

Celigo™ with Plate Stacker

Specification Sheet

Includes:

- Celigo™ Image Cytometer
- Power Supply and Power Cord
- PC Laptop with Windows OS
- Celigo Software (pre-installed on Laptop)
- USB 2.0 and USB 3.0 Connector Cables
- Automated Plate Stacker*
- Satellite Workstation
- FCS Express Image Software by De Novo Software

* Automated Plate Stacker includes Power Supply/Power Cord, Serial Cable for computer connection, Compressor (with Power Supply/Power Cord), and Airline.

Imaging

Cell Size 5 - 80 microns

Performance:

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:

(without Desktop)

Weight (Celigo only): 117 lbs (53 kg)

Weight (Stacker only): 47 lbs (21 kg)

Electrical Power: 100-240 VAC, 50/60 Hz, 600 W, 8.4A (max). Five electrical outlets needed for instrument and stacker

Fuse Rating: 10A SLO-BLO, 250 VAC 3AB (Celigo only)

Dimensions: Width: 20" (50.8 cm)
(Celigo only) Height: 17" (43.5 cm)
Depth: 25" (63.5 cm)*

* Depth: 29.5" (72.39 cm) if plate holder is extended

Dimensions: Width: 44" (112 cm)
(Celigo and Height: 40" (102 cm)
Plate Stacker) Depth: 25" (63.5 cm)*

Temperature and Humidity Requirements:

Operating Temperature: 15°C to 25°C

Operating Humidity: 10% to 90% RH, non-condensing

Shipping and Storage Temperature: -18°C to 65°C

Shipping and Storage Humidity: 10% to 90% RH, non-condensing

Manufacturer

Celigo

Revvity, Inc.
360 Merrimack Street, Suite 200
Lawrence, MA 01843 USA

Plate Stacker

Kbiosystems Limited
Units 5 to 10 Paycocke Close
Basildon, Essex, UK, SS143HS

- PC Specifications:** (Desktop Included)
- Model: Dell Precision Tower 3660 XE
 - OS: Windows 10 IoT Enterprise LTSC
 - CPU: Intel® i7-12700 25 MB Cache, 12 Core
 - RAM: 64 GB DDR5 ECC
 - Boot SSD: 512 GB 2.5" SATA SSD
 - Data Storage: 4TB 3.5" SATA SSD
 - GPU: NVIDIA T1000 4GB
 - Monitor: Dell 24" 1920 x1080 Full HD
 - Network: Additional 1Gbit PCI-e NIC
 - USB 3.0 Port (Quantity of 5 Minimum)
 - Microsoft SQL Server 2019 SW License
 - Celigo USB License Dongle
 - 3 Years Dell ProSupport Warranty

Configurations:

	<i>Optics</i>	<i>Applications</i>
BF	Brightfield (BF) Counting	Basic Application Suite
Celigo 4B Config	BF, 377/470 (Blue), 483/536 (Green), 531/629 (Red)	Basic Application Suite
Celigo 4C Config	BF, 377/470 (Blue), 483/536 (Green), 531/629 (Red)	Classic Application Suite
Celigo 5C Config	BF, 377/470 (Blue), 483/536 (Green), 531/629 (Red), 628/688 (Far Red)	Classic Application Suite Celigo Virology Suite – Requires additional purchase of Virology Suite License

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

- **377/470:** DAPI, Total and Dead Cell Nuclear Blue, Hoechst
- **483/536:** AlexaFluor® 488, Annexin V-FITC, AO (Acridine Orange), Calcein AM, Caspase-3 488, CFSE, CMFDA, GFP (Green Fluorescent Protein), Total and Dead Cell Nuclear Green
- **531/629:** AlexaFluor® 568, PE (R-Phycoerythrin), PI (Propidium Iodide), RFP (Red Fluorescent Protein), Total and Dead Cell Nuclear Red
- **628/688:** AlexaFluor® 647, Draq5®, Total and Dead Cell Nuclear Far Red

Basic Application Suite includes:

Cell Counting

- Direct Cell Counting
- Cell Viability
- Dead + Total
- Live + Dead
- Live + Dead + Total
- Live + Total

Confluence

Confluence 1

- Confluence 1+2
- Confluence 1+2+3
- Confluence 1+2+3+4
- Confluence 1+2+3+4+5

Colony

- Colony 1
- Colony 1+2 (Merge)
- Colony 1+2+3 (Merge)
- Colony 1+2+3+4 (Merge)
- Colony 1+2+3+4+5 (Merge)
- Colony 1+2+3+4 + Mask
- Colony 1+2+3 + Mask
- Colony 1+2 + Mask
- Colony 1 + Mask
- Single Colony Verification

Basic Application Suite includes (Continued):

Expression Analysis

- Target 1
- Target 1+2
- Target 1+2 (Merge)
- Target 1+2+3
- Target 1+2+3 (Merge)
- Target 1+2+3+4
- Target 1+2+3+4 (Merge)
- Target 1+2+3+4+5
- Target 1+2+3+4+5 (Merge)
- Target 1+2+3+4 + Mask
- Target 1+2+3 + Mask
- Target 1+2 + Mask Target 1 + Mask

Tumorsphere

- Tumorsphere 1
- Tumorsphere 1+2+3+4 + Mask
- Tumorsphere 1+2+3 + Mask
- Tumorsphere 1+2 + Mask
- Tumorsphere 1 + Mask

Classic Application Suite includes:

All Basic Applications

Cell Secretion

- Cell + Secretion

DNA Synthesis

- Synthesis + Total

Embryoid Body

Invasion and Migration

- Tumorsphere Invasion
- Tumorsphere Migration
- Wound Healing

Plate Reader

- Whole Well 1
- Whole Well 1+2
- Whole Well 1+2+3
- Whole Well 1+2+3+4
- Whole Well 1+2+3+4+5

PS Externalization

- PS Ext + Dead + Total
- PS Ext + Total

Spheroid Analysis

- Zonal Tumorsphere 1+2 + Mask
- Zonal Tumorsphere 1 + Mask

Virology Suite includes:

(Requires an additional purchase; not included in Basic or Classic Application Suites)

Foci Counting

Inverted Plaque

Plaque Counting

Viral Titration

Technical Diagram:

