

Celigo™

Specification Sheet

Includes:	<ul style="list-style-type: none"> • Celigo™ Image Cytometer • Power Supply and Power Cord • PC Desktop with Windows OS • Celigo Software (Pre-installed on Laptop) • USB 2.0 and USB 3.0 Connector Cables 	Available Accessories:	<ul style="list-style-type: none"> • Automated Plate Stacker • Satellite Workstation • FCS Express™ Image Software by De Novo Software
------------------	---	-------------------------------	---

Imaging Performance:	Cell Size	5 - 80 microns
	Conc. Range:	1.0 x10 ⁵ - 1.5 x 10 ⁷ 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:	117 lbs (53 kg)	Dimensions:	Width: 20" (50.8 cm) Height: 17" (43.5 cm) Depth: 25" (63.5 cm)*
	Electrical Power:	100-240 VAC, 50/60 Hz, 600 W, 8.4A (max)		
	Fuse Rating:	10A SLO-BLO, 250 VAC 3AB		* Depth: 29.5" (72.39 cm) if plate holder is extended.

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

PC Specifications: (Desktop Included)	<ul style="list-style-type: none"> • 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) 	<ul style="list-style-type: none"> • 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