

Cellometer™ Ascend

Specification Sheet

- Includes:**
- Cellometer™ Ascend Automated Cell Counter with integrated touchscreen computer and dual-fluorescence optic channels
 - Matrix Software
 - Two Ports on front of instrument (one USB A and one USB C)
 - Three Ports on back of instrument (two USB A and one Ethernet)
 - Power Supply and Power Cord
 - Phone/online applications support during set-up

Imaging Performance: Cell Size 5 - 80 microns
 Conc. Range: 2.0 x 10⁴ - 2.0 x 10⁷ cells/mL 2.5x Magnification / 1um/pixel resolution

Brightfield imaging, fluorescent imaging, and pattern-recognition software to quickly and accurately decluster, identify, and count individual cells.

Instrument Specifications: Weight: 24 lbs (10.5 kg) Dimensions: Width: 9.9" (25 cm)
 Depth: 13" (32.8 cm)
 Height: 18.7" (47.5 cm)

Input to Power Adapter: 100-240 VAC, 50/60 Hz, 1.5A
 Output to Instrument: 12 VDC, 10A

Display: Touchscreen (10.1") with advanced touch capabilities

Fluorescence Optics: **Excitation/Emission: 470 nm/534 nm**
 Example Fluorophores:

- Acridine Orange (AO)
- Calcein AM
- CFDA
- GFP (Green Fluorescent Protein)
- SYTO™ 9
- SYTO™ 13

Excitation/Emission: 531 nm/655 nm
 Example Fluorophores:

- Propidium Iodide (PI)
- Ethidium Bromide (EtBr)