










## Incucyte® Live-Cell Analysis 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, CX3	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	SX5, CX3	4768
Incucyte® Annexin V Orange Dye		One vial: 100 tests	SX5, CX3	4759
Incucyte® Caspase-3/7 Green Dye		One vial: 20 µL	SX1, S3, SX5, CX3	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, CX3	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	SX5, CX3	4846

### Proliferation

Product	Color	Quantity	Compatibility	Cat. No.
Incucyte® Nuclight Red Lentivirus 2.0 (puro)		One vial: 0.6 mL	SX1, S3, SX5 (Green/Red Optical Module)	BA-04887
Incucyte® Nuclight Green Lentivirus 2.0 (puro)		One vial: 0.6 mL	SX1, S3, SX5, CX3	BA-04888
Incucyte® Nuclight Red Lentivirus 2.0 (bleo)		One vial: 0.6 mL	SX1, S3, SX5 (Green/Red Optical Module)	BA-04889
Incucyte® Nuclight Green Lentivirus 2.0 (bleo)		One vial: 0.6 mL	SX1, S3, SX5, CX3	BA-04890
Incucyte® Nuclight Orange Lentivirus 2.0 (puro)		One vial: 0.6 mL	SX5, CX3	BA-04891
Incucyte® Nuclight NIR Lentivirus (puro)		One vial: 0.2 mL	SX5, CX3	4805
Incucyte® Nuclight Rapid Red Dye 2.0		One vial: 50 µg	SX1, S3, SX5 (Green/Red Optical Module)	BA-04903
Incucyte® Nuclight Rapid Green Dye		One vial: 50 µg	SX1, S3, SX5, CX3	BA-04902
Incucyte® Nuclight Rapid Orange Dye		One vial: 50 µg	SX5, CX3	BA-04904
Incucyte® Nuclight Rapid NIR Dye		One vial: 50 µL	SX5, CX3	4804










### Cell Cycle

Product	Color	Quantity	Compatibility	Cat. No.
Incucyte® Cell Cycle Green/Red Lentivirus 2.0		One vial: 0.2 mL	SX1, S3, SX5 (Green/Red Optical Module)	BA-04892
Incucyte® Cell Cycle Green/Orange Lentivirus 2.0		One vial: 0.2 mL	SX5, CX3	BA-04889

Ordering reagents is now much easier!  
Visit our online store: [www.sartorius.com/incucyte-shop](http://www.sartorius.com/incucyte-shop)



## Immune Cell Killing, Activation & Proliferation

Product	Color	Quantity	Compatibility	Cat. No.
Incucyte® Annexin V Green Dye		One vial: 100 tests	SX1, S3, SX5, CX3	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	SX5, CX3	4768
Incucyte® Annexin V Orange Dye		One vial: 100 tests	SX5, CX3	4759
Incucyte® Caspase-3/7 Green Dye		One vial: 20 µL	SX1, S3, SX5, CX3	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, CX3	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	SX5, CX3	4839







## Scratch Wound Cell Migration & Invasion

Product	Quantity	Compatibility	Cat. No.
Incucyte® Imagelock 96-well Plates	50 plates	SX1, S3, SX5, CX3	BA-04857
	10 plates		BA-04856
Incucyte® 96-well Woundmaker Tool	One wounder	SX1, S3, SX5, CX3	4563
Incucyte® Cell Migration Kit	One kit	SX1, S3, SX5, CX3	BA-04858

## Chemotaxis

Product	Quantity	Compatibility	Cat. No.
Incucyte® Clearview 96-well Plates	1 plate	SX1, S3, SX5, CX3	4582
	10 plates		4648
	50 plates		4599
Incucyte® Clearview 96-well Reservoir Plates	10 plates	SX1, S3, SX5, CX3	4600
	100 plates		4601

