

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.

From our workhorse, the Incucyte® S3, to our economical SX1, to the versatile SX5, and to the new, groundbreaking CX3 with spinning disk confocal capabilities for 3D models, we deliver real-time insights to simplify progress for labs of any size.

Key Incucyte® Specifications

		CX3	SX5	S3	SX1
Capacity and Vesselware Compatibility (1) 3 microplate trays included for SX5, S3 and CX3. Purchase separately for SX1. (2) Trays available for purchase separately. See our latest vessel compatibility list for full details.	Interchangeable Trays	3	3	3	1
	Microplate Capacity	6	6	6	2
	Microplate Compatibility	6, 12, 24, 48, 96, and 384 well (1)			
	Flask Compatibility	T-25, T-50, T-75, T-100, T-175, T-225 (2)			
	Other Labware Compatibility	35 mm, 60 mm, 100 mm, 150 mm dishes, and chamber slides (2)			
Optical System	Total Fluorescence Channels	3	Up to 5	2	2
	• Green/Orange/NIR Optical Module	Laser module	included	N/A	N/A
	• Green/Red Optical Module	N/A	optional	included	included
	• Metabolism Optical Module	N/A	optional	N/A	N/A
	Label-Free, HD Phase Imaging	■	■	■	■
	Objectives	4x, 10x, 20x included			4x, 10x, or 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	> 250 L required	>200 L required	>200 L recommended	>200 L recommended
	Incubation to 42 °C	■	■	■	■
Computing and Storage	Operating System	Windows 11 IoT Enterprise LTSC	64-bit Windows 10		
	Controller Data Storage	43.6 TB	27.3 TB	16.4 TB	16.4 TB
	Expandable RAID Storage available for purchase separately	87.2 TB	32.7 TB	32.7 TB	32.7 TB
	Controller Memory	64 GB	64 GB	48 GB	48 GB
Software Module Compatibility All software modules sold separately.	Incucyte® AI Cell Health	■	■	■	■
	Incucyte® Organoid Analysis	■*	■	■	■
	Incucyte® Cell-by-Cell Analysis	■	■	■	■
	Incucyte® Spheroid Analysis	■	■	■	■
	Incucyte® Scratch Wound Analysis	■*	■	■	■
	Incucyte® Chemotaxis Analysis	■*	■	■	■
	Incucyte® Neurotrack Analysis	■*	■	■	■
	Incucyte® Neuronal Activity Analysis	■*	■		
	Incucyte® ATP Analysis		■		
	Incucyte® Angiogenesis Analysis	■*	■	■	■

*Coming soon with future software iterations

Incucyte® Reagents, Consumables and Software

Sartorius provides a comprehensive selection of reagents, consumables, and software designed to meet the specific needs of your applications, enabling long-term, kinetic live-cell imaging and analysis. With optimized, turnkey solutions, you can easily obtain precise and

reliable measurements of cell health, movement, and function, seamlessly linking phenotype and function to pathological processes. Accessing biologically relevant information has never been simpler!

Find Your Solution to Meet Your Application Needs:

Cell Health

Proliferation

Viability

Apoptosis

Cytotoxicity

Mitochondrial Membrane Potential

ATP Metabolism

Cell Function

Immune Cell Killing

Antibody Internalization

Live-Cell Immunocytochemistry

Phagocytosis

NETosis

Neuronal Activity

Cell Movement and Morphology

Chemotaxis Migration and Invasion

Scratch Wound Migration and Invasion

Immune Cell Activation and Proliferation

Neurite Outgrowth

Assays for 3D Models

Spheroid Growth

Spheroid Invasion

Spheroid Immune Cell Killing

Organoid Assay

Cell Monitoring and Workflows

Cell Culture QC

Organoid QC

EV Uptake

Key Advantages

- Gather valuable data using sensitive, non-perturbing reagents for long-term, live-cell studies.
- Improve productivity with lab-tested protocols and intuitive, purpose-built software.
- Capture data-rich insights from every sample using fluorescent reagent combinations.
- Support visualization and automate cell movement studies with Incucyte® consumables.

Germany

Sartorius Lab Instruments GmbH & Co. KG
Otto-Brenner-Straße 20
37079 Göttingen
Phone +49 551 308 0

USA

Sartorius Corporation
3874 Research Park Drive
Ann Arbor, MI 48108
Phone +1 734 769 1600

Orders EMEA

Email: orderhandling.lps-ne@sartorius.com

Orders USA

Email: LPS.OPM.NA@sartorius.com

 **Find out more:** [sartorius.com/incucyte](https://www.sartorius.com/incucyte)

 **For questions, email:** AskAScientist@sartorius.com