

Cellaca PLX

Specification Sheet

- Includes:**
- Cellaca™ PLX High-throughput Image Cytometer
 - Power Supply and Power Cord
 - PC Laptop with Windows OS
 - Matrix™ Software (pre-installed on Laptop)
 - FCS Express® from De Novo Software by Dotmatics® (pre-installed on Laptop)
 - 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 x10⁵ – 1.0 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: 44 lbs (20 kgs) 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 (SW Minimum Requirements):**
- 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 3.0 Port
 - SQL Server Standard SW License
- Laptop Included*

Fluorescence Optics:

| Excitation LEDs (nm) | Emission Filters* (nm) |
|----------------------|-------------------------|
| 370 | 452, 534, 605, 655, 692 |
| 475 | 534, 605, 655, 692 |
| 531 | 605, 655, 692 |
| 628 | 692 |

*Bandpass, Center Wavelength

Commonly Used Fluorescent Dyes and their Excitation LED/Emission Filter Combinations:

- **370/452:** DAPI, Dead Blue, Hoechst
 - **475/534:** Annexin V-FITC, AO (Acridine Orange), Calcein AM, Caspase-3 488, CMFDA, 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)
 - **628/692:** APC (Allophycocyanin), RubyDead
-