



Automated cell counters for single cell sequencing.

Scientists in single cell analysis choose Revvity cell counters

- Low sample volume
- Accurate cell and nuclei counts of cell lines and primary samples
- Reliable counting of isolated nuclei from messy samples
- Automated assessment of cell clumps
- Broad dynamic range for concentration and viability measurement

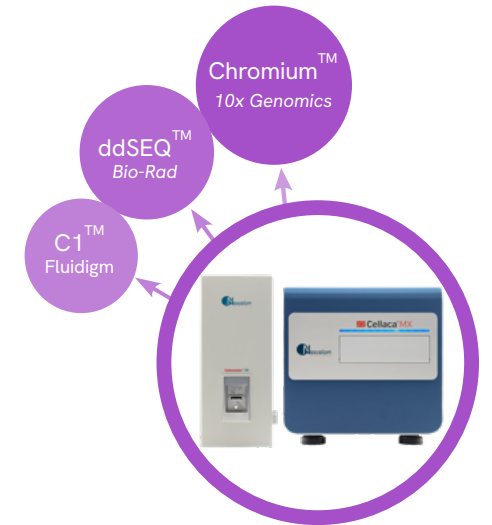
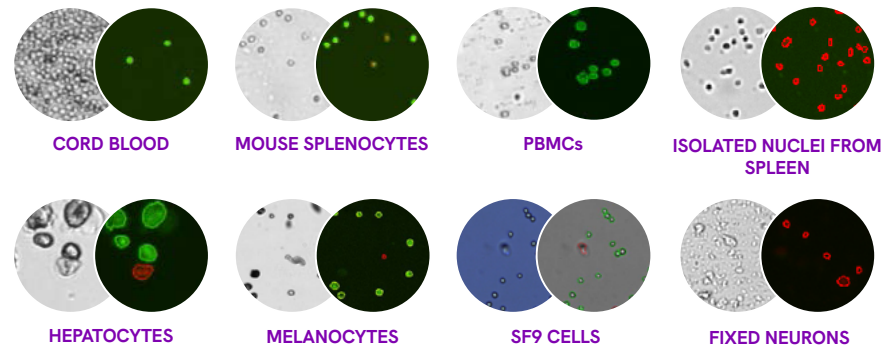
When cell counting results are critical

The field of single-cell analysis is advancing rapidly and continues to provide novel insights into diverse applications, such as gene expression dynamics and disease pathogenesis.

Accurate cell counts are critical in sample preparation for single-cell sequencing.

Accuracy from cell lines to primary samples — including clumpy cells

The Cellometer® K2 and Cellaca® MX cell counters can be customized to handle a variety of cell types (primary cells, tumor digest, insect cells, cell lines, fragile cells, clumpy cells and more) at low or high concentrations.

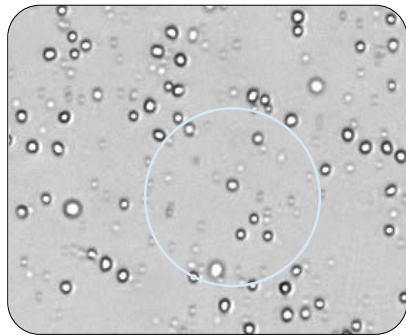


Low sample volume, complete counts

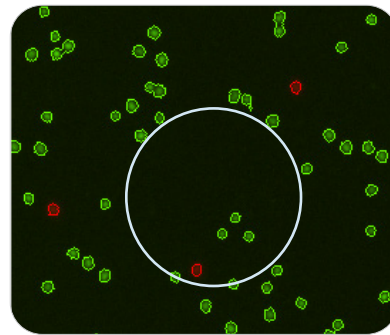
Cell samples in single cell analysis are precious. The Cellometer and Cellaca systems use only 10 µl and 25 µl respectively for accurate counts.

Dual fluorescent staining counts only nucleated cells

The brightfield image on the left shows the combination of nucleated cells, red blood cells, and platelets present in the sample. The red blood cells are not visible in the fluorescent image on the right, only the live (green) and dead (red) nucleated cells are counted.



| Brightfield



| AO/PI

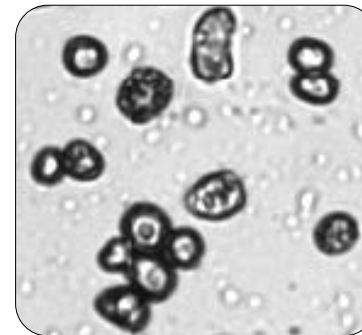
User proven performance

"The Cellometer K2 cell counter helped us significantly with reducing processing time, saving time for cell and nuclei counting, and also more accurate cell counts." - Dr. Christine Cheng, Boston University

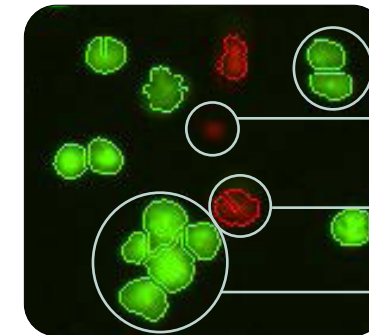
"After testing this cell counter and comparing it to other counting methods, we have seen that the Cellca MX produces tighter/accurate counts and saves us time." - Kimberly Sandy, Pfizer

Specialized algorithm for clumpy cells

The fluorescent image (far right) shows bright green acridine orange (AO) positive hepatocytes declustered by the Cellometer algorithm. Red circled hepatocytes are PI positive (dead) while free nuclei are not counted.



| Brightfield



| AO/PI

Live AO positive

Free nuclei are not counted

Dead, PI positive

Declustered clumpy cells accurately counted

Custom applications

- Viability for Immune Cells[†]
- Viability for "Fragile Cells"^{††}
- RBC+ Platelet Contamination Assay[†]
- QA/QC Pre Viability Using "Do Not Cluster"
- Trypan Blue Count for SF9 Cells
- Viability for Pigmented Cells[†]
- Total Nuclei Count[†]

[†] Using acridine orange / propidium iodide (AO/PI) ^{††} Using Calcein AM / propidium iodide (CAM/PI)

^{*} Using propidium iodide (PI)

Request a free demo!

revvity