

Vi-CELL BLU Analyzer Data Exports and Offline Analysis Instructions

Andrew Wu, Application Scientist, Beckman Coulter Life Sciences [AWU01@BECKMAN.COM]



Introduction

This tech note provides examples of several types of data available for export from the Vi-CELL BLU analyzer and a method to export and import zipped, encrypted sample data for offline analysis. Please refer to the IFU for more detailed information.

Exports can be obtained from six main locations:

- 1. Home screen
 - a. Sample data
 - i. Single sample (.csv)
 - ii. Multiple sample (.csv)
 - b. Loadable template from Create Sample Set screen
 - i. Encrypted sample template (.dat)
- 2. Reports \rightarrow Logs screen
 - a. Audit log (.csv)
 - b. ample activity log (.csv)
 - c. System error log (.csv)
 - d. Scheduled logs
 - i. Audit log (.csv)
 - ii. System error log (.csv)
 - e. Concentration slope log (.csv)
- 3. Quality control screen
 - a. Single sample (.csv)
 - b. Multiple sample (.csv)
- 4. Reports \rightarrow Results screen
 - a. Completed run summary report (.pdf)
 - b. Run results report (.csv and .pdf)
 - c. QC report (.pdf)
 - d. Cell types report (.pdf)
 - e. Instrument status report (.pdf)
 - f. Scheduled data exports
 - i. Encrypted zipped files (.zip)
 - ii. Unencrypted zipped files (.zip)
- 5. Settings → Storage screen
 - a. Configuration template (.cfg)
 - i. The import configuration is enabled for new instruments; once a sample is run, the button will be disabled.
 - b. Encrypted zipped files (.zip)
- 6. PostgreSQL database
 - a. Software dependent formats
 Accelerating Answers | 2

Examples of available data exports:

1. Sample data from home screen:

ScoutUIMasterP	age							- 🗆 X
=	Vi-CELL BLU	\odot		11/21/2024 7:4 Idle	16 AM	93	🧓 🏖 v	
+	Sample Template	Sample ID Sample	Cell Type BCI Default	Dilu v 1	tion Wash Normal (Тад 200 µL) 🔻	• •	
◆ 10/24/20◆ 10/24/20		VI-CELL Sample	set (10-24-2024) Ilitiple Sa	imple (c	sv)			Complete
Status Completed	Pos Samp 02 Sample	le ID Total (x10^ cells/mL 4.6	6) Viable (x10^6) cells/mL 4.6	Viability Total (%) Cells 99.1 12586	Avg Diam (μm) 15.46 asda	ell Type Dilution	Wash Normal (200 µL)	Tag
€ 10/24/20	4 1	Vi-CELL Sample	e Set (10-24-2024)					Complete
© 10/24/21	Single sam		e Set (10-24-2024) e Set (10-24-2024)					Complete
 ♥ 10/21/20 ● 8/4/2020 	24	VI-CELL Sample Vi-CELL Sample	≥ Set (10-22-2024) ≥Set 1					Complete Complete

Single sample summary results (single page CSV icon):

Sample ID	Images for analysis	Cell type	Minimum Diameter (µm)	Maximum Diameter (µm)	Images	Cell sharpness
Example1	100	Yeast	3	20	100	4
Part number	Lot number	Reagent	Expiration	In service date	Effective expiration	1
C06019	836140	Cleaning Agent	5/10/2045 0:00	6/24/2024 0:00	9/22/2024 0:00	
C06019	836140	Disinfectant	5/10/2045 0:00	6/24/2024 0:00	9/22/2024 0:00	
C06019	836140	Buffer Solution	5/10/2045 0:00	6/24/2024 0:00	9/22/2024 0:00	
C06019	836140	Trypan Blue	5/10/2045 0:00	6/24/2024 0:00	9/22/2024 0:00	
Image#	Cell count	Viable cells	Total (x10^6) cells/mL	Viable (x10^6) cells/mL	Viability (%)	Average diameter (µm)
1	235	231	8.52	8.37	98.3	6.97
2	242	239	8.77	8.66	98.8	6.88
3	223	219	8.08	7.94	98.2	7.02

Sample set summary results (multiple page CSV icon):

