

Get your  
counting time  
back.



revvity

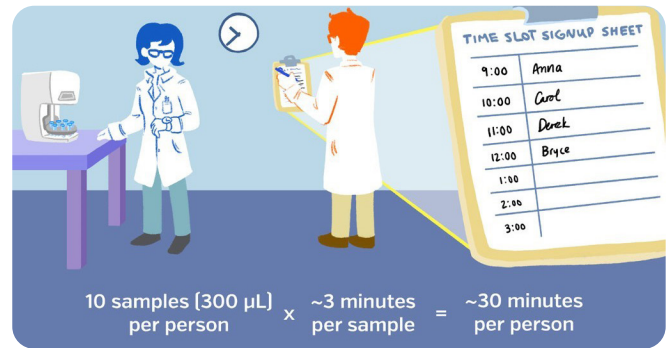
Cellaca<sup>®</sup> MX high-throughput cell counter

# Forget cell counting as you know it

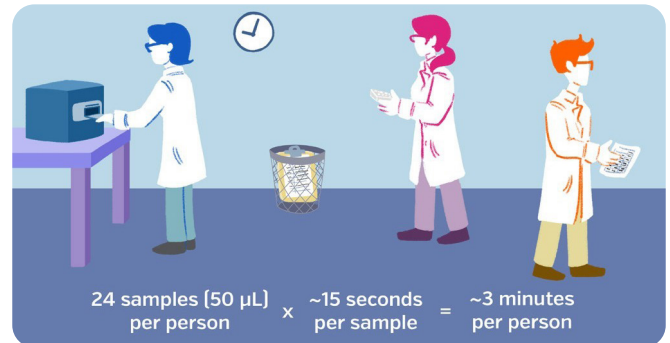
Cellaca MX changes everything. Count 24 samples in 2.5 minutes or less — skip the sign-up sheet and the long waits. Get your all-day, multi-scientist experiments done in 1.5 hours with one pair of hands. Analyze low-volume primary cell samples, and count only the cells you're looking for.

Cellaca MX lets you do high-speed, high-throughput counting.

- Analyze cell lines or primary cells like PBMCs, T cells and tumor digests
- Batch up to 24 samples at a time for cell count, viability by Trypan blue or AO/PI
- Assess sample quality and counts with higher accuracy using image cytometry
- Move cells on to next steps faster to minimize variability in results



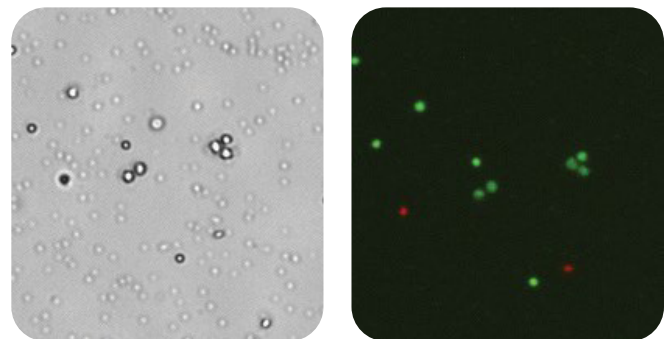
Carousel counter



Cellaca MX

## Count the right cells every time

Cellaca MX gives you the flexibility to run assays that fit your sample and experimental conditions best. Use brightfield to check viability. Switch to fluorescence to get specific details on cell health or analyze heterogeneous populations.<sup>1</sup> Cellaca MX identifies only your cells of interest — tissue debris and red blood cells won't artificially change the count. Move on to next steps knowing the results you'll get there are based on an accurate count.



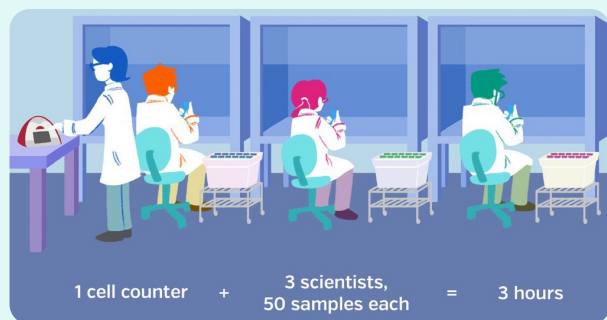
Brightfield and fluorescent images of PBMCs. In brightfield, white blood cells, platelets, and red blood cells are seen. Nuclear fluorescent staining using both green and red channels accurately identifies only the white blood cells.

