノードのプロット表示

選択したノードまたはバブル内の細胞についてプロットで表示して確認することができます。

- ① ノードもしくはバブルを選択します
- ② 画面左の機能ツール Events Plot を開きます。
- ③ Plot type:プロットの種類を選択します。
- ④ Y-Channel:Y軸のパラメータを選択します。
- ⑤ X-Channel: X 軸のパラメータを選択します。

Pop-Out New Plot をクリックすると、プロット図を新しいウィンドウで表示できます。



3-5-6バブル内の細胞(イベント)の出力

バブル内のイベントデータを出力し、新たな Experiment を作成することができます。SPADE 解析による細胞集団をバブルとして出力し、ドットプロットやヒートマップなどで再解析することができます。

① 画面左機能ツールの Actions を開きます。

- ② Export の Select Bubbles To Export より、出力したいバブルを選択します。
- ③ Export to New Experiment をクリックします。
- ④ バブル内のイベントを新規の Experiment へ出力します。



3-6CITRUS

3-6-1 CITRUS とは

CITRUS (cluster identification, characterization, and regression)とは、相関または予測の アルゴリズムを用い、統計的な群間比較です。群(グループ)間の有意差やバイオマーカー探索を抽出 するのに役立ちます。

CITRUSは2つの工程で構成されています。最初に群の識別なしにすべてのサンプルで SPADE に似た クラスタリングを行います。次に各クラスタリングについて、群間で有意差があるものを相関または予測として 抽出します。抽出に用いるアルゴリズムは SAM、PAMR、LASSO です。

各群3サンプル(n=3)以上で CITRUS 自体の実行は可能ですが、解析結果として用いるには n=8 以上を強く推奨します。

3-6-2CITRUS 設定·実行

- ① デブリ・ダブレット・死細胞除去や対象の細胞集団を絞るゲーティングを前処理として済ませておき ます。
- ② Experiment メニューバー> Advanced analyses>CITRUS > New analysis を選択します。
- ③ New SPADE analysis name:に名前を入力します。
- ④ 画面が切り替わります。
- ⑤ Population ボックスの Choose を押し、対象の細胞集団を選択します。

Actions	🚹 Illustrations	🌱 Sample Tags	🖢 SPADE 🦳 🗳 vis	NE 🖉 CITR	US 🏠 Gating	
∂ CI1	rRUS Analysis Set	Copy Settin	gs Status:	New	Run CITRUS Analysis	
Analys	is Name: test 🧷 💳				解析名を変更	
Popu Singl Unse CD3- CD3- CD3- (CD3 (CD3) Cells Unga	ulation () Choose e cells elected population: CD20+18 Cell CD33+Monocyte :Tcell +:Tcell ted 7 present	File Grouping Cf II (3) unstim (3) Unassigned (1) 6 files in 2 group	Clustering Cluste	choose Setup choose Setup stering	Population:解れ 集団(ゲート名) File Grouping: ループ分け Clustering Chen レーションを規定する 択	斤対象となる細胞 を選択 サンプルファイルをグ anels : サブポピュ るパラメーターを選
File-l	nsation (i)	•			適用コンペンセーショ	コンの選択
Associa	ation Models (j) gnificance Analysis of earest Shrunken Cent -Penalized Regressio	[:] Microarrays (SAM) - roid (PAMR) - Predici n (LASSO via GLMNE	Correlative tive T) - Predictive		使用するモデルを選 ※判断つかない場合 果から検討 ・SAMー相関 ・PAMRー予測 ・LASSO via GLN 群のみ)	訳 合は全部選択し結 METー予測(2

- ⑥ File Grouping ボックスの Choose を押します。
- ⑦ Add file group(s)ボタンを押し、グループ名をカンマ区切りで入力します。
- ⑧ グループ名が表に追加されるので、属するファイルのラジオボタンを ON にします。グループ名とファイル名が部分一致すると自動で割当てられます。間違いがないかも確認します。
- ⑨ Clustering channels ボックスの Choose を押し、クラスタリングに使用するパラメータを選択します。
- Association Models を選択します。
 - SAM-相関
 - PAMR-予測
 - LASSO via GLMNET-予測(2群のみ)
- ⑪ Cluster Characterization で、各クラスターの何をもとに差異を検討するかを選択します。
 - Abundance:存在比率
 - Medians:パラメータ強度(選択するとリストが表示されるので、そこから選択)
- ② Event Sampling で、CITRUS にインプットする細胞のサンプリング方法を選択します。
- ③ Equal:各ファイルから同数(最少のファイルに揃えます)
- ⑭ Max per file: ファイルからの数を揃えない
- ⑤ Event sampled per file で各ファイルから用いる細胞数を入力します。Est. total events to be clustered に総細胞数が表示されます。