Sample ID	Cell count	Viable cells	Total (x10°6) cells/mL	Viable (x10^6) cells/mL	Viability (%)	Average diameter (µm)
Example1	22310	21720	8.09	7.87	97.4	6.88
Example2	101990	93128	36.97	33.75	91.3	5.97
Example3	100404	93765	36.39	33.98	93.4	6.02
Example2 Example3	101990 100404	93128 93765	36.97 36.39	33.75 33.98	91.3 93.4	

Sample set file from sample loading screen:

ScoutUIMasterPa	je					- 0
Create Samp	le Set					Vi-CELL
		Sample Set (11-21-2024)				Cave/Load Sample Set Templates
		Sample ID Sample	Cell Type BCI Default	Dilution T	Wash Normal (200 µL)	File Syster, Save/Load
Carous	el		Pos	Sample ID	Cell Type	
05 06 07 08 09		24 23 21 20 19 18 18 17 16 14				
			Ca	ncel Add		
		C				

2. Reports \rightarrow Logs screen exports:

	ada Paga				- 0	×		E Louilliterellige				
	Vi-CELL BLU	11/21/2824 84	E AM	🔒 🎧	Vi-CELL	-		VI-CELL BLU	\odot	11/21/2024	0.01 AM	- 🕄 💮 😂
									2			
	VPB							Eginesare Ecogr				
	bucostaske				6			Audit log	Audit log			
								Sample activity log	Date time	User norme	Event type	Description
		M-GRL ▼ 11/14/2024	11/21/20	24 113				System ense log 🧿	10/20/2021 7:36 /44	v+ctu	er Qugin	Cursule logis OK
	rts 🚺	Provide Antonio Antonio					N (1997)	Concentration slope history log	11/21/28/4 2.36 AV		ext.accpusing tand	A Cap using standard concernation e
		well BOVIAL						Scheduled log exports			except	
		Conments	_						10/0/014 236 200	Michiel	est india	Control Mode
									11/0/2004 9/33 AM	WHEEL	er cogn	Consider login OK
						11					a stoor contrat ionie t	
		Cel court		Cell type					11/0/2021 9:33 AM		est of Bisenscie	Office Mode
									11/5/2004 8:39 AM	VECEL	ext, logis	Console logis OK
		U vale ter tour		C resnertos come					11/5/2024 8:39 AM		evt_accpusingstand	A-Cup using standard concentration a
		Average dameter (sm)		Analysis by					110/004 815 44		except aut officiencois	Cilling Mode
		Average viable clameter (un		Reasalysis by					11/5/2021 0-20 AM	Victur	and leavin	Computer basis Off
Image: Section of the section of t				-		-			11/5/2004-028 AM		ext.ecopoingstand	A Cup using standard concernation :
											and an an an and a sheet	
averte 11/2/2020 (S) To Rouse (S)		ELINGUS E V-CILI BIU Preuts Elogs	•	est0.01		4/3/382	с Sa 820 км Фе			6 8	— 6 х Фили —	
		Vector BUL Participation State Sala Anton Sala Anton Sala Sala Anton Sala Anton Sala Sala Anton Sala Anton Sala Sala Anton Sala Sala Anton Sala Sala Anton Sala Sala An	C Stangular log The Stangular Stang Stangular Stangular St		Canada Ca	and the second s		terd © ∑terda ■		@ @	Bran -	
		Sectors Burger			General Policy General Policy General			Dere Vielen D		a	⊕ xou =	
		Vector Bill			General P			Per d		()	e vai -	
		Autor CLU NUL O Rendo CLU NUL Autor C Autor C	C Supported large The support		American A A A A A A A A A A A A A A A A A A A	An American State		Next Control of the second se		8 •		

Audit log:

Date time	User name	Event type	Description
			Sample: Sample
			Cell Type: BCI Conc Beads
			Position: F-1
			Tag: QC: 2M
9/27/2024 10:29	factory_admin	evt_sampleresultcreated	
			Sample: Sample
			Cell Type: BCI Conc Beads
			Position: E-1
			Tag: QC: 2M
9/27/2024 10:28	factory_admin	evt_sampleresultcreated	
			Sample: Sample
			Cell Type: BCI Conc Beads
			Position: D-1
			Tag: QC: 2M
9/27/2024 10:26	factory_admin	evt_sampleresultcreated	