## Immunocytochemistry

Product	Color	Quantity	Compatibility	Cat. No.
Incucyte® Mouse IgG1 Fabfluor-488 Antibody Labeling Kit		One kit	SX1, S3, SX5, CX3	4745
<ul style="list-style-type: none"> <li>■ Incucyte® Mouse IgG1 Fabfluor-488 Dye</li> <li>■ Incucyte® Opti-Green Background Suppressor</li> </ul>				
Incucyte® Mouse IgG2a Fabfluor-488 Antibody Labeling Kit		One kit	SX1, S3, SX5, CX3	4743
<ul style="list-style-type: none"> <li>■ Incucyte® Mouse IgG12a Fabfluor-488 Dye</li> <li>■ Incucyte® Opti-Green Background Suppressor</li> </ul>				
Incucyte® Mouse IgG2b Fabfluor-488 Antibody Labeling Kit		One kit	SX1, S3, SX5, CX3	4744
<ul style="list-style-type: none"> <li>■ Incucyte® Mouse IgG12b Fabfluor-488 Dye</li> <li>■ Incucyte® Opti-Green Background Suppressor</li> </ul>				
Incucyte® Mouse IgG1 Fabfluor-555 Antibody Labeling Kit		One kit	SX5, CX3	BA-04873
<ul style="list-style-type: none"> <li>■ Incucyte® Mouse IgG1 Fabfluor-555 Dye</li> <li>■ Incucyte® Opti-Orange Background Suppressor</li> </ul>				
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

## Neuronal Activity

Product	Color	Quantity	Compatibility	Cat. No.
Incucyte® Neuroburst Orange Lentivirus	■	One vial: 2 mL	SX5, CX3	4736
Incucyte® Neuroactive Orange Kit	■	One kit	SX5, CX3	4761

## Neurite Outgrowth: Monoculture & Co-culture

Product	Color	Quantity	Compatibility	Cat. No.
Incucyte® rCortical Neurons	n/a	One vial: 2 x 10 <sup>6</sup> cells	SX1, S3, SX5, CX3	4753
Incucyte® rAstrocytes	n/a	One vial: 2 x 10 <sup>6</sup> cells	SX1, S3, SX5, CX3	4586

## 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	SX5, CX3	4812
Incucyte® Mouse IgG1 Fabfluor-pH Red Antibody Labeling Dye	■	One vial: 50 µg	SX1, S3, SX5 (Green/Red Optical Module)	4723
Incucyte® Mouse IgG1 Fabfluor-pH Orange Antibody Labeling Dye	■	One vial: 50 µg	SX5, CX3	BA-04885
Incucyte® Mouse IgG2a Fabfluor-pH Red Antibody Labeling Dye	■	One vial: 50 µg	SX1, S3, SX5 (Green/Red Optical Module)	4750
Incucyte® Mouse IgG2a Fabfluor-pH Orange Antibody Labeling Dye	■	One vial: 50 µg	SX5, CX3	BA-04886
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 <ul style="list-style-type: none"> <li>■ Incucyte® CytoATP Lentivirus</li> <li>■ Incucyte® CytoATP Non-binding Control Lentivirus (puro)</li> </ul>	■	One kit	SX5 (Metabolism)	4772
Incucyte® MMP Orange Reagent Kit <ul style="list-style-type: none"> <li>■ MMP Orange Reagent</li> <li>■ FCCP</li> <li>■ Oligomycin A</li> </ul>	■	One kit	SX5	4775
Incucyte® Caspase-3/7 Orange Dye for Metabolism	■	One vial: 20 µL	SX5 (Metabolism)	4776

## 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	SX5, CX3	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, CX3	4616

## EV Uptake Assay

Product		Quantity	Compatibility	Cat. No.
Incucyte® ExoFluor Green EV Labeling Kit		One kit	SX1, S3, SX5, CX3	BA-04877
Incucyte® ExoFluor Red EV Labeling Kit		One kit	SX1, S3, SX5 (Green/Red Optical Module)	BA-04900
Incucyte® ExoFluor NIR EV Labeling Kit		One kit	SX5, CX3	BA-04901

### 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

 For further information, visit  
[sartorius.com](https://www.sartorius.com)