









## Incucyte® Live-Cell Analysis System Reagents and Consumables

Use Incucyte® Reagents and Consumables with your live-cell analysis system for a turnkey solution to a broad range of cell health, movement, function, and other applications. Improve efficiency while maintaining cell

health with Incucyte Reagents that are non-perturbing to cell proliferation and morphology. Find your solution to meet your application needs!

### Apoptosis & Cytotoxicity

Product	Color	Quantity	Compatibility	Cat. No.
Incucyte® Annexin V Green Dye		One vial: 100 tests	SX1, S3, SX5	4642
Incucyte® Annexin V Red Dye		One vial: 100 tests	SX1, S3, SX5 (Green/Red Optical Module)	4641
Incucyte® Annexin V NIR Dye		One vial: 100 tests	S3 for Neuroscience, SX5	4768
Incucyte® Annexin V Orange Dye		One vial: 100 tests	S3 for Neuroscience, SX5	4759
Incucyte® Caspase-3/7 Green Dye		One vial: 20 µL	SX1, S3, SX5	4440
Incucyte® Caspase-3/7 Red Dye		One vial: 20 µL	SX1, S3, SX5 (Green/Red Optical Module)	4704
Incucyte® Cytotox Green Dye		Five vials: 5 µL each	SX1, S3, SX5	4632
Incucyte® Cytotox Red Dye		Five vials: 5 µL each	SX1, S3, SX5 (Green/Red Optical Module)	4633

### Proliferation & Cell Cycle

Product	Color	Quantity	Compatibility	Cat. No.
Incucyte® Nuclight Green Lentivirus (puro)		One vial: 0.2 mL	SX1, S3, SX5	4624
		One vial: 0.6 mL		4475
Incucyte® Nuclight Green Lentivirus (bleo)		One vial: 0.2 mL	SX1, S3, SX5	4626
		One vial: 0.6 mL		4477
Incucyte® Nuclight Red Lentivirus (puro)		One vial: 0.2 mL	SX1, S3, SX5 (Green/Red Optical Module)	4625
		One vial: 0.6 mL		4476
Incucyte® Nuclight Red Lentivirus (bleo)		One vial: 0.2 mL	SX1, S3, SX5 (Green/Red Optical Module)	4627
		One vial: 0.6 mL		4478
Incucyte® Nuclight Orange Lentivirus (puro)		One vial: 0.2 mL	S3 for Neuroscience, SX5	4771
Incucyte® Nuclight NIR Lentivirus (puro)		One vial: 0.2 mL	S3 for Neuroscience, SX5	4805
Incucyte® Nuclight Rapid Red Dye		One vial: 50 µL	SX1, S3, SX5 (Green/Red Optical Module)	4717
Incucyte® Nuclight Rapid NIR Dye		One vial: 50 µL	S3 for Neuroscience, SX5	4804
Incucyte® Cell Cycle Green/Red Lentivirus (puro)		One vial: 0.6 mL	SX1, S3, SX5 (Green/Red Optical Module)	4779
Incucyte® Cell Cycle Green/Orange Lentivirus (puro)		One vial: 0.6 mL	SX5	4809

\*Pre-labeled Nuclight cell lines are also available for purchase. Please visit [shop.incucyte.com](http://shop.incucyte.com) for more information

## Spheroid

Product	Color	Quantity	Compatibility	Cat. No.
Incucyte® Cytolight Green Lentivirus (puro)	■	One vial: 0.6 mL	SX1, S3, SX5	4481
Incucyte® Cytolight Red Lentivirus (puro)	■	One vial: 0.6 mL	SX1, S3, SX5 (Green/Red Optical Module)	4482

## Immune Cell Killing, Activation & Proliferation

Product	Color	Quantity	Compatibility	Cat. No.
Incucyte® Annexin V Green Dye	■	One vial: 100 tests	SX1, S3, SX5	4642
Incucyte® Annexin V Red Dye	■	One vial: 100 tests	SX1, S3, SX5 (Green/Red Optical Module)	4641
Incucyte® Annexin V NIR Dye	■	One vial: 100 tests	S3 for Neuroscience, SX5	4768
Incucyte® Annexin V Orange Dye	■	One vial: 100 tests	S3 for Neuroscience, SX5	4759
Incucyte® Caspase-3/7 Green Dye	■	One vial: 20 µL	SX1, S3, SX5	4440
Incucyte® Caspase-3/7 Red Dye	■	One vial: 20 µL	SX1, S3, SX5 (Green/Red Optical Module)	4704
Incucyte® Cytolight Rapid Green Dye	■	One vial: 100-1,000 tests	SX1, S3, SX5	4705
Incucyte® Cytolight Rapid Red Dye	■	Five vials: 100-1,000 tests each	SX1, S3, SX5 (Green/Red Optical Module)	4706

