

ALL IN ONE: FROM RAW DATA TO P-VALUE

Summary visualization and statistical tests now right in your platform

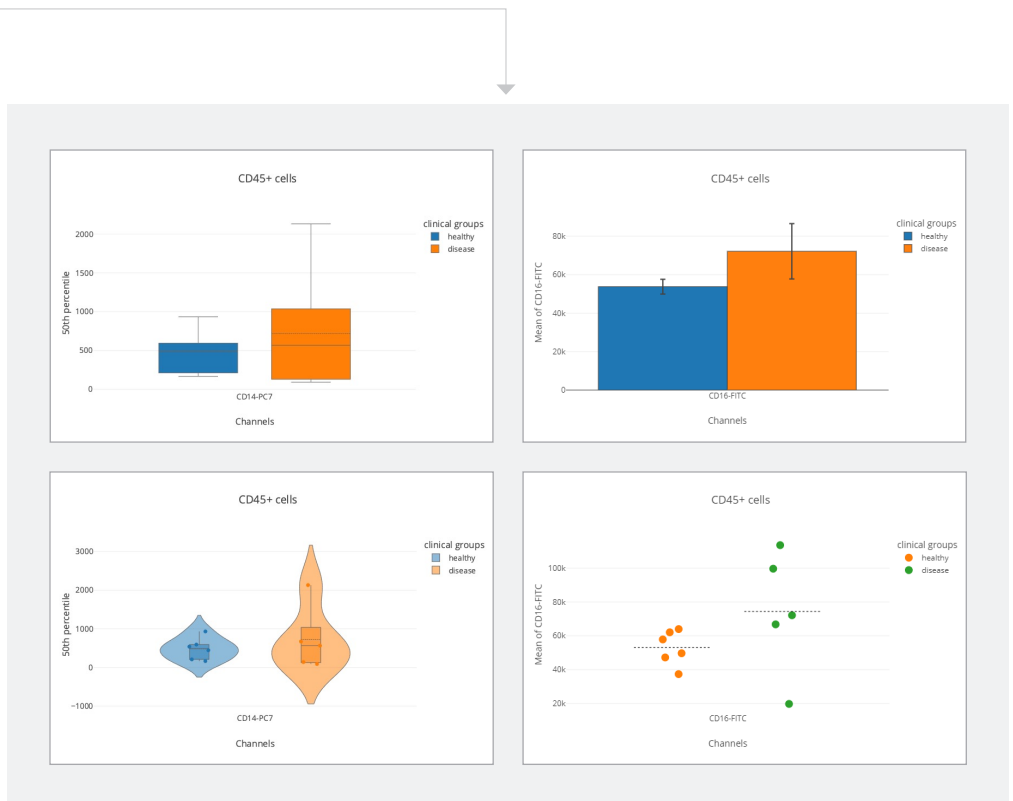
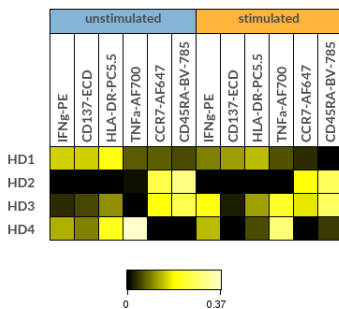
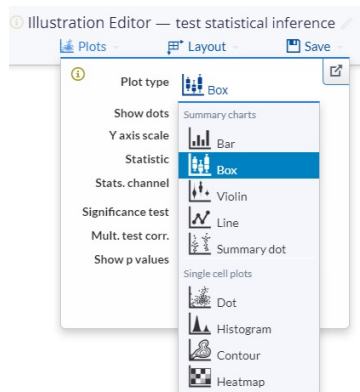


Tired of having to copy-paste percentage or mean values between different software to test for statistical significance? If you need a tool that provides a quick overview of your results, identifying statistically relevant differences between samples, everything within your analysis software, join the Cytobank platform and move from event-level data to summary statistics in few clicks.

Visually Explore Differences Between Groups

Bar charts, Box plots, Violin plots, Line and Summary dot plots and Dimension-Annotated Heatmaps, now available directly in the Illustration Editor, will allow you to easily summarize your meaningful results.

