

Cellometer® Auto T4

Specification Sheet

- Includes:**
- Cellometer Auto T4 Automated Cell Counter
 - Cellometer Software
 - Power Supply and Power Cord
 - USB 2.0 Connector Cable
 - Phone/online applications support during set-up

- Available Accessories:**
- PC Laptop

Imaging Performance:

Cell Size 5 – 60 microns (most mammalian cells)
 Conc. Range: 10⁵ – 10⁷ cells/ml

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

Instrument Specifications:

Weight: 10.4 lbs (4.7 kg)

Dimensions: Width: 3.5" (8.9 cm)
 Depth: 4.2" (10.7 cm)
 Height: 12.6" (32.0 cm)

Input to Power Adapter: 100-240 VAC, 50/60 Hz, 0.5A
 Output to Instrument: 12 VDC, 1.5A

- PC/Laptop Minimum Requirements:**
 (If purchasing instrument without PC Laptop)
- Windows 10
 - 2.2 GHz Processor or higher
 - 4 GB RAM (8 GB RAM or higher recommended)
 - Integrated Graphics Card
 - 1024 x 768 pixels Display Resolution
 - USB 2.0 Port