<sup>1</sup> Accurate measurement of peripheral blood mononuclear cell concentration using image cytometry to eliminate RBC-induced counting error, Chan. et.al, J Immunol Methods, 388:25-32, 2013.

# Get your time back

Save time and resources every time with Cellaca MX.

Experiment: 50 mice, 3 samples per mouse, 150 samples



Single slide counter



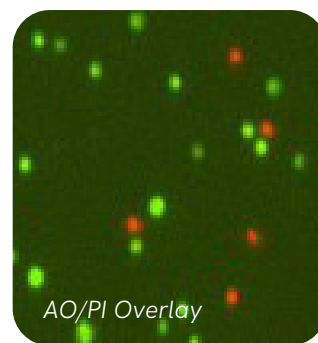
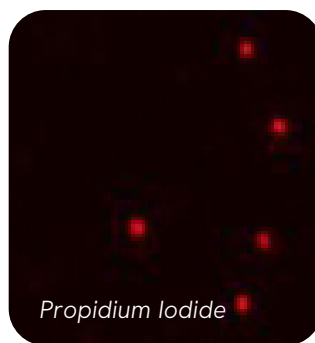
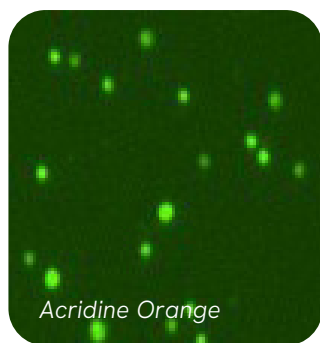
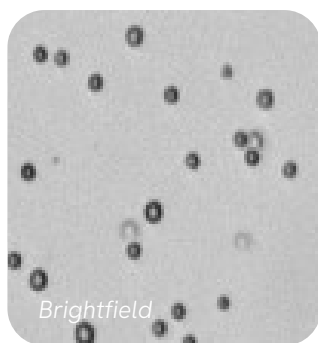
Cellaca MX

Single slide counter	Cellaca MX
4 researchers	1 researcher
1 count/run	24 counts/run
12 hours hands-on time	1.5 hours hands-on time

**Cellaca MX gets the project done 10.5 hours faster, and saves the time of 3 researchers.**

## Analyze primary cells

Getting accurate information on cell health is next to impossible with other systems as they often need 350  $\mu$ L of sample or more. Cellaca MX only needs 25  $\mu$ L, which makes analysis of tumor biopsy tissue, mouse splenocytes, tumor infiltrating lymphocytes (TILs) and more completely possible. You'll also have enough sample left for downstream analyses.



Human PBMC sample stained with Acridine Orange/Propidium iodide (AO/PI) and imaged using brightfield and fluorescence on Cellaca MX.

# Everything you need for your assays

Cellaca MX comes with everything you need to start counting on day one, including software, consumables, and reagents. As your experimental needs change, choose from a wide range of dyes that are tested and validated for use on Cellaca MX.

## System specifications

Description	Specification
Samples per run	24
Counting speed per sample	2-17 sec (assay dependent)
Volume per sample	25-200 $\mu$ L
Imaging methods	Brightfield, Trypan Blue, Fluorescence
Sample types	Cultured cell lines, blood samples, primary cell samples, digested tissue cells, isolated and purified cells
Compatible dyes and fluorophores	Hoechst, DAPI, PE, APC, FITC, CFDA, Calcein AM, 7AAD
Excitation LED	365, 470, 527 and 620 nm
Emission filters (bandpass, center wavelength)	452, 534, 605, 655 and 692 nm
Plate format	SBS Format
Automation ready	Available API for Robotic Integration
Dimensions and weight	13 in x 13 in x 16 in (33 cm x 33 cm x 41 cm) 42 lb (19 kg)

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



revvity

Revvity, Inc.  
940 Winter Street  
Waltham, MA 02451 USA  
[www.revvity.com](http://www.revvity.com)

For a complete listing of our global offices, visit [www.revvity.com](http://www.revvity.com)  
Copyright ©2023, Revvity, Inc. All rights reserved.

1202468

[www.revvity.com](http://www.revvity.com)