

Incucyte® S3 Live-Cell Analysis System



Operating Systems:	64-bit Windows 10			Network Connection:	Ethernet 10 100 1000 Mbps		
Objectives:	4X 10X 20X			Power:	100-240 VAC 50 60 Hz 3.5 Amps max @ 120 V 1.75 Amps max @240 V		
Image Resolution:	4X: 2.82 µm/pixel 10X: 1.24 µm/pixel 20X: 0.62 µm/pixel			Controller Dimensions (H X W X D):	5.5" X 17.0" X 22.7" 14.0 cm X 43.2 cm X 57.7 cm		
Exported Image Formats:	JPEG, PNG, TIFF, MetaMorph Multi Dimensional File v2.0, Meta Imaging Series TIFF (Single Multi-Plane) v1			Controller Weight:	41 lbs 18.6 kg		
Exported Movie Formats:	WMV AVI MPEG-4			Microscope Dimensions (H X W X D):	12.6" X 17.7" X 18.8" 32.0 cm X 45.0 cm X 47.8 cm		
Fluorescence Excitation Emission Dual Color Module 4614:	Channel	Excitation	Emission	Controller Operating Environmental Conditions:	0° to 33° C 5% to 90% RH Non-Condensing		
	Green	441-481 nM	503-544 nM	Microscope Operating Environmental Conditions:	0° to 42° C 5% to 95% RH Non-Condensing		
	Red	567-607 nM	622-704 nM				
* Internal Data Storage:	27.3 TB RAID Hard Drive 64 GB RAM						

* Reflects specifications for instruments manufactured after May 31, 2021

Learn more at

www.sartorius.com/incucyte

North America

Phone +1 734 769 1600, ext. 1

Email orders.US07@sartorius.com

Europe

Phone +44 (0) 1763 227400

Email euorders.UK03@sartorius.com

Japan

Phone +81 3 6478 5202

Email orders.US07@sartorius.com

China

Phone +86 21 6878 2300

**Rest of Asia Pacific and other
countries around the world:**

Phone +65 6872 3966



For further contacts, visit

www.sartorius.com