

Cellometer® Auto 2000

Specification Sheet

- Includes:**
- Cellometer Auto 2000 Automated Cell Counter with integrated touchscreen computer and dual-fluorescence optic channels
 - Cellometer Software
 - Two USB 2.0 Ports and one Ethernet Port
 - Power Supply and Power Cord
 - Phone/online applications support during set-up

Imaging Performance: Cell Size 5 - 300* microns
 Conc. Range: 10⁵ - 10⁷ cells/ml

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

* Cellometer CHT4-PD300 slides are required for cells > 80 microns in diameter.

Instrument Specifications: Weight: 26 lbs (11.8 kg) Dimensions: Width: 11.1" (28.2 cm)
 Depth: 12.8" (32.4 cm)
 Height: 13.4" (34.0 cm)

Input to Power Adapter: 100-240 VAC, 50/60 Hz, 1.5A
 Output to Instrument: 12 VDC, 7.5A

Display: 10.4" Touchscreen

Fluorescence Optics: Excitation/Emission: 470 nm/535 nm
 Example Fluorophores:

- Acridine Orange (AO)
- Calcein AM
- CFDA
- SYTO™ 9
- SYTO™ 13

Excitation/Emission: 540 nm/605 nm
 Example Fluorophores:

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