ctions 🛛 🗽 Illustrations 🔷 🌱 Sample Tags	🔶 SPADE	🗳 visn	IE 🖉 CITRUS	🏠 Gating	
	Settings	Status: N	ew Rur	CITRUS Analysis	
Cluster Characterization ()	of the experiment.	f Z Edit sca	各クラスターのどち を選択 Abundance:ネ Medians:パラン 解析に寄与する・ Equal:各ファイ ント数に揃える) ・Max per file:	らをもとにして差異を 子在比率 メータ強度(パラメー イベントのサンプリング ルから同数(最少ご 各ファイルの最大数	 -タを選択 ブ方法 ファイルの1 (数を揃
Additional Settings	٥				
Events sampled per file	5000	í	各ファイルから用い	るイベント数を入力)
Est. total events to be clustered	30000		**=+* -*** /-/= 0		
Minimum cluster size (%)	5 (i)	ĥ	坑計七テルに使用	ヨタるクラスターの最近	小イヘント
Est. minimum cluster size	1500	3	交差検証の標本	データ分割数を設定	と(予測せ
Cross Validation Folds	1 (i)	,	レのみ使用)		
False Discovery Rate (%)	1 (i)		海隍性が生じえる	変の期待値を設定	Þ
Normalize Scales	- (j	1	高陽 圧が上しる 低くするとモデルに	開いられるパラメータ	⊑ 9数が制限
Plot Theme	Light •	÷	チェックすると、各 」、標準偏差が1	パラメータ全イベント ⁻ . になるようノーマラィ	での平均を ′ズ

- Minimum cluster size (%)で、CITRUS の最初のステップのクラスタリングでのクラスターの最
 小数(%を変えると下の Ext. minimum cluster size に下図が表示)を設定します。
- ① Cross Validation Folds で、交差検証の標本データ分割数を設定します(予測モデルでのみ 使用)。解析時間に大きく影響するので、検討では1を、論文使用では10を推奨します。
- 18 False Discovery Rate(%)で、FDR 閾値を設定します(PAMR でのみ使用)。
- ⑨ Normalize Scales では、チェックを入れると各パラメータの全イベントでの平均をゼロ、標準偏差が1になるようにノーマライズします。
- 20 Plot Theme で、結果グラフの背景を白か黒にできます。
- 上部の Run CITRUS Analysis ボタンを押します。
- 22 画面上部にステータスバーが表示されたら、Cytobankを終了しても大丈夫です。
- 23 完了すると結果のリンクが登録のメールアドレスに送られてきます。

3-6-3 CITRUS 結果の確認

- ① CITRUS 完了通知メールのリンクをクリックすると CITRUS 画面が開きます。
- ② Download results をクリックして、結果をダウンロードします。
- ③ Zip フォルダがダウンロードされるので、解凍します。
- ④ フォルダの中には以下の3つが入っています。



- Results フォルダ:結果
- Supporting_files: R やアルゴリズム関連
- Citrus_run_xxx_info.txt: CITRUS 設定情報
- ⑤ Results フォルダを開くと以下の2つのファイルと各モデルの結果フォルダが入っています。モデルフォ ルダは、設定での選択によって作成されます。また、群間に有意差のない場合はフォルダ自体が 作成されません。
- markerPlotsAll.pdf: CITRUS 最初の工程のクラスタリングマップ(1ページにまとめてある)
- markerPlots.pdf: CITRUS 最初の工程のクラスタリングマップ (パラメータごとにページ分かれている)
- sam_results:SAMの結果
- pamr_results:PAMRの結果