Any changes you make to your gating strategy will automatically propagate through to your summary charts.



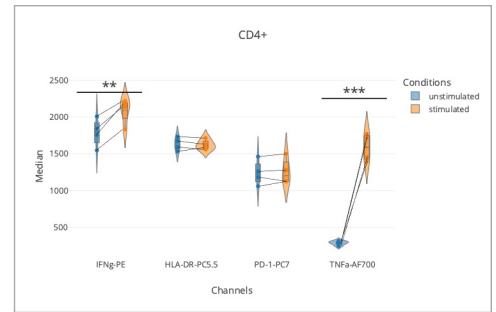
Continued on next page

Assess Statistical Significance Directly In Your Illustration

You can now access Student's T test, ANOVA, Wilcoxon signed-ranks, Mann-Whitney and Kruskal-Wallis statistical tests to evaluate p-values for significant differences among groups. Don't wait until analysis is complete to check if that difference is significant!

Easy-to-interpret graphs and a comprehensive report table will allow you to get the most out of your results.

	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-ranks test	1	2		✓
Kruskal wallis H test	1	> 2		
One-way analysis of variance	1	> 2	✓	
Two-way analysis of variance	2	> 1	✓	

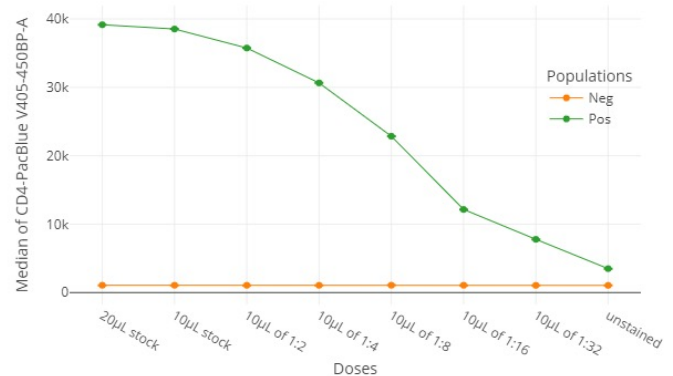


Population	Channel	Condition A	Condition B	Mean A	Mean B	SD A	SD B	n A	n B	p	Significance p-value
CD4+	IFNg-PE	unstimulated	stimulated	1788	2089	165.7	151.1	4	4	0.005766	**
CD4+	HLA-DR-PC5.5	unstimulated	stimulated	1634	1622	77	59.31	4	4	0.6361	ns
CD4+	PD-1-PC7	unstimulated	stimulated	1241	1259	146.4	153.4	4	4	0.544	ns
CD4+	TNFa-AF700	unstimulated	stimulated	287.8	1588	24.15	163	4	4	0.000802	***

Easily Titrate Your Antibodies and Calculate Stain Index

Determining the correct concentration or titer of antibody to use in your experiment is a critical quality control parameter. Summary Charts and our new Stain Index visualization enable you to directly see how stain index and fluorescence intensities change.

Annotate your files with Reagent and Dosage tags, gate on positive and negative populations and build your Illustration. No exporting steps are required.



The Cytobank platform is a cloud-based, single-cell analysis solution for high-dimensional data, visual data management and controlled sharing.



Get Personalized Support

Our Application Specialists at Cytobank_Support@beckman.com
Your local sales representative at Cytobank_Sales@beckman.com



Access online training resources

Support Articles at support.cytobank.org
Videos and more in the Cytobank Learning Center on beckman.com



Access machine learning – assisted analysis with no need to code



Upload any type of data in a tabular format



Securely store & easily share data



Organize data & create auditable workflows



For research use only. Not for use in diagnostic procedures. © 2021 Beckman Coulter and Beckman Coulter Life Sciences. All rights reserved. Beckman Coulter, the stylized logo, and the Beckman Coulter product and service marks mentioned herein are trademarks or registered trademarks of Beckman Coulter, Inc. in the United States and other countries. All other trademarks are the property of their respective owners.

For Beckman Coulter's worldwide office locations and phone numbers, please visit "Contact Us" at beckman.com

21.10.3004.FLOW