Sample activity log after selecting single sample:

User name	Date time	Sample ID	Cell type	Analysis type	Status
factory_admin	9/27/2024 10:10	Sample	2M (BCI Conc Beads)	Viable (TB)	Completed
factory_admin	9/27/2024 10:10	Sample	2M (BCI Conc Beads)	Viable (TB)	Completed
factory_admin	9/27/2024 10:10	Sample	2M (BCI Conc Beads)	Viable (TB)	Completed

System error log:

Date time	User name	Error code	Description
	hai sarvica		
10/31/2024 15:32	DCI_Service	2164590081	[Warning] Instrument - Precondition - Plate present: Not met
40/04/000444.00		0004 405005	
10/31/2024 14:29		2231435265	[Warning] Fluidics - General: Nightly clean cycle skipped
10/31/2024 14:29		2198011907	[Warning] Reagents - Reagent pack - Main bay: Empty

Concentration slope log:

Date time	User name	Concentration slope type	# of consumables	Consumable	Assayvalue (x10^6)/mL	Lot number	Expiration	Slope	Intercept
12/31/2017 19:00	factory_admin	cal_Concentration	3	Factory Default	2	n/a	1/1/2100	368	(
12/31/2017 19:00	factory_admin	cal_Concentration	3	Factory Default	4	n/a	1/1/2100	368	(
12/31/2017 19:00	factory_admin	cal_Concentration	3	Factory Default	10	n/a	1/1/2100	368	(
12/31/2017 19:00	factory_admin	cal_Size	3	Factory Default	2	n/a	1/1/2100	0.53	-5.53
12/31/2017 19:00	factory_admin	cal_Size	3	Factory Default	4	n/a	1/1/2100	0.53	-5.53
12/31/2017 19:00	factory_admin	cal_Size	3	Factory Default	10	n/a	1/1/2100	0.53	-5.53

Scheduled audit and system error log:

Date time	User name	Event type	Error code	Location	Description
9/27/2024 10:43	Scheduled data exports	evt_Instrumentdataexported	-	Audit	Exported: 1. Result of Sample: <sample> User : <factory_admin> Timestamp : <sep 14:14:41="" 2024="" 27="" utc=""> Celltype : <bci beads="" conc="">. 2. Result of Sample:</bci></sep></factory_admin></sample>
9/27/2024 10:29	factory_admin	evt_sampleresultcreated	-	Audit	Sample: Sample Cell Type: BCI Conc Beads Position: F-1 Tag: QC: 2M
9/27/2024 10:28	factory_admin	evt_sampleresultcreated	-	Audit	Sample: Sample Cell Type: BCI Conc Beads Position: E-1 Tag: QC: 2M
9/27/2024 10:26	factory_admin	evt_sampleresultcreated	-	Audit	Sample: Sample Cell Type: BCI Conc Beads Position: D-1 Tag: QC: 2M
9/27/2024 10:25	factory_admin	evt_sampleresultcreated	-	Audit	Sample: Sample Cell Type: BCI Conc Beads Position: C-1 Tag: QC: 2M
9/27/2024 10:20	factory_admin	evt_sampleresultcreated	-	Audit	Sample: Sample Cell Type: BCI Conc Beads Position: B-1 Tag: QC: 2M
9/27/2024 10:15	factory_admin	evt_sampleresultcreated	-	Audit	Sample: Sample Cell Type: BCI Conc Beads Position: A-1 Tag: QC: 2M
9/27/2024 10:08	factory_admin	evt_qcontrolcreate	-	Audit	Quality Control Name: 2M Cell Type: "BCI Conc Beads" Lot information: 123 Expiration Date: 2024-Sep-27 Assay value: 2.010 Plus/Minus: 10.0 %
9/27/2024 10:05	factory_admin	evt_login	-	Audit	Console login OK
9/27/2024 10:04		evt_offlinemode	-	Audit	Offline Mode
9/27/2024 10:03		evt_offlinemode	-	Audit	Offline Mode
9/26/2024 14:41	factory_admin	evt_login	-	Audit	Console login OK
9/26/2024 14:41		evt_offlinemode	-	Audit	Offline Mode
9/26/2024 14:34		evt_offlinemode	-	Audit	Offline Mode