- glmnet_results:LASSO via GLMNETの結果	
▶ CITRUS_New_CITRUS_results ● ● ● ● 新しいフォルダー ・ CITRUS_New_CIT 整理 ライブラリに追加 共有 新しいフォ (二) 名前	
☐ citrus_run_51264_info.txt	定等の情報
📜 supporting_files アルゴリズムや R の	関連ファイル
a le results 结果	
 ・ CITRI を理 ・ ライブラリ(・ ・ ・	JS_New_C) results) に追加 ~ 共有 ~ 新しい
🗙 🔶 名前	クラスタリングマップ(まと
The second secon	^{NotsAll.pdf} クラスタリングマップ(別フ
📜 📜 sam_res	sults SAM の結果
amnet ∎	^{ssults} PAMR の結果
	LASSO via GLMNET の結果

3-6-4 CITRUS 結果ファイル

markerPlots.pdfまたはmarkerPlotsAll.pdfを開きます。クラスターツリーの上のパラメータについて、各クラスターの強度が色で示されています。各クラスターがどのような細胞集団であるかを確認することが出来ます。

※CIRUS のクラスターツリーには矢印があり、親→子クラスターとなっています。親クラスターは子ク ラスターを含んでいます。



- ② 各モデルの result ファイルを開きます。以下の5種類のファイルが入っています。
- clusters….pdf-各クラスター内の細胞の表現型を確認するために使用します。各クラスター の表現型はバックグラウンドの表現型とオーバーレイされたヒストグラムで表示されています。クラ スターが赤色、全細胞が青色で示されており、ピークでノーマライズされています。



- featurePlots….pdf-グループ間で有意差のあるクラスターのツリー上の位置を確認するため に使用します。抽出されたクラスターは赤色で示され、親子クラスターの場合は囲われていま す。



- features….pdf-グループ間で有意差のあるものについて、その違いをボックスプロットで表示しています。
 - CITRUS 設定で cluster characterization を「Medians」にしていた場合:プロット 上部のクラスター番号の記載パラメータにおける違いを図示しています。グラフのスケール は、データスケーリング(変換値)に一致します。
 - CITRUS 設定で cluster characterization を「abundance」にしていた場合:プロット上部のクラスターにおける細胞数の違いを示します。値は、CITRUS インプット細胞の総数に対するクラスター内細胞数の割合(0から1の間)で表示されます(例えば、あるサンプルの総細胞数が1,000で、クラスター内の細胞数が100である場合、値は0.1となります)。

注意:abundanceの場合のスケールがLog10と表示されている場合、不具合ですので無視してください。



- features….csv- features….pdf の数値情報のスプレッドシートです。
 注意: Medians 結果の値は CITRUS 解析を開始した時点でのスケール設定で変換されています。
- ModelErrorRate.pdf-予測モデル(PAM, LASSO)の結果のクオリティ調査に使用します。 SAM は相関法であり、予測モデルではないため、このファイルは作成されません。



- ③ また、これら5種類のファイルについて、モデルによって3種類の結果が出力されています。
- 相関モデル (SAM)
 - ・ …fdr_0_01…: FDR を1%としたときの結果
 - ・ …fdr_0_05…: FDR を 5%としたときの結果
 - ・ …fdr_0_10…: FDR を10%としたときの結果
- 予測モデル (PAMR、LASSO)
 - ・ …cv_min…: 交差検証で最もエラー率の低くなる候補数の時の結果
 - ・ …cv_1se…: 交差検証で最も低いエラー率よりも標準偏差が1高い時の結果で、一般的に新規のデータで予測検討する際に採用されます。
 - ・ …cv_fdr_constrained…: (PAMR のみ)設定した FDR 以下のすべての候補

3-6-5 CITRUS 結果ファイルの見方

- (予測モデル PAMR、LASSO の場合) ModelErrorRAte.pdfを開き、cv.1se または cv.minの Model Cross Validation Error Rate(Y 値)を確認します。2 群間の場合、この 値が約 50%だと完全ランダム、3 群の場合は約 66%です。
- ② Features….pdfを開き、違いのあるクラスターの数やその度合いを把握します。
- ③ featurePlots….pdfを開き、違いのあるクラスターがクラスタリングマップのどの位置にあるかを把 握します。
- ④ markersPlots.pdfを開き、違いのあるクラスターが、どのような細胞集団であるかを把握します。Clusters….pdfも参照します。

