



revvity

# Increase your productivity with high-throughput cell counting.

## Cellaca<sup>®</sup> MX High-throughput automated cell counter

### Speed

Count up to 24 samples in only 48 seconds with Trypan blue exclusion, and count complex samples with fluorescence in as little as 3 minutes for 24 samples. The Cellaca<sup>®</sup> MX is one of the fastest cell counters on the market.

### Small sample volume

The small sample size - only 25  $\mu$ L - required by the Cellaca MX counter makes it possible to readily analyze precious bioprocessing or cell line development samples and enables the same sample source to be used for additional downstream analysis.

### Consistency

High sample to sample and instrument consistency.



### Automation

Standardized API Automation compatibility with automation-ready counting plates.

### Flexibility

Cellaca MX has structured and adaptable software with customizable applications, dashboards, and data templates along with LIMS compatibility for seamless data flow with over 200 cell libraries and 19 different assays.

Number of samples	Amount of time
<b>Cellaca MX trypan blue</b>	
24	48 seconds
480	30 minutes
<b>Other cell counter</b>	
12	36 minutes
480	24 hours

Feature	Cellaca MX BF	Cellaca MX FL2	Cellaca MX FL5	Competitor platform	Competitor legacy platform
Channels	Brightfield	Brightfield, Green, Red	Brightfield, Blue, Green, Red, Far Red	Brightfield *video imaging through flow	Brightfield *video imaging through flow
Number of channels	1	4	14	1	1
Commonly used dyes	Trypan blue	Trypan blue, AO/ PI, Calcein AM/PI	Trypan blue, AO/ PI, Hoechst, DAPI, CFDA, Calcein AM, 7AAD	Trypan blue	Trypan blue
Counting speed	2 sec per sample (24 samples - 48 seconds)	2 - 17 sec per sample (24 samples - 48 sec with bright field or 3 min with fluorescence)	2 - 17 sec per sample (24 samples - 48 sec with bright field or 3 min with fluorescence)	3 min per sample (24 samples - 72 min) * note they list in seconds and it makes it seem minimal	FAST Mode* < 80 sec per sample (24 samples - 32 min) Normal mode 130 sec p/sample (24 samples - 52 min)
Volume	25 µL - 100 µL sample volume 50 µL - 200 µL total well volume	25 µL - 100 µL sample volume 50 µL - 200 µL total well volume	25 µL - 100 µL sample volume 50 µL - 200 µL total well volume	500 µL	200 µL Normal Mode 170 µL FAST Mode*
Suitable for complex primary samples and small immune cells	No	Yes	Yes	No	No
Robotics compatible	Yes	Yes	Yes	Yes (custom configuration eliminates 21 CFR Part 11 compliance due to SW validation)	Yes (custom-quoted, but not implemented yet)
Regulatory	21 CFR Part 11	21 CFR Part 11	21 CFR Part 11	21 CFR Part 11	21 CFR Part 11
Computer	Windows 10	Windows 10	Windows 10	External PC, Win 7 OS	External PC, Win 7 OS

"The ease of use and fast, high-throughput capabilities definitely make the Cellaca® MX the best cell counter on the market, even over the Beckman Vi-CELL which I've used for the past decade."

- Jasmine Ta, Eli Lilly

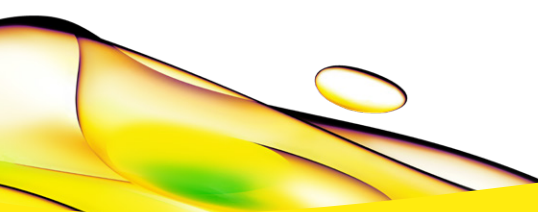
"We were rapidly able to use the Cellaca MX for pipeline PBMC immunology assays integrating both flow cytometry and single-cell genomics platforms. . . . The ability to run a batch of 24 samples in a single plate has been integrated into our data analysis streams. Solid, reliable cell counter with excellent support from the manufacturer."

- Julian Reading, Allen Institute for Immunology

"The Cellaca MX high-throughput cell counter has streamlined the development efforts of our cell therapy product. The instrument combines the accuracy we require for cell viability measurements using the dual live/ dead fluorescent dyes with the automation needed for intensive characterization studies. The time savings for our team with the increased speed and automation will easily justify the cost of the instrument."

- Will Heaton, DiscGenics

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



revvity