

Cellometer® K2

Specification Sheet

Includes:	<ul style="list-style-type: none"> • Cellometer K2 Automated Cell Counter • Matrix Software or Cellometer Software • Two Standard Fluorescence Optics Modules • Power Supply and Power Cord • USB 2.0 Connector Cable • Phone/online applications support during set-up 	Available Accessories:	<ul style="list-style-type: none"> • PC Laptop* <p>* Recommended for Matrix Software.</p>
------------------	---	-------------------------------	--

Imaging Performance:	<p>Cell Size 5 - 300** microns</p> <p>Conc. Range: 10⁵ - 10⁷ cells/ml</p>	<p>** Cellometer CHT4-PD300 slides are required for cells > 80 microns in diameter.</p>
<p>Brightfield imaging, fluorescent imaging, and pattern-recognition software to quickly and accurately decluster, identify, and count individual cells.</p>		

Instrument Specifications:	<p>Weight: 23 lbs (10.4 kg)</p> <p>Input to Power Adapter: 100-240 VAC, 50/60 Hz, 1.0A</p> <p>Output to Instrument: 12 VDC, 3.34A</p>	<p>Dimensions:</p> <p>Width: 6.0" (15.2 cm)</p> <p>Depth: 8.5" (21.6 cm)</p> <p>Height: 14.0" (35.6 cm)</p>
-----------------------------------	--	---

	<u>Cellometer Software</u>	<u>Matrix Software</u>
PC/Laptop Minimum Requirements: (If purchasing instrument without PC Laptop)	<ul style="list-style-type: none"> • Windows 10 • Intel® i5 (1.6 - 4.40 GHz) Processor • 10 Core • 4 GB RAM • Integrated Graphics Card • 1080p Display Resolution • 500 GB Hard Drive • USB 2.0 Port 	<ul style="list-style-type: none"> • Windows 10 • Intel® i7 (2.10 - 4.80 GHz) Processor • 16 Core, 24 Threads • 32 GB RAM • NVIDIA RTX A2000 8GB GDDR6 • 1080p Display Resolution • 1 TB+ Hard Drive • USB 2.0 Port

Available Fluorescence Optics Modules:	<p>VC-535-403</p> <p>Excitation/Emission: 470nm/535nm</p> <p>Example Fluorophores:</p> <ul style="list-style-type: none"> • Acridine Orange (AO) • CFDA • FITC 	<p>VC-660-503</p> <p>Excitation/Emission: 540nm/660nm</p> <p>Example Fluorophores:</p> <ul style="list-style-type: none"> • Propidium Iodide (PI) • Ethidium Bromide (EtBr) • 7-AAD
---	---	--