例)刺激なしと LPS 刺激ありの 2 群で、median & SAM モデルを用いた FDR0.01 の CITRUS 結果: 6 候補が抽出され、その features….pdf を見ると大きく異なる候補が 1 つ、cluster 29993 における pp38 パラメータに着目した。クラスター番号と Feature Plot….pdf から、着目クラスターのクラスタリングマップ上の位置を把握し、これを marker Plots.pdf と照合すると CD33+の集団であることが分かった。



3-6-6 CITRUS 結果から見出した候補クラスターの詳細解析・グラフ化

抽出された候補クラスターを、各ファイル各クラスターとして fcs ファイルにし Cytobank の新規 Experiment に出力することで、詳細解析やグラフ化することができます。 作成された新規 Experiment で、Cytobank の他の機能と組み合わせて様々な作図や検討ができる 他、csv ファイルでエクスポートすると他のソフトウエアでも扱うことができます。

- ① CITRUS 結果から、候補クラスターを抽出しクラスター番号をメモします。
- ② CITRUS 結果画面を表示します。
- ③ Export clusters to new experiment ボタンを押します。
- ④ IDs of clusters to export で、メモした候補クラスター番号にチェックを入れます。

- ⑤ 全細胞の情報も一緒に出力したい場合は、Original files to included でファイルを選択します。
- ⑥ 実行ボタンを押します。



4各種機能

4-1ダウンロード・エクスポート

4-1-1 Experiment アーカイブ (acs) のダウンロード

Experiment 丸ごとをファイル化してダウンロードします。

サーバーの異なる Cytobank で Experiment を移動させる場合や、Cytobank のライセンス契約を一 定期間(90 日間)以上停止する場合、不要な Experiment を端末ローカルで保管する場合に使 用します。

<u>Export Experiment to ACS</u>もご覧ください。

4-1-2統計計算結果・サンプルタブ・ゲート情報・Experiment内 FCS ファイル&添付ファイル・ Illustrations で表示している統計表・Illustrations で表示している細胞情報の FCS/text フ ァイルのエクスポート

Export をご覧ください。

4-2 Experiment の複製

<u>Clone 機能</u>にて Experiment を複製できます。

4-2-1 Experiment の削除

Experimentの削除は2段階になっており、ゴミ箱にいれる段階と完全消去となっています。 ゴミ箱内の Experiment は30日後に自動的に完全消去されます。

4-2-1-1ゴミ箱に入れる削除

Experiment manager で削除する Experiment を選択し、 Delete ボタンを押します。

Experiment Manager		Actions fo	or 1 selected experiment	Apply label	Archive	<u>च</u> Delete	Onselect		
New experiment		Q All e	Q All experiments						
Q All	27		Name				P		
Mine	4	2 1	🔅 SPADE demo (C	Clone)			6		
🤽 Shared with me	10		2012-12 Healthy	Human PBMC w	/ 26 Surface	Markers fro	m .		
🔅 Cytobank curated	15		Fluidigm (Clone)				1		
Public	1		Data for public				3		
â Trash	0		o Data loi pablio				0		

または、対象の Experiment で Actions>Delete を選択します。

	PBMC (fluorescence)	4
<	Experiment Actions Q View Summary	γ
~	🚳 Cloning 🔹 🕨	
	🔁 Import	
2	🔁 Export	3M 3M
	🗓 Reset	one
0	💼 Delete	F.
0	Edit Reagents [admin only]	3
(I)	Course(c)	

4-2-2ゴミ箱に入れた Experiment を戻す・完全消去する

Experiment Manager の Trash を開き、Experiment を選択して Remove from trash (ゴミ 箱から戻す) または Permanently delete (完全消去) をクリックします。

Experiment Manager		Ac	tions fo	r 1 selected experiment	Remove from trash	A Permanently delete	3 Unselect
New experiment		ŵ	Tras	h			
Q AII	27			Name			P
dine Mine	2		1	🔅 SPADE demo (C	Clone)		6
🤽 Shared with me	10						
Cytobank curated	13						
Public	1						
💼 Trash	2						

※アカウントのデータ空き容量は Permanently delete 実行後に増加します(処理にしばらく時間が かかります)。ゴミ箱に入れた状態では空き容量は増えません。

※完全消去(Permanently delete)したデータは復活しませんので十分ご注意ください。

4-3 FCS ファイル詳細情報

メニューバーActions>View summary の一番下 FCS Files に表示されている各ファイルの details をクリックすると、FCS ファイルのメタデータキーワードなど詳細を確認することができます。

