

Cellaca™ MX

Specification Sheet

- Includes:**
- Cellaca™ MX High-throughput Cell Counter
 - Power Supply and Power Cord
 - PC Laptop with Windows OS
 - Matrix Software (pre-installed on PC Laptop)
 - USB 3.0 Connector Cable
 - Phone/online applications support during set-up

Imaging Performance: Cell Size 5 - 80 microns
 Conc. Range: 1.0 x10⁵ - 1.5 x 10⁷ cells/ml (recommended)

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

Instrument Specifications: Weight: 42 lbs (19.1 kg) Dimensions: Width: 13" (33.0 cm)
 Depth: 16" (40.6 cm)
 Height: 13" (33.0 cm)

Input to Power Adapter: 100-240 VAC, 50/60 Hz, 1.5A
 Output to Instrument: 12 VDC, 7.5A

- PC Specifications:** (Laptop Included)
- Windows 10 IoT Enterprise 64 bit or higher
 - Intel® i7 Core™ i7-10750H 2.1 GHz Processor or higher
 - 6 Cores or higher
 - 16 GB RAM or higher
 - NVIDIA RTX A1000 Graphics Card or equivalent/better
 - 1 TB+ Hard Drive (total of C: + D: drives)
 - USB 3.0 Port
 - SQL Server Standard SW License

Fluorescence Optics:**Cellaca MX FL5****Excitation LEDs: 370, 475, 525, and 628 nm****Emission Filters: 452, 534, 605, 655, and 692 nm
(bandpass, center wavelength)***Commonly Used Excitation LED/Emissions Filter Combinations:*

- **365/452:** DAPI, Hoechst
- **470/534:** Annexin V-FITC, AO (Acridine Orange), Calcein AM, Caspase-3 488, CFDA, GFP (Green Fluorescent Protein)
- **527/655:** 7-AAD, PI (Propidium Iodide), RFP (Red Fluorescent Protein)

Cellaca MX FL2**Excitation LEDs: 475 and 527 nm****Emissions Filters: 534 and 655 nm
(bandpass, center wavelength)***Commonly Used Compatible Dyes and Assays:*

- Annexin V-FITC
- AO/PI (Acridine Orange/Propidium Iodide)
- Calcein AM
- Caspase-3 488

Cellaca MX BF**Excitation LEDs: N/A****Emissions Filters: N/A***Commonly Used Compatible Dye:*

- Trypan Blue
-