

# Cellaca<sup>®</sup> PLX

## Specification Sheet

- Includes:**
- Cellaca PLX Image Cytometry System
  - Power Supply and Power Cord
  - PC Laptop with Windows OS
  - Matrix Software (pre-installed on PC Laptop)
  - FCS Express Flow Software from De Novo Software (pre-installed)
  - PLX Slide Holder
  - USB 3.0 Connector Cable
  - Phone/online applications support during set-up

**Imaging Performance:** Cell Size: 5 – 80 microns  
 Conc. Range: 1.0 x10<sup>5</sup> – 1.5 x 10<sup>7</sup> 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
  - Intel<sup>®</sup> i7 (2.10 – 4.80 GHz) Processor
  - 16 Core, 24 Threads
  - 32 GB RAM
  - NVIDIA RTX A2000 8GB GDDR6 Graphics Card
  - 1080p Display Resolution
  - 1 TB+ Hard Drive
  - USB 3.0 Port
  - SQL Server Standard SW License

---

**Fluorescence  
Optics:****Excitation LEDs: 370, 475, 531, and 628 nm****Emission Filters: 452, 534, 605, 655, and 692 nm  
(bandpass, center wavelength)***Commonly Used Fluorescent Dyes and their  
Excitation LED/Emission Filter Combinations:*

- **365/452:** DAPI, Dead Blue, Hoechst
  - **470/534:** Annexin V-FITC, AO (Acridine Orange), Calcein AM, Caspase-3 488, CFDA, Dead Green, GFP (Green Fluorescent Protein), KIRAVIA Blue 520™
  - **531/605:** PE (R-Phycoerythrin), RFP (Red Fluorescent Protein)
  - **531/655:** 7-AAD, PI (Propidium Iodide)
  - **620/692:** APC (Allophycocyanin), RubyDead
-