

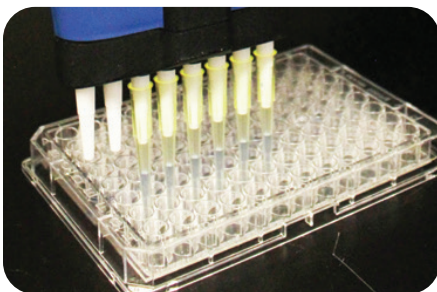


revvity

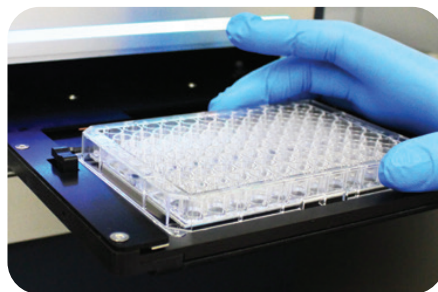
Increase your productivity with high-throughput cell counting.

Cellaca[®] MX high-throughput cell counter

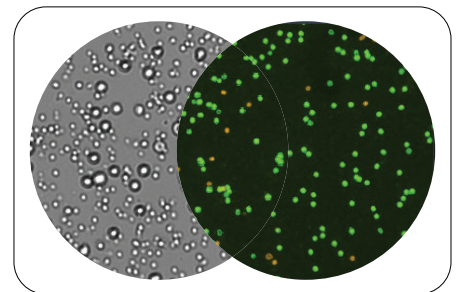
- High-throughput cell counting and analysis with brightfield and two fluorescent channels for use with multiple cell types and lines (including primary and messy samples)
- Count up to 24 samples with fluorescence in 3 minutes or less using a plate-based format
- Low loading volume conserves precious primary samples for additional downstream analysis in bioprocessing or cell line development workflows
- Automation integration ability with optional API
- Optional 21 CFR part 11 module for compliance with regulatory requirements



1. Pipette sample



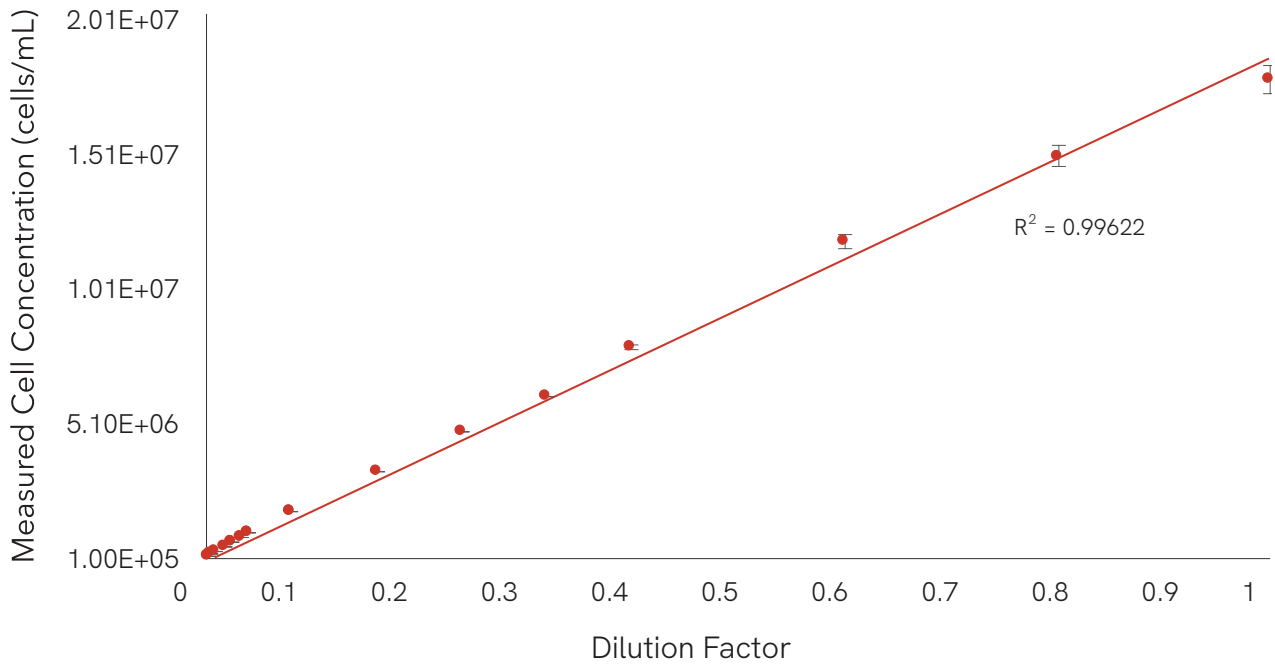
2. Insert plate and count



3. Get images and data

For research use only. Not approved for diagnostic or therapeutic use.