3. Quality controls screen:

🗈 Scou	tUIMasterPag	e						- 0	×		
≡		Vi-CELL BLU	\odot	11/21/	2024 10:02 AM Idle		🔫 🧓 🏖 Vi-CELL -				
Qı	uality c	ontrols		Historical 🔊	Quality contro	ols 📄 Image 🕻	Graphs 🔤		>		
	Name	Cell type	Assay parameter	Lot number	Assay Multiple	e Samples (csv)	expiration	Comments			
		BCI Conc Beads	Total (x10^6) cells/mL	124	Single	e Sample (csv)	10/25/2024		٦		
		BCI Viab Beads	Total (x10^6) cells/mL	746	50.00	+/- 10 %	10/24/2024				
		BCI Conc Beads	Total (x10^6) cells/mL	120	2.07	+/- 10 %	2/25/2021				
		BCI Conc Beads	Total (x10^6) cells/mL	121	4.03	+/- 10 %	2/14/2021				
		BCI Conc Beads	Total (x10^6) cells/mL	118	10.03	+/- 10 %	9/12/2020				
		BCI Conc Beads	Total (x10^6) cells/mL	117	0.49	+/- 10 %	6/20/2020				

Sample set summary results (multiple page CSV icon):

Sample ID	Cell count	Viable cel	Total (x10	Viable (x1	Viability (Average d	Average v	i Average c	Average vi	i Average ce	Average b	Bubble co	Cluster co	Images fo	Cell type
Sample	575	0	0.21	0	0	10.15	0	0.98	0	5	131	0	0	100	2M (BCI Conc Beads)
Sample	575	0	0.21	0	0	10.16	0	0.98	0	5	131	0	0	100	2M (BCI Conc Beads)
Sample	576	0	0.21	0	0	10.15	0	0.98	0	5	131	0	0	100	2M (BCI Conc Beads)

Single sample summary results (single page CSV icon):

Sample ID	Images for analysis	Cell type	Minimum Diameter (µm)	Maximum Diameter (µm)	Images	Cell sharpness	Minimum	Decluster	Aspiration	Viable sp	Viable sp	Mixing cy	cAnalysis date/time
Sample	10	2M (BCI Conc Beads)	2.5	12	100	17	0.75	High	3	50	1	3	9/27/2024 10:14
Part number	Lot number	Reagent	Expiration	In service date	Effective expiration								
C06019	65432	1 Trypan Blue	10/27/2024 0:00	9/27/2024 0:00	10/27/2024 0:00								
Image#	Cell count	Viable cells	Total (x10^6) cells/mL	Viable (x10^6) cells/mL	Viability (%)	Average diameter (µm)	Average vi	Average c	i Average vi	Average c	Average b	Bubble co	Cluster count
1		6 (0.22	0	0	10.05	0	0.98	0	6	130	0	1 0
2		4 (0.15	0	0	10.52	0	0.99	0	4	130	0	1 0
3		9 (0.33	0	0	9.82	0	0.98	0	9	131	0	1 0
4		4 0	0.15	0	0	9.69	0	0.98	0	4	132	0	0
5		6 (0.22	0	0	10.18	0	0.98	0	6	132	0	0

4. Reports \rightarrow Results screen:

≡	Vi-CELL BL	U	2		11/	14/2024 10:29 AM Idle		300	🔠 factory_admin 📔 🗕
ۍ ۱		ogs							
8		<u> </u>							
X (Quality controls		Instrument status						PDF
o e s			Print title Vi-CELL BLU	v1.4.4.1					
as N. P	Reports		Comments						
0 :									
ہا ھ									
 •									