## Scratch Wound Cell Migration & Invasion

Product	Quantity	Compatibility	Cat. No.
Incucyte® Imagelock 96-well Plates	50 plates	SX1, S3, SX5	4379
	10 plates		4806
Incucyte® 96-well Woundmaker Tool	One wounder	SX1, S3, SX5	4563
Incucyte® Cell Migration Bundle	One kit	SX1, S3, SX5	4493
Incucyte® Cell Migration   Invasion Bundle	One kit	SX1, S3, SX5	4474
Incucyte® Cell Invasion Accessories	One kit	SX1, S3, SX5	4444

## Chemotaxis

Product	Quantity	Compatibility	Cat. No.
Incucyte® Clearview 96-well Plates	1 plate	SX1, S3, SX5	4582
	10 plates		4648
	50 plates		4599
Incucyte® Clearview 96-well Reservoir Plates	10 plates	SX1, S3, SX5	4600
	100 plates		4601

## Immunocytochemistry

Product	Color	Quantity	Compatibility	Cat. No.
Incucyte® Mouse IgG1 Fabfluor-488 Antibody Labeling Dye	■	One vial: 50 µg	SX1, S3, SX5	4745
Incucyte® Mouse IgG2a Fabfluor-488 Antibody Labeling Dye	■	One vial: 50 µg	SX1, S3, SX5	4743
Incucyte® Mouse IgG2b Fabfluor-488 Antibody Labeling Dye	■	One vial: 50 µg	SX1, S3, SX5	4744

**Antibody Internalization**

Product	Color	Quantity	Compatibility	Cat. No.
Incucyte® Human Fabfluor-pH Red Antibody Labeling Dye	■	One vial: 50 µg	SX1, S3, SX5 (Green/Red Optical Module)	4722
Incucyte® Human Fabfluor-pH Orange Antibody Labeling Dye	■	One vial: 50 µg	S3 for Neuroscience, SX5	4812
Incucyte® Mouse IgG1 Fabfluor-pH Red Antibody Labeling Dye	■	One vial: 50 µg	SX1, S3, SX5 (Green/Red Optical Module)	4723
Incucyte® Mouse IgG2a Fabfluor-pH Red Antibody Labeling Dye	■	One vial: 50 µg	SX1, S3, SX5 (Green/Red Optical Module)	4750
Incucyte® Mouse IgG2b Fabfluor-pH Red Antibody Labeling Dye	■	One vial: 50 µg	SX1, S3, SX5 (Green/Red Optical Module)	4751
Incucyte® Rat Fabfluor-pH Red Antibody Labeling Dye	■	One vial: 50 µg	SX1, S3, SX5 (Green/Red Optical Module)	4737

**NEW! Metabolism**

Product	Color	Quantity	Compatibility	Cat. No.
Incucyte® CytoATP Lentivirus Reagent Kit ■ Incucyte® CytoATP Lentivirus ■ Incucyte® CytoATP Non-binding Control Lentivirus (puro)	■	One kit	SX5 (Metabolism)	4772
Incucyte® MMP Orange Reagent Kit ■ MMP Orange Reagent ■ FCCP ■ Oligomycin A	■	One kit	S3 for Neuroscience, SX5	4775
Incucyte® Caspase-3/7 Orange Dye for Metabolism	■	One vial: 20 µL	S3 for Neuroscience, SX5 (Metabolism)	4776

**Neuronal Activity**

Product	Color	Quantity	Compatibility	Cat. No.
Incucyte® Neuroburst Orange Lentivirus	■	One vial: 0.2 mL	S3 for Neuroscience, SX5	4736
Incucyte® Neuroactive Orange Kit	■	One kit	S3 for Neuroscience, SX5	4761

