

Cellometer K2

Specification Sheet

Includes:	<ul style="list-style-type: none"> • Cellometer™ K2 Fluorescent Cell Counter • Cellometer™ Software or Matrix™ 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>
-----------------------------------	--	---

PC Specifications (SW Minimum Requirements):	<p><u>Cellometer Software</u></p> <ul style="list-style-type: none"> • Windows® 11 • Intel® Core™ i5 (1.6 - 4.40 GHz) Processor or higher • 10 Core or higher • 4 GB RAM or higher • Integrated Graphics Card • 1080p Display Resolution • 500 GB Hard Drive • USB 2.0 Port or higher 	<p><u>Matrix Software (Laptop Included)</u></p> <ul style="list-style-type: none"> • Windows 11® LTSC 24H2 or higher • Intel® Core™ i7 (2.10 - 5.30 GHz) Processor or higher • 20 Core, 28 Thread or higher • 32 GB RAM or higher • NVIDIA RTX 2000 Ada 8 GB Graphics Card or equivalent/better • 2 TB Hard Drive (total of C: + D: drives) • USB 2.0 Port or higher
---	---	---

Available Fluorescence Optics Modules:	<p>VC-535-403</p> <p>Excitation/Emission: 475 nm/534 nm</p> <p>Example Fluorophores:</p> <ul style="list-style-type: none"> • Acridine Orange (AO) • CFDA • FITC 	<p>VC-660-503</p> <p>Excitation/Emission: 542 nm/655 nm</p> <p>Example Fluorophores:</p> <ul style="list-style-type: none"> • Propidium Iodide (PI) • Ethidium Bromide (EtBr) • 7-AAD
---	---	--