FCS Files – 11					
File Name	Sample Name	Sample Tage	Papel	Evente	Sizo
File Name	Sample Name	Sample Tays	Fallel	Events	3126
pbmc_lrs005_il10.fcs 💿 details ┥	IL10	PBMC_LRS005	Panel 1	325374	22.3 MB
pbmc_lrs005_il6.fcs @ details	IL6	PBMC_LRS005	Panel 1	326838	22.4 MB

4-4 FCS ファイル以外のデータを Cytobank ヘインポート (DROP 機能)

DROP 機能では、フローサイトメーターが出力する FCS ファイルのデータ部分と同様、列に対象(シング ルセル)・行に測定数値(発現タンパク質)の形式となっているスプレッドシート(.csv または.tsv)ファイ ルのデータを Cytobank ヘインポートし解析を可能とします。

データ例	種類	プロット/統計	viSNE	SPADE	CITRUS
Cytometry, RNA-Seq, segmented image features	シングルセル 連続変数	Yes	Yes	Yes	Yes
RNA-Seq, Protein expression, clinical	バルク	Yes	Yes	Yes	
features, cell population features	連続変数				
Bulk DNA, clinical features	バルク 順序変数	Yes	Yes		

Premium Cytobank ではインポートできる行数は 100、Enterprise Cytobank では 818 です。

- ① 以下を確認し、.csv または.tsv ファイルを用意します。
- 列がイベント情報(1細胞の情報/1サンプルの情報)・行が測定数値(蛍光強度/発現遺 伝子)
- 行名(パラメータ名)に重複がない
- 非数字情報は削除するか数字情報へ変換します(例:性別は1,2,3とスコア化)。
- 行が 100 以上あるデータの場合、ソートや PCA で行数を減らしてください。
- ② New experiment をクリックします。
- ③ ファイルアップロード画面で、CSV または TSV ファイルをドラッグ&ドロップします。

Add More				
Name	\$	Size	† Type	\$
kidneyCancers_TCGA_RNAseq_f support_tutorial.csv	pkm-	10.3 MB	$csv\toFCS$	\otimes
You can still drag and drop files 🏸	 Convert all csv files to analysis (FCS) files Upload File 	10.33 MB	1 files	

- ④ Upload File ボタンをクリックします。
- ⑤ インポートする範囲を選択します。一般的にはパラメータ部分がグレーになるように指定します。除きた い行は行番号を Indices of columns to skip に入力します。

Select data matrix start										
Select the first (upper left) number in your data matrix. This position will be applied to all files. If you do not see the start of the numerical data matrix, edit your file to remove some of the headers and try again. Ensure you've reviewed our data formatting requirements [2]										
Indices of columns to s	Indices of columns to skip: 1,5,23, etc.									
Learn more about this dataset: http://blog.cytobank.org /2017/08/02/u										
Sample	MT- CO3_fpkm	MT- CO1_fpkm	RN7SL2_fpkm	MT- ND4_fpkm	MT- CO2_fpkm	MT- RNR2_fpkm	MT- ATP6_fpkm	MT- CYB_fp		
TCGA1	6721.620589	7265.666105	19.58601257	8771.732779	5718.798011	7708.309919	5821.272308	3690.2		
TCGA2	15558.84999	15176.78872	44.17243377	11808.18688	9220.986543	8712.944802	8982.356166	7757.3		
TCGA3	10463.71562	10828.1279	108.9822284	7162.489243	5872.184764	8324.959206	4618.079125	4637.4		
TCGA4	25930.08591	17911.69963	164.8555006	26260.24462	16582.45011	10166.54967	18332.3363	15309		
TCGA5	8779.025841	7539.677283	47.38836637	7922.356054	7627.864235	6563.617226	5743.976888	5083.6		
TCGA6	7849.529267	10172.33364	14.88718874	9145.638264	7952.284599	4673.18427	6033.879801	3495.2		

⑥ コンバートしたファイルの概要を確認し、Open experiment ボタンを押します。

4-5アカウントの確認・設定変更

画面右上の My profile ボタン(Username)をクリックすると、設定画面(account settings)ボ タンと現在のデータ使用状況(Experiments の数と使用データ容量)、ログアウトボタンが表示されま す。