**Neurite Outgrowth: Monoculture & Co-culture**

Product	Color	Quantity	Compatibility	Cat. No.
Incucyte® Neurolight Red Lentivirus (Synapsin)	■	Two vials: 0.45 mL each	SX1, S3, SX5 (Green/Red Optical Module)	4807
Incucyte® Neurolight Orange Lentivirus (Synapsin)	■	Two vials: 0.45 mL each	S3 for Neuroscience, SX5	4808
Incucyte® Neuroprime Red Kit	■	One kit	SX1, S3, SX5 (Green/Red Optical Module)	4585
Incucyte® Neuroprime Orange Kit	■	One kit	S3 for Neuroscience, SX5	4760
Incucyte® rCortical Neurons	n/a	One vial: 2 x 10 <sup>6</sup> cells	SX1, S3, SX5	4753
Incucyte® rAstrocytes	n/a	One vial: 2 x 10 <sup>6</sup> cells	SX1, S3, SX5	4586

## Phagocytosis

Product	Color	Quantity	Compatibility	Cat. No.
Incucyte® pHrodo® Red Cell Labeling Kit	■	One kit	SX1, S3, SX5 (Green/Red Optical Module)	4649
Incucyte® pHrodo® Orange Cell Labeling Kit	■	One kit	S3 for Neuroscience, SX5	4766
Incucyte® pHrodo® Red E. coli Bipoarticles®	■	One vial: 2 mg	SX1, S3, SX5 (Green/Red Optical Module)	4615
Incucyte® pHrodo® Green E. coli Bipoarticles®	■	One vial: 2 mg	SX1, S3, SX5	4616
Incucyte® pHrodo® Red Zymosan Bipoarticles®	■	One vial: 1 mg	SX1, S3, SX5 (Green/Red Optical Module)	4617
Incucyte® pHrodo® Green Zymosan Bipoarticles®	■	One vial: 1 mg	SX1, S3, SX5	4618
Incucyte® pHrodo® Red S. aureus Bipoarticles®	■	One vial: 2 mg	SX1, S3, SX5 (Green/Red Optical Module)	4619
Incucyte® pHrodo® Green S. aureus Bipoarticles®	■	One vial: 2 mg	SX1, S3, SX5	4620

## Angiogenesis

Product	Color	Quantity	Compatibility	Cat. No.
Incucyte® Primekit	■	One kit	SX1, S3, SX5	4452
Incucyte® Primekit VEGF/Suramin Kit	■	Two vials	SX1, S3, SX5	4437
Incucyte® Primekit bFGF/Suramin Kit	■	Two vials	SX1, S3, SX5	4438
Incucyte® Primekit Assay Media Kit	■	One bottle: 125 mL	SX1, S3, SX5	4541
Incucyte® Cytolight Green HUVEC	■	One vial: 1.7 x 10 <sup>6</sup> cells	SX1, S3, SX5	4453
Incucyte® Cytolight Green Lentivirus	■	One vial: 0.6 mL	SX1, S3, SX5	4513

Ordering reagents is now much easier!

Visit our online store:

[shop.incucyte.com](http://shop.incucyte.com)

## Sales and Service Contacts

For further contacts, visit  
[www.sartorius.com](http://www.sartorius.com)

### Essen BioScience, A Sartorius Company

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

E-Mail: [AskAScientist@sartorius.com](mailto:AskAScientist@sartorius.com)

#### North America

Essen BioScience Inc.  
300 West Morgan Road  
Ann Arbor, Michigan, 48108  
USA  
Telephone +1 734 769 1600  
E-Mail: [orders.US07@sartorius.com](mailto:orders.US07@sartorius.com)

#### Europe

Essen BioScience Ltd.  
Units 2 & 3 The Quadrant  
Newark Close  
Royston Hertfordshire  
SG8 5HL  
United Kingdom  
Telephone +44 1763 227400  
E-Mail:  
[euorders.UK03@sartorius.com](mailto:euorders.UK03@sartorius.com)

#### APAC

Essen BioScience K.K.  
4th Floor Daiwa Shinagawa North  
Bldg.  
1-8-11 Kita-Shinagawa  
Shinagawa-ku, Tokyo  
140-0001  
Japan  
Telephone: +81 3 6478 5202  
E-Mail: [orders.US07@sartorius.com](mailto:orders.US07@sartorius.com)

Specifications subject to change without notice.

© 2020. All rights reserved. Incucyte, Essen BioScience, and all names of Essen BioScience products are registered trademarks and the property of Essen BioScience unless otherwise specified. Essen BioScience is a Sartorius Company. Publication No.: 8000-0741-A00

Status: 08 | 2020