Completed run summary report (select up to 3 parameters):

		v	i-CELL BLU	J v1.4.4.	1		
BECKMAN COULTER Life Science	l 5					Serial number	: LocalInstrumentDe 9/27/2024 10:35:39 A
			Completed r	un summar	У		
Sample ID	Analysis date	Total (x10^6) cells/mL	Viable (x10^6) cells/mL	Viability (%)	Cell type	Analysis by	Signature
Sample	9/27/2024 10:29:49 AM	0.42	0.00	0.0	2M (BCI Conc Beads)	factory_admin	
Sample	9/27/2024 10:28:26 AM	0.42	0.00	0.0	2M (BCI Conc Beads)	factory_admin	
Sample	9/27/2024 10:26:50 AM	0.42	0.00	0.0	2M (BCI Conc Beads)	factory_admin	
Sample	9/27/2024 10:25:15 AM	0.21	0.00	0.0	2M (BCI Conc Beads)	factory_admin	
Sample	9/27/2024 10:19:55 AM	0.21	0.00	0.0	2M (BCI Conc Beads)	factory_admin	
Sample	9/27/2024 10:14:41 AM	0.21	0.00	0.0	2M (BCI Conc Beads)	factory_admin	
yeastfull	8/7/2024 4:18:48 PM	36.39	33.98	93.4	Yeast	bci_service	
yeasthungry	8/7/2024 4:16:50 PM	36.97	33.75	91.3	Yeast	bci_service	
yeastold	8/7/2024 4:14:53 PM	8.09	7.87	97.4	Yeast	bci_service	

Run results report:

From open sample (folder icon), with multiple page CSV icon selected:

Sample ID	Cell count	Viable cells	Total (x10^6) cells/mL	Viable (x10^6) cells/mL	Viability (%)	Average diameter (µm)	Average viable diameter (µm)
Sample	575	0	0.42	0	0	10.17	0
Sample	575	0	0.42	0	0	10.17	0
Sample	574	0	0.42	0	0	10.16	0
Sample	576	0	0.21	0	0	10.15	0
Sample	575	0	0.21	0	0	10.16	0
Sample	575	0	0.21	0	0	10.15	0

Accelerating Answers | 7

PDF icon after selecting sample (1st annotated image checked):



QC report after selecting one QC name:



Cell types report:

				Vi-C	ELL BL	U v1.4	.4.1					
Serial number : LocalInstrumentDe Life Sciences 9/27/2024 10:39:46 A Cell types User name factory_admin												
Cell type	Minimum diameter (µm)	Maximum diameter (µm)	Images	Cell sharpness	Minimum circularity	Decluster degree	Viable spot brightness (%)	Viable spot area (%)	Aspiration cycles	Mixing cycles	Concentration Adjustment Factor (%)	
BCI Default	1.00	50.00	100	7.0	0.10	Medium	55.0	5.0	3	3	0.0	
Mammalian	6.00	30.00	100	7.0	0.10	Medium	55.0	5.0	3	3	0.0	
Insect	8.00	50.00	100	7.0	0.10	Medium	55.0	5.0	3	3	0.0	
Yeast	3.00	20.00	100	4.0	0.10	High	45.0	2.0	3	3	0.0	
BCI Viab Beads	5.00	25.00	100	22.0	0.50	Medium	50.0	1.0	3	3	0.0	
BCI Conc Beads	2.50	12.00	100	17.0	0.75	High	50.0	1.0	3	3	0.0	
BCI L10 Beads	5.00	15.00	100	22.0	0.50	Medium	50.0	1.0	3	3	0.0	
Type0	1.00	50.00	100	7.0	0.10	Medium	55.0	5.0	3	3	0.0	
Type1	1.00	50.00	100	7.0	0.10	Medium	55.0	5.0	3	9	0.0	
Type2	1.00	50.00	100	7.0	0.10	Medium	55.0	5.0	9	9	0.0	
Туре3	1.00	50.00	100	7.0	0.10	Medium	55.0	5.0	9	3	0.0	
Туре3 - Сору	25.00	50.00	100	7.0	0.10	Medium	55.0	5.0	9	3	0.0	
Mammalian - 6x mix	6.00	30.00	100	7.0	0.10	Medium	55.0	5.0	3	6	0.0	

Instrument status report:

