



# We've raised the bar for cell counting.

## Cellometer Ascend automated cell counter

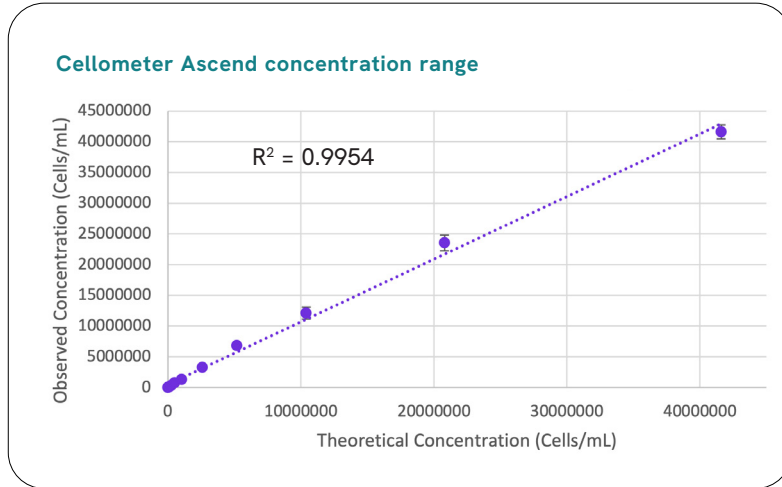
Introducing the Cellometer™ Ascend™ automated cell counter, the newest innovation in our renowned, trusted portfolio of cell counting instruments.

- **Cell counts and viability information:** dual fluorescence and brightfield imaging assist with increasing accuracy for nuclei and cell counting
- **Auto-count upon slide load:** image capture begins immediately allowing rapid size, count, concentration, and viability calculations
- **Count up to eight samples at a time:** choose between a three- or eight-chamber slide
- **Multiple fields of view:** capture up to eight images per sample to increase the accuracy of counts
- **Analysis of various sample types:** including cultured cell lines, PBMCs, primary cells, whole blood, and bone marrow
- **Small sample volume:** as little as 10 µl of cell sample needed
- **Broad concentration range:** count samples from  $2 \times 10^4$  –  $4 \times 10^7$  cells/mL
- **Advanced auto-focus:** slide or image-based focus, adjusted for your particular sample and cell type
- **Bi-directional (x, y) stage movement:** to capture higher sample volume, especially for samples with low cell concentration
- **Favorites feature:** customize experiment settings and save commonly used assays
- **21 CFR Part 11 ready:** optional add-on that includes an audit trail, user access control, and digital signature



## Confidence elevated

The versatile Cellometer Ascend can analyze various cell sample types - both primary cell samples as well as cultured cells, at a wide range of concentrations. In particular, the system has been optimized to analyze low concentration samples for applications such as single cell sequencing and nuclei counting.



Graph of results for cell concentration dynamic range on the Cellometer Ascend. Data set was taken on a concentration series of cultured Jurkat cell line. Samples from  $2 \times 10^4$  -  $4 \times 10^7$  cells/mL can be counted without further dilution. The CV% at each concentration was below 10%.

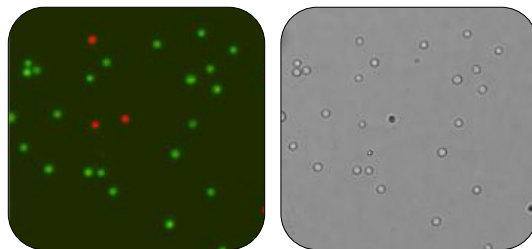
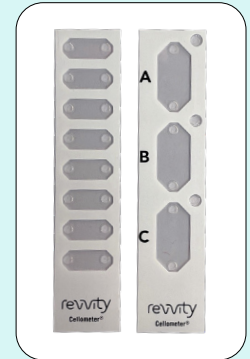


The Cellometer Ascend counter is operated via Matrix software. The 10" screen provides clear images of your cell sample and shows how it was counted.

## Optimized reagents and consumables

The dual-fluorescence stain combination AOPI (acridine orange - propidium iodide) is the preferred method for viability analysis of primary cell samples that may contain debris or red blood cell contamination.

The Cellometer Ascend's single-use counting slides, with either 3 or 8 counting chambers, bring throughput flexibility to the benchtop.



Left: AO/PI  
Right: Brightfield



revvity