Total Jurkat Cell Concentration
Linear Fit Between 1.8×10^5 to 1.8×10^7 cells/mL

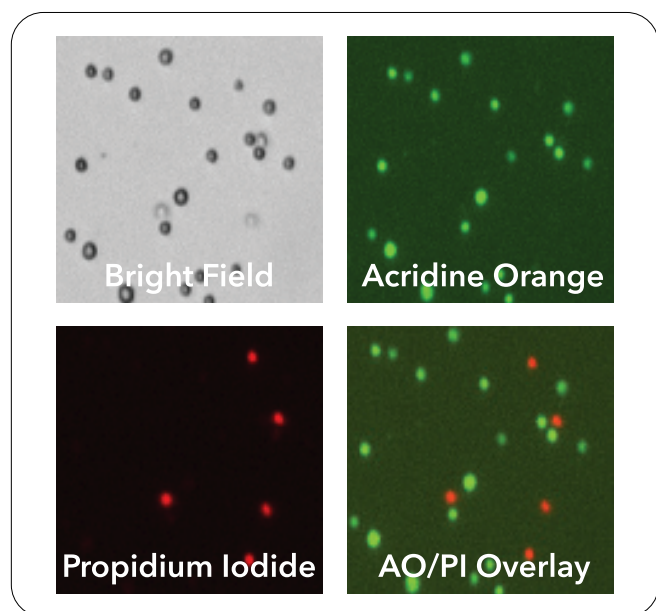


Mean cell concentration cells/mL (n=9)	%CV
1.8E+05	0.7%
2.3E+05	0.4%
3.2E+05	0.4%
5.0E+05	0.3%
6.9E+05	0.3%
8.5E+05	0.5%
1.0E+06	0.5%
1.9E+06	0.1%
3.4E+06	0.3%
4.9E+06	0.2%
6.3E+06	0.2%
8.2E+06	0.1%
1.2E+07	0.7%
1.6E+07	1.9%
1.8E+07	2.4%

System specifications	
Dimensions	13 in x 13 in x 16 in (33 cm x 33 cm x 41 cm)
Weight	42 lbs (19 kg)
Excitation LED	365, 470, 527 and 620 nm
Emission Filters (bandpass, center wavelength)	452, 534, 605, 655 and 692 nm
Volume per Sample	50 -200 μ L
Counting Speed per Sample	2 sec - 17 sec (assay dependent)

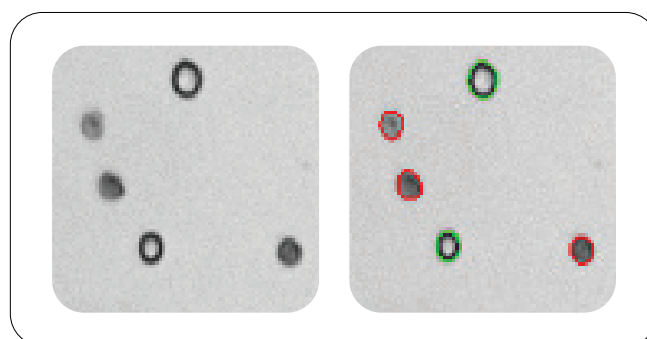
Compatible dyes and fluorophores				
Hoechst	GFP	APC	CFDA	CFP
DAPI	PI	JC-1	Calcein AM	YFP
AO	PE	FITC	RFP	7AAD

Counting of human PBMCs stained with AO/PI

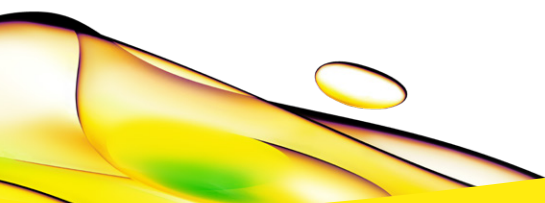


A human PBMC sample stained with acridine orange/propidium iodide (AO/PI) and imaged on the Cellaca MX system.

Counting of CHO cells stained with Trypan blue



CHO cells stained, imaged and automatically analyzed by the Cellaca MX system. Live cells are outlined in green while dead, Trypan blue (TB) cells are outlined in red



revvity