

Cellometer® X2

Specification Sheet

Includes:	<ul style="list-style-type: none"> • Cellometer X2 Automated Cell Counter • 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
------------------	--	-------------------------------	---

Imaging Performance:	<p>Cell Size 0.5* – 15 microns</p> <p>Conc. Range: 5 x 10⁵ – 10⁷ cells/ml</p>	<p>* Cellometer CHT4-CD025 slides are required for cells < 20 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>	Dimensions:	<p>Width: 6.0" (15.2 cm)</p> <p>Depth: 8.5" (21.6 cm)</p> <p>Height: 14.0" (35.6 cm)</p>
-----------------------------------	--	--------------------	--

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
--	--

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 • SYTO™ 9 	<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
---	--	--