

# Cellometer® Spectrum

## Specification Sheet

<b>Includes:</b>	<ul style="list-style-type: none"> <li>• Cellometer Spectrum Image Cytometer</li> <li>• Cellometer Software</li> <li>• Two Fluorescence Optics Modules</li> <li>• Power Supply and Power Cord</li> <li>• USB 2.0 Connector Cable</li> <li>• Phone/online applications support during set-up</li> </ul>	<b>Available Accessories:</b>	<ul style="list-style-type: none"> <li>• PC Laptop (includes FCS Express Flow Software from De Novo Software)</li> </ul>
------------------	--	-------------------------------	--

<b>Magnification Selection:</b>	<ul style="list-style-type: none"> <li>• 5x (e.g., most mammalian cells)</li> <li>• 10x (e.g., yeast, platelets, algae)</li> </ul>
---------------------------------	--

<b>Imaging Performance:</b>	<u>5x Magnification</u>	<u>10x Magnification</u>
Cell Size:	5 - 300* microns	0.5 - 15** microns
Conc. Range:	10 <sup>5</sup> - 10 <sup>7</sup> cells/ml	10 <sup>5</sup> - 10 <sup>7</sup> cells/ml

Brightfield imaging, fluorescent 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.

\*\* Cellometer CHT4-CD025 slides are required for cells < 20 microns in diameter.

<b>Instrument Specifications:</b>	<b>Weight:</b> 24.0 lbs (10.9 kg)	<b>Dimensions:</b> Width: 6.0" (15.2 cm) Depth: 8.5" (21.6 cm) Height: 14.0" (35.6 cm)
	<b>Input to Power Adapter:</b> 100-240 VAC, 50/60 Hz, 1.0A	
	<b>Output to Instrument:</b> 12 VDC, 3.34A	

<b>PC/Laptop Minimum Requirements:</b> (If purchasing instrument without PC Laptop)	<ul style="list-style-type: none"> <li>• Windows 10</li> <li>• Intel® i5 (1.6 - 4.40 GHz) Processor</li> <li>• 10 Core</li> <li>• 4 GB RAM</li> <li>• Integrated Graphics Card</li> <li>• 1080p Display Resolution</li> <li>• 500 GB Hard Drive</li> <li>• USB 2.0 Port</li> </ul>
--	--

---

<b>Fluorescence Optics Modules, Standard:</b>	<b>S1-534-470</b> <b>Excitation/Emission: 475 nm/534 nm</b> <i>Example Fluorophores:</i> <ul style="list-style-type: none"><li>• Acridine Orange (AO)</li><li>• GFP (Green Fluorescent Protein)</li><li>• Annexin V-FITC</li></ul>	<b>S1-655-527</b> <b>Excitation/Emission: 525 nm/655 nm</b> <i>Example Fluorophores:</i> <ul style="list-style-type: none"><li>• Propidium Iodide (PI)</li><li>• Ethidium Bromide (EB)</li><li>• 7-AAD</li></ul>
<b>Fluorescence Optics Modules, Available:</b>	<b>S1-605-527</b> <b>Excitation/Emission: 525 nm/605 nm</b> <i>Example Fluorophores:</i> <ul style="list-style-type: none"><li>• Propidium Iodide (PI)</li><li>• Ethidium Bromide (EtBr)</li><li>• R-phycoerythrin (PE)</li><li>• RFP (Red Fluorescent Protein)</li></ul> <b>S1-594-470</b> <b>Excitation/Emission: 475 nm/594 nm</b> <i>Example Fluorophores:</i> <ul style="list-style-type: none"><li>• Algae (Chlorophyll A and B)</li></ul>	<b>S1-692-620</b> <b>Excitation/Emission: 628 nm/692 nm</b> <i>Example Fluorophores:</i> <ul style="list-style-type: none"><li>• Allophycocyanin (APC)</li><li>• Alexa Fluor® 647</li><li>• Cy5™</li><li>• SYTOX™ Red</li></ul> <b>S1-452-365</b> <b>Excitation/Emission: 370 nm/452 nm</b> <i>Example Fluorophores:</i> <ul style="list-style-type: none"><li>• DAPI</li><li>• Hoechst</li><li>• ViaStain™ Dead Cell Nuclear Blue</li></ul>

---