



Specification Comparison of Vi-CELL XR Analyzer and Vi-CELL BLU Analyzer

The Vi-CELL BLU analyzer leverages the key performance features of the Vi-CELL XR analyzer but incorporates many design improvements that our customers have requested over the years. The specifications listed below are for informational purposes only and subject to change. Please check beckman.com for most current information.



Comparison of system specifications

Criteria	Vi-CELL XR Analyzer	Vi-CELL BLU Analyzer
Single sample analysis time	180 sec	130 sec Normal Mode 90 sec Fast Mode
Carousel loading	12-position carousel	24-position carousel
Microplate loading	-	Standard multi-well microplates
Automation support	Automation Ready	Automation Ready
Cell Viability Method	Trypan blue dye exclusion	Trypan blue dye exclusion
Cell Size Range	3-70 μm	2-60 μm
Focus System	Autofocus	Autofocus
Image Transfer	Firewire	USB 3.0
Camera Sensor Size	1280x960	2048x2048
Captured Image Size	882x661 pixels	2048x2048
Analyzed Image Dimensions	832x611 μm	1000x1000 μm
Effective Resolution	1 μm per pixel	0.5 μm per pixel
Imaged volume per frame	43 nL	65 nL

Criteria (Continue)	Vi-CELL XR Analyzer	Vi-CELL BLU Analyzer
Minimum Sample Volume	500 µL	200 µL Normal Mode 170 µL Fast Mode
Cell Density Range	5x10 ⁴ – 1x10 ⁷ cells/mL	5x10 ⁴ – 1.5x10 ⁷ cells/mL
Counting Accuracy	± 10%	± 10%
Counting repeatability	± 10%	± 5%
Reagent Pack	200 activities per pack	310 activities per pack
Facilitates 21 CFR part 11 compliance	Yes	Yes
Data formats	TXT, MS Excel	CSV, PDF, Encrypted Binary
Touchscreen interface	Depends on external PC	Integrated
Computer System	External PC, Win 7 OS	Integrated PC, Win 10 OS, 64 bit
Computer System Specifications	Varies depending on configuration ordered	i7-6700 Intel CPU 4 cores at 3.4 GHz, 16 GB DDR4-2400 RAM, 500 GB Samsung EVO 850 SSD
Connectivity	Via PC system	4 x USB 3.0 ports, Gigabit Ethernet
Height	44.5 cm (17.5 in) excluding PC	45 cm (16.5 in)
Width	38 cm (15 in) excluding PC	54 cm (21 in)
Depth	41 cm (16 in) excluding PC	42 cm (18 in)
Weight	11.3 Kg (25 lb) excluding PC	28 Kg (63 lb)
Power	Power: 50 watts (65 watts max) Voltages: 100V, 120V, 220V, or 240 V50/60 Hz	Power: 50 watts (65 watts max) Voltages: 100V, 120V, 220V, or 240 V50/60 Hz

Recommended Control Beads

C09145 Vi-CELL BLU 50% Viability Control

C09147 0.5M Concentration Control, 0.5x10⁶/mL

C09148 2M Concentration Control, 2x10⁶/mL

C09149 4M Concentration Control, 4x10⁶/mL

C09150 10M Concentration Control, 10x10⁶/mL

6602796 L10 Standard, nominal 10 µm, Latex Particle (NIST Traceable) 1 x 15 mL



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

© 2024 Beckman Coulter, Inc. 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.

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

2023-GBL-EN-102551-v2

