










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	4633
Incucyte® Cytotox Red Dye		Five vials: 5 µL each	SX1, S3, SX5 (Green/Red Optical Module)	4632
Incucyte® Cytotox NIR Dye		One vial: 100 µL	S3 for Neuroscience, SX5	4846

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 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: 15 µg	SX1, S3, SX5	4705
Incucyte® Cytolight Rapid Red Dye	■	Five vials: 50 µg	SX1, S3, SX5 (Green/Red Optical Module)	4706
Incucyte® Cytolight Rapid Orange Dye	■	One vial: 1 mg	S3 for Neuroscience, SX5	4839

Scratch Wound Cell Migration & Invasion

Product	Quantity	Compatibility	Cat. No.
Incucyte® Imagelock 96-well Plates	50 plates	SX1, S3, SX5	BA-04857
	10 plates		BA-04856
Incucyte® 96-well Woundmaker Tool	One wounder	SX1, S3, SX5	4563
Incucyte® Cell Migration Kit	One kit	SX1, S3, SX5	BA-04858
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 Kit	■	One kit	SX1, S3, SX5	4745
■ Incucyte® Mouse IgG1 Fabfluor-488 Dye				
■ Incucyte® Opti-Green Background Suppressor				
Incucyte® Mouse IgG2a Fabfluor-488 Antibody Labeling Kit	■	One kit	SX1, S3, SX5	4743
■ Incucyte® Mouse IgG12a Fabfluor-488 Dye				
■ Incucyte® Opti-Green Background Suppressor				
Incucyte® Mouse IgG2b Fabfluor-488 Antibody Labeling Kit	■	One kit	SX1, S3, SX5	4744
■ Incucyte® Mouse IgG12b Fabfluor-488 Dye				
■ Incucyte® Opti-Green Background Suppressor				
Incucyte® Mouse IgG1 Fabfluor-594 Antibody Labeling Dye	■	One vial: 50 µg	SX1, S3, SX5 (Green/Red Optical Module)	4844
Incucyte® Mouse IgG2a Fabfluor-594 Antibody Labeling Dye	■	One vial: 50 µg	SX1, S3, SX5 (Green/Red Optical Module)	BA-04863



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





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 ⁶ cells	SX1, S3, SX5	4753
Incucyte® rAstrocytes	n/a	One vial: 2 x 10 ⁶ cells	SX1, S3, SX5	4586

NEW! Kinase Activity

Product	Color	Quantity	Compatibility	Cat. No.
Incucyte® Kinase Akt Green/Red Lentivirus (puro)		One vial: 0.2 mL	SX1, S3, SX5 (Green/Red Optical Module)	BA-04868

Phagocytosis

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

North America

Sartorius Corporation
565 Johnson Avenue
Bohemia, NY 11716
USA
Phone +1 734 769 1600

Europe

Sartorius UK Ltd.
Longmead Business Centre
Blenheim Road
Epsom
Surrey, KT19 9QQ
United Kingdom
Phone +44 1763 227400

Asia Pacific

Sartorius Japan K.K.
4th Floor, Daiwa Shinagawa North Bldg.
1-8-11, Kita-Shinagawa 1-chome
Shinagawa-Ku
Tokyo 140-0001
Japan
Phone +81 3 6478 5202

Specifications subject to change without notice.

© 2021. All rights reserved. Incucyte and all names of Sartorius products are registered trademarks and the property of Sartorius AG and/or one of its affiliated companies.
Incucyte-Reagents-Consumables-Flyer-en-L-8000-0741-C00-Sartorius Status: 11 | 2021