

# Introducing the Incucyte® SX5 Live-Cell Analysis System

More Colors. More Insights. More Possibilities.

## Leading the Way With Living Cells

See more information in every sample and explore more applications. Leverage up to 5 different fluorescence channels, up to 3 at a time, for long term timelapse experiments.

## Go Where Your Research Takes You

Study complex immune-tumor cell interactions, synaptic activity in neuronal co-cultures, metabolism in cancer cells, and much more—with a single platform.

## Protect Your Cells

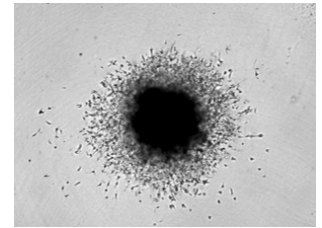
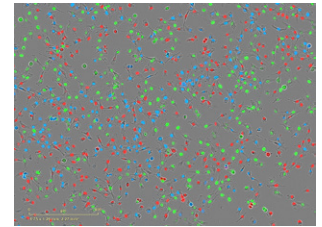
Patent-pending 3-color optical module includes a long wavelength, low phototoxicity Near IR channel and reagents designed for long term live-cell experiments.

## Improve Productivity

Enjoy walk-away convenience as images are automatically acquired and analyzed in microplate format, up to six in parallel.



The Incucyte SX5 Live-Cell Analysis System offers more channels, more reagents and more purpose-built software for more applications—allowing you to derive deeper, physiologically relevant information about your cells. Never miss powerful insights again, with the Incucyte SX5 Live-Cell Analysis System, Software, Reagents, and Consumables.



## Dedicated to Living Cells

- Up to 5 different fluorescence channel options
- Multiplex HD Phase with up to 3 fluorescence channels at a time (Green/Orange/Near IR)
- 4x, 10X, and 20X lenses on an automated turret
- Purpose-built software modules, reagents and consumables for turnkey applications

## Support for Multiple Users

- Support for 3 interchangeable vessel trays and over 600 vessels, up to 6 microplates in parallel
- Remote, networked access with unlimited, free licenses

Learn more at

[www.sartorius.com/incucyte](http://www.sartorius.com/incucyte)

E-mail: [orders.US07@sartorius.com](mailto:orders.US07@sartorius.com)

**North America:** +1 734 769 1600, ext. 3

**Europe:** +44 7515 947101

**APAC:** +81 3 6478 5202

See how the Incucyte is driving research forward at

[www.essenbioscience.com/publications](http://www.essenbioscience.com/publications)

## See What You Can Do With the Incucyte Live-Cell Analysis System!

### Cell Health & Proliferation

Proliferation & Cell Counting  
Cell Cycle  
Apoptosis  
Cytotoxicity  
Viability  
Mitochondrial Membrane Potential NEW!  
ATP Metabolism NEW!

### Cell Function

Immune Cell Killing  
Antibody Internalization  
Immunocytochemistry  
Phagocytosis  
Neurite Dynamics  
Neuronal Activity  
Angiogenesis

### 3D Cell Models

Spheroid Growth & Viability  
Spheroid Invasion

### Cell Movement & Morphology

Chemotaxis Migration & Invasion  
Scratch Wound Migration & Invasion

Specifications subject to change without notice.

© 2020, Essen BioScience, Inc., part of the Sartorius Group. All Rights Reserved. Incucyte and all names of Incucyte products are registered trademarks and the property of Essen BioScience unless otherwise specified. Incucyte is a Sartorius brand. Printed in the EU or US on paper bleached without chlorine. Publication No.: 8000-0737-A00 Version 1 | 2020 | 04