	Vi-CELL BLU v1.4.4.1						Vi-CELL BLU v1.4.4.1							
BECKMAN COULTER Life Sciences		Serial numbe	r : LocalInstrumentDefa 9/27/2024 2:42:43 PM	6	BECKM COULTER Life Scien	AN ces			Se	rial number : L 9/:	ocalinstrumentDefa 27/2024 2:42:43 PM			
				L	ast concent	tration slo	pe date	12/31/2017						
	Instrument st	atus report					ssav value	Expiration date	Lot num	ber				
							(x10^6)/mL	Expiration date	201110					
About instrument				2	M		2.00	1/1/2100		-				
Vi-CELL BLU software version	on 1.4.4.1	Serial number	LocalInstrumentDef	4	м		4.00	1/1/2100		-				
Backend software version	1.4.4.25503		a	1	OM		10.00	1/1/2100		-				
Firmware version	N/A	No. of samples lifetir	me 6											
UI software version	1.4.4.1	Instrument health	OK											
Image analysis software ver	sion 4.3.8			Stor	age									
Syringe pump firmware ver	sion N/A			Ш.,										
Camera firmware version	N/A			2	06.7 GB fre	e of 920.0	GB							
_				•)ther	708.5 G	B							
Reagents					ata	2.2 GB								
Container	Trypan Blue			E	xport	2.4 GB								
Part number	4													
Lot number	C10819			LOW	/ level co	ntrois								
Effective expiration	10/27/2024				Optics			Sensor status st	ate	Motor status	state			
Uses total	300				Brightfie	Id LED	20.00	Carousel	•	Radius home	•			
Uses remaining	300						20000	Tube		These house				
Remaining warte tube trav	120				nosition	locus	30000	Tube		Theta nome				
capacity	120				Elow cell	death	0.06	Reagent door		Probe home	•			
					(mm)	uepui	0.00	Reagent pack	•	Focus home	•			
Users										Reagent upper	•			
1. factory_admin										Reagant lower	17			
										Reagent lower	н			
Cell types					Status :	state		Voltage		Temperature				
1. BCI Default 2	. Mammalian	3. Insect	4. Yeast		Caroural	/06 wall		+ 3 31/	2 21 V	Control board	25.2%			
5. BCI Viab Beads 6	. BCI Conc Beads	7. BCI L10 Beads	8. Type0		carouse	/ so wen		+5.54	5.51 4	control board	25.2 C			
9. Type1 1	0. Type2	11. Type3	12. Type3 - Copy		Valve		F	+5V Sensor	5.02 V	CPU	25.3°C			
13. Mammalian - 6x mix 1	4. Mammalian - 9x mix	15. Mammalian - 1x mix	16. Yeast_Custom		Syringe		0	+5V Circuit	4.98 V	Optical case	25.1°C			
-	Сору				Probe		0	+12V	11.67 V					
17. Yeast - VSA test					Motor th	eta	1789564	+24V	24.01 V					
Analysis type							560							
ruary so type					Motor ra	dius	0							
1. Viable (TB)					Motor fo	cus	-1							
Concentration slope					Motor re	agent	1150000							
User name	factory_admin			ll v ,	ctive	Inactiv	. X (Jake							
Slope	368.00			" "	CUVE	- mactiv	e ee Unki							
Intercept	0			Com	nents									

Scheduled encrypted and unencrypted exports (Settings>Reports):

Note: Unencrypted files do not have image annotations and cannot be uploaded to the offline software.

	Vi-CELL BL	U	\odot	9/30/2	024 12:38 PM Idle	120	, 🛛 🚨 facto	ry_admin –
ស្		ogs						
3	Cell types							
\mathbf{M}	Quality controls		Schedu	uled data exports				
¢ °	Settings			🖌 🧪 Edit 🔲 🔟	elete			
== 	Reports		Enabled	Schedule Name	Comments		Export Filename (.zip)	Repeat
\odot	Sign out			Scheduled Encrypted Export	Encrypted Export Example		Encrypted Export	Once
Ð	Lock			Scheduled Unencrypted Export	Unencrypted Export Example		Unencrypted Export	Once
₽								

Input screen for scheduled data export options:

Vi-CELL BLU	0
Schedule Name	Enabled
Comments	optional description of schedule
Export to	path to destination folder
Export filename	
Encrypt export?	Export CSVs and non-encrypted images
Notify	email address to send confirmation of export
Repeat	Once
Export on	9/30/2024 15 at 0 : 00 24 Hour 🔻
Export samples	From 9/30/2024 15 To 9/30/2024 15
	Additional data filters 📀
	Save Cancel

5. Settings \rightarrow Storage screen:

🌣 Settings 🛛 🔽 Securi	ty 🥃 Storage
Storage 187.3 GB free of 920.0 GB Other Data Export	Clean up/Expc User name F All Sample ID
System backup Export configuration	

