



Cydem VT Automated Clone Screening System

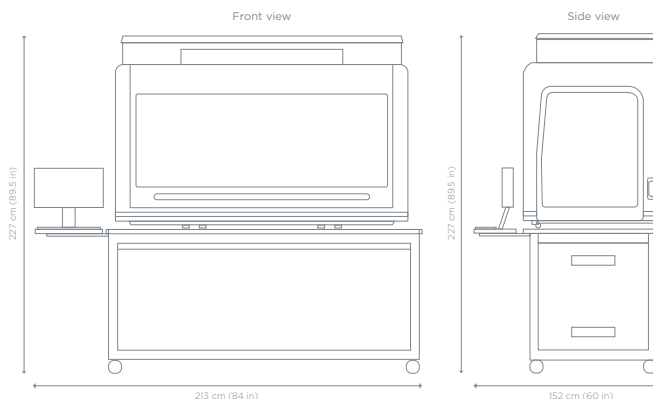
Technical Instrument Parameters

Specifications	Description
Combined size (As installed - Instrumentation, HEPA filter, Workstation cart with monitor)*	- Width: 213 cm (84 in) - Depth: 152 cm (60 in) - Height: 227 cm (89.5 in) with door closed - Height: 232 cm (91.5 in) with door open
Instrument chassis (Cydem VT liquid handler)	- Width: 170 cm (67 in) - Depth: 81 cm (32 in) - Height: 112 cm (44 in) with door closed
Workstation cart size**	- Width: 183 cm (72 in) - Depth: 81 cm (32 in) - Height: 84 cm (33 in)
Weight (As installed - Instrumentation, HEPA filter, Workstation cart with monitor)	630 kg (1389 lbs) with reagents and consumables 581 kg (1281 lbs) without reagents and consumables

* Monitor and mounting arm positioned at a 45° angle.

** Workstation cart is always included in the Cydem VT system.

System Dimensions



The Cydem VT Automated Clone Screening System combines multiple trusted Beckman Coulter technologies into one advanced platform that empowers lab managers with a far more effective, high-throughput method for pool and top clone screening in cell line development—offering increased confidence that promising clones don't get overlooked in the process of consistently selecting the best.

- High-throughput clone and minipool screening in cell line development
- Automated, walk-away operation
- pH-controlled fed-batch cultivation of 96 parallel cultures with individual gassing
- On-deck cell count, viability and antibody titer analysis
- Fully integrated system
- All results in one software

Environmental Requirements

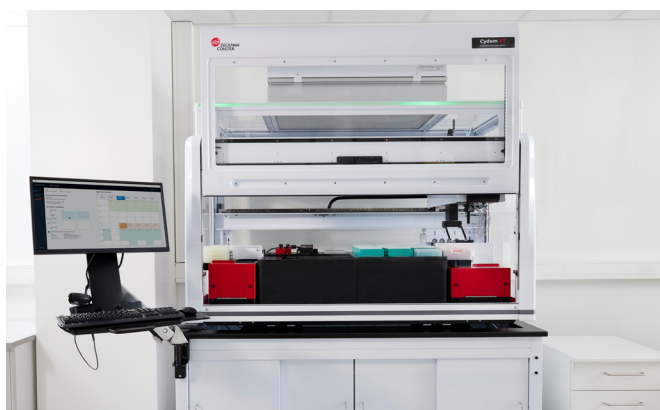
Specifications	Description
Electrical requirements	- Europe (230 V-option): Each Rated 16A. Requires two 230 VAC, 16A CEE 7/7 outlets (same or separate 16A service) - North America (115 V-option): Each Rated 16A. Requires two 120 VAC, 16A NEMA 5-20 outlets (same or separate 20A service)
Fluid requirements:	- DI water, degassed - Isopropyl alcohol - Reagent packs (cell health and titer) 0.5 M NaOH
Gassing requirements	- CO₂, O₂, N₂ 1 - 1.5 bar inlet pressure - Connection fitting to 4 mm OD tubing for all 3 gases
Access space	92 cm (36 in.) access all around the instrument during service required 122 cm (48 in.) in front of the door required at all times
Floor	Typical lab floor with no major slope

Part numbers	Description
D03895	Cydem VT Automated Cell Culture System, 115 V
D03896	Cydem VT Automated Cell Culture System, 230 V
C84309	Cydem VT System Cell Culture Standard Experiment Kit
C86430	Cydem VT System IgG Titer Standard Experiment Kit

Operating Conditions

Specifications	Description
Operating temperature	19 - 25 °C
Humidity	20 - 85 % rH (non-condensing) @ 30 °C
Altitude	Up to 2000 m (3562 ft)

Part numbers	Description
C86225	Cydem VT System IgG Titer Extended Experiment Kit
C84310	Cydem VT System Cell Health Test Kit
C87046	Human IgG Standard Kit, 10 mg
C88594	Cydem VT System Consumables Kit



From one clone can arise a therapy that helps millions of people. From one lab manager can flow a string of selection successes that elevates the team and speeds discovery of the next big cure. It's not imagination. It's more a reality than ever.

The Cydem VT system combines multiple trusted technologies into a standalone system. It's a single platform that allows user-independent cell growth, and antibody titer measurement to accelerate the selection of the best clones.

Discover the difference you can make with an instrument like no other at beckman.com/cydemvt.



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

©2025 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-104324-v5

