

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

Incucyte® SX1 Supported Applications and Objectives

	Application	Compatible Objectives		
		4X	10X	20X
Cell Health & Proliferation	Confluence	■	■	■
	Cell-by-Cell Analysis Adherent*		■	■
	Cell-by-Cell Analysis Non-Adherent*			■
	Apoptosis	■	■	■
	Cytotoxicity	■	■	■
	Label-Free Viability (AI Cell Health Analysis)*		■	■
	Single Spheroid*	■	■	
	Multi Spheroid*	■		
	Organoid QC*	■		
	Organoid Assay*	■		
Cell Function	Cell Cycle†	■	■	■
	Immune Cell Killing	■	■	■
	Antibody Internalization†		■	■
	Phagocytosis		■	■
	NETosis	■	■	■
	Angiogenesis*	■	■	
	Live-Cell Immunocytochemistry†		■	■
Cell Movement and Morphology	Kinase Akt Assay	■	■	■
	Chemotaxis*		■	
	Scratch Wound*	■	■	■
	Immune Cell Clustering & Proliferation†	■	■	■
	Neurite Outgrowth*		■	■
Cell Culture Workflows	Advanced Label-Free Classification Analysis*		■	■
	Cell Culture QC	■	■	■
	Dilution Cloning	■		
	Whole Well	■		
	Reporter Genes†	■	■	■
	Stem Cell Monitoring & Reprogramming	■		
	Transfection Efficiency†	■	■	■

* Requires software module

† Additional software available for enhanced analysis

North America

Sartorius Corporation
300 West Morgan Road
Ann Arbor, Michigan 48108
USA
Phone +1734 769 1600
E-Mail: ips.opm.na@sartorius.com

Europe

Sartorius UK Ltd.
Longmead Business Centre
Blenheim Road
Epsom
Surrey, KT19 9QQ
United Kingdom
Phone +44 1763 22740
E-Mail: euorders.UK03@sartorius.com

Asia Pacific

Sartorius Singapore Pte. Ltd.
30 Pasir Panjang Road,
#06-32, Mapletree Business City
Singapore 117440
Phone : + 65 6872 3966
Email: sartoriusap@sartorius.com



Find out more: www.sartorius.com/incucyte



For questions, email: AskAScientist@sartorius.com