

Incucyte[®] SX5 Live-Cell Analysis System



Operating System:	64-bit Windows 10
Included Objectives:	4X / 10X / 20X
Image Resolution:	4X: 2.82 um/pixel 10X: 1.24 um/pixel 20X: 0.62 um/pixel
Exported Image Formats:	JPEG, PNG, TIFF, MetaMorph Multi Dimensional File v2.0, Meta Imaging Series TIFF (Single / MultiPlane) v1
Exported Movie Formats:	WMV AVI MPEG-4
Internal Data Storage:	27.3 TByte RAID Hard Drive 64 GByte RAM
Network Connection:	Ethernet 10/100/1000 Mbps

Power:	100-240 VAC 50/60Hz 3.5 Amps max @ 120V 1.75 Amps max @ 240V
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
Controller Weight:	36.5 lbs 16.6 Kg
Microscope Dimensions (H X W X D):	13.43" X 17.7" X 18.7" 34.1 cm X 45.0 cm X 47.4 cm
Microscope Weight:	51 lbs 23.1 Kg
Controller Operating Environmental Conditions:	0° to 33° C 5% to 90% RH Non-Condensing
Microscope Operating Environmental Conditions:	0° to 42° C 5% to 95% RH Non-Condensing
Incubator Specification:	200L minimum internal volume

Optical Modules

	Channel	Excitation	Emission
G/O/NIR (standard)	Green Orange NIR	453-485 nm 546-568 nm 648-674 nm	494-533 nm 576-639 nm 685-756 nm
G/R (optional: Cat. No. 4832)	Green Red	441-481 nm 567-607 nm	503-544 nm 622-704 nm
Metabolism (optional: Cat. No. 4820)	ATP	473-498 nm 524-550 nm	565-591 nm
	Orange NIR	524-550 nm 648-674 nm	565-591 nm 685-756 nm

Incucyte® SX5 Application Specifications

	Application	G/O/NIR (Included with SX5)	Metabolism (Optional Cat. No. 4820)	G/R (Optional Cat. No. 4832)
Cell Health and Proliferation	Confluence	■	■	■
	Cell-by-Cell Analysis Adherent*	■	■	■
	Cell-By-Cell Analysis Non-adherent*	■	■	■
	Apoptosis†	■	■	■
	Cytotoxicity	■		■
	Single Spheroid*	■	■	■
	Multi-Spheroid*	■	■	■
	Cell Cycle†	■		■
	ATP Metabolism*		■	
	Mitochondrial Membrane Potential*	■	■	
Cell Function	Immune Cell Killing†	■	■	■
	Antibody Internalization†	■	■	■
	Phagocytosis	■	■	■
	NETosis	■		■
	Neuronal Activity*	■		
	Angiogenesis*	■		■
	Live-Cell Immunocytochemistry†	■		■
Cell Movement and Morphology	Chemotaxis*	■	■	■
	Scratch Wound*	■	■	■
	Immune Cell Clustering and Proliferation†	■	■	■
	Neurite Outgrowth*	■		■
Cell Culture Workflows	Cell Culture QC	■	■	■
	Dilution Cloning	■	■	■
	Whole Well	■	■	■
	Reporter Genes†	■	■	■
	Stem Cell Monitoring and Reprogramming	■	■	■
	Transfection Efficiency†	■		■

* Requires optional software module

† Additional software available for enhanced analysis

Sales and Service Contacts

For further contacts, visit
www.sartorius.com/Incucyte

 facebook.com/essenbioscience/

 linkedin.com/company/essen-bioscience/

 twitter.com/EsSENBioScience

North America

Phone: +1734 769 1600, ext. 1

E-Mail: orders.US07@sartorius.com

Europe

Phone +44 (0) 1763 227400

E-Mail: euorders.UK03@sartorius.com

Asia-Pacific

Phone: +81 3 6478 5202

E-Mail: orders.US07@sartorius.com