



revvity

Cell counting made simple.

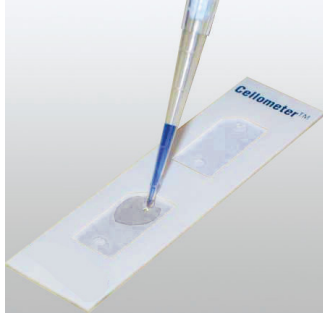
Cellometer® Auto T4 brightfield cell counter

Collect more data, more quickly,
more easily and more accurately than
traditional manual cell counting

- **Cell line analysis:** Obtain cell images, concentration, viability, and cell size in **< 30 seconds**
- **Automated data management:** Archive sample images and auto-save counting and viability results to a secure network
- **Visual confirmation of results:** View cell morphology in real time to ensure only cells of interest are being counted
- **Accurate cell counting:** Pattern-recognition software accurately counts cells in clumps, identifies irregular cell shapes, and eliminates debris from cell counts
- **Cell size analysis:** Overlay cell size histograms to analyze changes in cell size over time. Count specific cell populations based on size



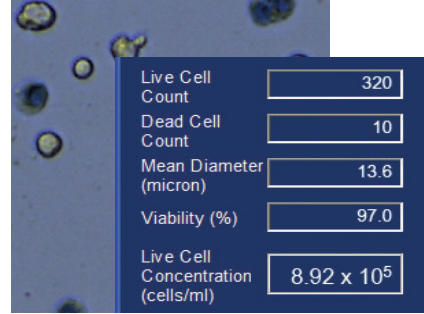
Simple cell counting and Trypan blue viability in seconds!



Pipette 20 μ L of sample



Insert disposable slide



View image & results in <30 seconds

GMP/GLP support module available for access and audit control

- Four different access level controls for user logging
- Audit trail by user
- Autolocking of system (inactivity timeout)
- Password aging reset requirement
- Auto printing option

Cellometer Auto T4 is very intuitive and works great! The best cell counter I have used!

- University of Colorado, Anschutz Medical Campus

The Cellometer Auto T4 is an excellent device that is easy to use and provides very accurate and consistent cell counts. It saves my lab a lot of time. Highly recommended!

- University of Pittsburgh Medical Center

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