Cytobank	Experin	ments Pro	jects	Admin	Help	Cytobar	nkSupport *
Experiment Mana	ger Q	۹.			Cytob Suppo setting	ank rt's Account gs	
• New experiment	t	🖴 Inbo	ĸ		support.accoun 305 experimen 184.6 GB used	it@cytobank.org ts	• Settings
Q All	1599		Name				
🖬 Mine	299				🕞 Log out		ister

4-5-1パスワードの変更

- ① My profile>Account settings (設定画面ボタン)をクリックします。
- ② My account 画面が開きます。

③ Authentication 欄の Update password 部分で New password 等を入力し、 Update Password ボタンを押します。

My account 🥒 Edit	
Authoritization	
Authentication	
Update Password	
Current Password	
New Password	
Confirm New Password	
Passwords should be 8 to 40 characters, v	vith at least one letter and one number.
Update Password	

4-5-2使用データ容量と内訳の確認

My account 画面の Experiment statistics セクションで使用中のデータ容量(使用率)、 Experiment 数、共有している Experiment 数などが確認できます。

My account 🧪 Edit	
Experiment statistics	
Experiment statistics	
Total data owned (i)	128 GB (26% used of 500 GB)
Total experiments owned (1)	311
Uploaded or cloned	197
Experiments shared with user	109
Shared directly	109
Shared only via projects	0

4-5-3ユーザーコネクションのリクエストとリストの確認

My account 画面のスクロール下、User connections セクションで、ユーザーコネクションを確立したい ユーザーへのリクエストとコネクションのあるユーザーリストが確認できます。

My accou	nt 🥒	Edit				
User connections						
Req	uest new u	ser cor	nnection			
use	r@email.com		🕀 Submit			
Exis	ting conne	ctions				
Use	er		Date connected	# shared experiments	# shared projects	
	· •	•••	• ••		2	-

4-5-4 登録 E-mail アドレスの変更

- ① My account 画面上部の Edit ボタンを押します。
- ② Edit my account details 画面が開きます。
- ③ Contact email ボックスに変更する E メールアドレスを入力します。※間違えないようにお気をつけ ください。
- ④ 必要であれば First name と Last name を変更することもできます。
 - ※ Username は変更できない仕様になっています。

60	Cytobank	Experiments	Projects		Help	
<	Edit my acc	count details				
		Gravatar	Personalize or change	your Gravatar.		
		First name	Beckman			
		Last name	Coulter]		
		Contact email	cytobank@beckman.com]		
			Use the address you want to re	 cceive Cytobank email (e.g. jane@companyx.com or john@institu	ition.org)	
		Privacy	🗹 Do not display my email :	address to other users		
		Location	Japan			
		Country	Select a country	~		
		Postal code]		
		Address]		
		City				
	Compa	ny/organization	ВСКК			
	Princ	cipal investigator		Invite a new user		
		Bio				
	Default	new experiment landing page	Set up new experiment	~		
		Default project	None	~		
	Er	mail notifications				
	Receive n	ews and updates from Cytobank				
			Update user settings			

⑤ Update user settings ボタンを押します。

4-5-5その他細かいアカウント設定の変更

Edit my account settings 画面では以下の設定変更も可能です。

- Default new experiment landing page:新規 Experiment を作成した時にデフォルト表示されるページを変更できます。
- Default project: ログインした時にデフォルトで表示される Project を設定できます。
- Email notification: チェックを外すと Cytobank からの通知メール (解析完了など)を解除す ることができます。
- Receive news and updates from Cytobank: チェックを外すとアップデートなどのお知らせを 送らないようにできます。

4-6ログアウト

画面右上の My profile ボタンをクリックし、Logout ボタンを押すとログアウトできます。 パスワード機密保持のため、共通のパソコンでお使いの場合は忘れずにログアウトしてください。また、他の 人も使用する可能性がある端末ではブラウザにパスワードを記憶させないようにご注意ください。

5バージョン対応表

Cytobank version	簡易マニュアル版
7.2	1.000
8.1	2.000
9.3	3.000
10.2	4.000
10.3	4.100

ベックマン・コールター株式会社

社:〒135-0063 東京都江東区有明3-5-7 TOC有明ウエストタワー 本

お客様専用 🚾 0120-566-730 🛛 03-6745-4704 e-mail bckk_ls_web@beckman.com URL https://www.beckman.jp