

Incucyte®

Empower Live-Cell Analysis
Inside Your Incubator



Product Information

Analyze your cells for days, weeks or even months as they sit stationary in the stable environment of your tissue culture incubator. With Incucyte's user-friendly interface and robust instrument portfolio, any cell biologist can gain dynamic insights into the health, morphology, movement and function of their cell models.

Incucyte® offers a range of systems to meet your needs and budget. From our workhorse, the Incucyte® S3, to our economical SX1, to the newest member, the frontier-breaking SX5, we deliver real-time insights to simplify progress for labs of any size.

Key Incucyte[®] Specifications

	Application	SX5	S3	SX1
Capacity and Vesselware Compatibility (1) 3 microplate trays included for SX5 and S3. Purchase separately for SX1. (2) Trays available for purchase separately. See our latest vessel compatibility list for full details.	Interchangeable Trays	3	3	1
	Microplate Capacity	6	6	2
	Microplate Compatibility	6, 12, 24, 48, 96, & 384 well (1)		
	Flask Compatibility	T-25, T-50, T-75, T-100, T-175, T-225 (2)		
	Other Labware Compatibility	35mm, 60mm, 100mm, 150mm dishes, and chamber slides (2)		
Optical System	Total Fluorescence Channels	Up to 5	2	2
	<ul style="list-style-type: none"> ▪ Green/Orange/NIR Optical Module 	included	N/A	N/A
	<ul style="list-style-type: none"> ▪ Green/Red Optical Module 	optional	included	included
	<ul style="list-style-type: none"> ▪ Metabolism Optical Module 	optional	N/A	N/A
	Label-Free, HD Phase Imaging	■	■	■
	Objectives	4X, 10X, 20X included		4X, 10X, 20X purchase separately
	Objective Automation	Automated turret		Manual setting in software to switch
Multi-user Support	Remote Network Capability	■	■	■
	Unlimited, Free User Licenses	■	■	■
	Independent Vessel Scheduling	■	■	N/A
	Multi-objective Scheduling	■	■	N/A
Incubation	All systems leverage the stable incubation of a tissue culture incubator			
	Incubator size	>200 L required	>200 L recommended	>200 L recommended
	Incubation to 42 °C	■	■	■
Computing and Storage	Operating System	64-bit Windows 10		
	Controller Data Storage	27.3 TB	27.3 TB*	27.3 TB*
	Expandable RAID Storage	Add an additional 32.7 TB of storage with Incustore [®]		
	Controller Memory	64GB	64 GB*	64 GB*

* Reflects specifications for instruments manufactured after May 31, 2021.

Key Incucyte® Specifications (Continued)

	Application	SX5	S3	SX1
Software Module Compatibility	Incucyte® Cell-by-Cell Analysis	■	■	■
All software modules sold separately.	Incucyte® Spheroid Analysis	■	■	■
	Incucyte® Scratch Wound Analysis	■	■	■
	Incucyte® Chemotaxis Analysis	■	■	■
	Incucyte® Neurotrack Analysis	■	■	■
	Incucyte® Neuronal Activity Analysis	■		
	Incucyte® Angiogenesis Analysis	■	■	■
	Incucyte® Organoid Analysis Software Module	■	■	■
	Incucyte® Advanced Label-Free Classification Analysis Software Module	■	■	■
	Incucyte® AI Cell Health Analysis Software Module	■	■	■
	Incucyte® ATP Analysis	■		

Incucyte® Application Compatibility

	Application	SX5	S3	SX1
Cell Health and Proliferation	Confluence	■	■	■
	Cell-by-Cell Analysis Adherent**	■	■	■
	Cell-By-Cell Analysis Non-adherent**	■	■	■
	Apoptosis†	■	■	■
	Cytotoxicity	■	■	■
	Single Spheroid**	■	■	■
	Multi-Spheroid**	■	■	■
	Organoid**	■	■	■
	Cell Cycle†	■	■	■
	Label-Free Viability (AI Cell Health Analysis)**	■	■	■
	ATP Metabolism**	■		
	Mitochondrial Membrane Potential**	■		

Incucyte[®] Application Compatibility (Continued)

	Application	SX5	S3	SX1
Cell Function	Immune Cell Killing [†]	■	■	■
	Antibody Internalization [†]	■	■	■
	Phagocytosis	■	■	■
	NETosis	■	■	■
	Neuronal Activity**	■		
	Angiogenesis**	■	■	■
	Kinase Akt Assay	■	■	■
	Live-Cell Immunocytochemistry [†]	■	■	■
Cell Movement and Morphology	Chemotaxis**	■	■	■
	Scratch Wound**	■	■	■
	Immune Cell Clustering and Proliferation [†]	■	■	■
	Advanced Label-Free Classification Analysis**	■	■	■
	Neurite Outgrowth**	■	■	■
Cell Culture Workflows	Cell Culture QC	■	■	■
	Dilution Cloning	■	■	■
	Whole Well	■	■	■
	Reporter Genes [†]	■	■	■
	Stem Cell Monitoring and Reprogramming	■	■	■
	Transfection Efficiency [†]	■	■	■

** Requires software module

[†] Additional software available for enhanced analysis

For more information, visit
www.sartorius.com/incucyte

North America

Sartorius Corporation
300 West Morgan Road
Ann Arbor, Michigan 48108
USA
Phone +1 734 769 1600
Email: lps.opm.na@sartorius.com

Europe

Sartorius UK Ltd.
Longmead Business Centre
Blenheim Road
Epsom
Surrey, KT19 9QQ
United Kingdom
Phone +44 1763 227400
Email:
orderhandling.lps.ne@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