Configuration export from storage screen file:

20240927_104744.cfg

6. PostgreSQL read-only database Detailed Results table (PGAdmin4):

ResultDate timestamp without time zone	ProcessingStatus smallint	TotCumulativeImages smallint	TotalCellsGP /	TotalCellsPOI 🖋	POIPopulationPercent double precision	CellConcGP double precision	CellConcPOI double precision	AvgDiamGP double precision
2024-09-27 14:15:37	0	1	4	0	0	147200	0	9.6893978119
2024-09-27 14:25:35	0	1	7	0	0	257600	0	10.391514097
2024-09-27 14:29:58	0	1	4	0	0	294400	0	10.524515152
2024-09-19 20:24:04	0	1	956	914	95.606694561	34649853.564	33127579.663	5.897874291
2024-09-27 14:25:36	0	1	5	0	0	184000	0	10.249420929
2024-08-07 20:16:50	0	1	1003	902	89.930209372	36353350.549	32692644.263	6.0237758771
2024-09-27 14:20:45	0	1	4	0	0	147200	0	9.6893978119
2024-09-27 14:26:58	0	1	6	0	0	441600	0	10.054876804
2024-09-27 14:28:35	0	1	4	0	0	294400	0	10.524515152
2024-09-19 20:24:01	0	1	1023	956	93.450635386	37078242.883	34649853.564	6.0469955773

Instructions for exporting zipped encrypted files from the storage screen and importing into the offline software for remote troubleshooting, archiving and analysis.

Sample data may be exported for troubleshooting and optimizing cell types, archiving sample runs and reviewing data on an off-instrument computer. This may be done using an encrypted zipped sample data export file. The exported file will include all cell types, historic data including annotated images, sample names, run parameters and user credentials, and is identical to what can be viewed on the instrument. Please refer to IFU appendix E for more information.

To export the zipped encrypted file:

1. Navigate to settings



2. Navigate to storage, filter by date (1), select samples (2), and click the export icon (3)

≡ vi	-CELL BLU	\bigcirc	9/27/2024 12:4 Idle	2 PM	🧓	actory_admin -
🔅 Setting	s 🥂 🤽 Security	Storage	about 🥪	? Help		
206.8 GB	free of 920.0 GB	Clean up/Expe User name All	ort From: 9/20/2024 15	To: 9/27/2024 15	S	00
Data		Sample ID	Date time	User name	Images	Cell type
Concert.		Sample	9/27/2024 10:29:49	AM factory_admin	100	2M (BCI Conc Beads)
Export	9	Sample	9/27/2024 10:28:26	AM factory_admin	100	2M (BCI Conc Beads)
		Sample	9/27/2024 10:26:50	AM factory_admin	100	2M (BCI Conc Beads)
		Sample	9/27/2024 10:25:15	AM factory_admin	100	2M (BCI Conc Beads)
System		Sample	9/27/2024 10:19:55	AM factory_admin	100	2M (BCI Conc Beads)
Export cont	iguration	Sample	9/27/2024 10:14:41	AM factory_admin	100	2M (BCI Conc Beads)
Import con	figuration					

3. Enter the number of images to export and the export location



The file will populate at the specified location. Encrypted sample sets are in .zip format.

Example of exported file:

Encrypted_20240927_104301.zip

Note: Removing the .zip and changing the format when renaming may not allow future import.

Instructions for importing the encrypted zipped file into the offline software:

- 1. Locate the import_data.bat file from C:\Instrument\Software.
- 2. Double-click to run.
- 3. Paste the zipped file location into the command prompt window.

The file will load into the software and once completed, the data will be available the next time the software is open. The Vi-CELL BLU offline software must be installed on the host computer.

Summary:

The Vi-CELL BLU analyzer allows versatile data handling options with various methods and formats for data exports and reports. Scheduled exports can simplify data archiving. The encrypted zipped sample data export option enables remote troubleshooting and data analysis on off-instrument computers. For additional assistance with data handling or the offline software, please reach out to Beckman Coulter Life Sciences.



Product is not verified or validated for use in diagnostic procedures.

©2025 Beckman Coulter, Inc. All rights reserved. Beckman Coulter, the stylized logo, and 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