

IncuCyte® Live-Cell Analysis System

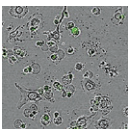
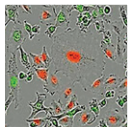
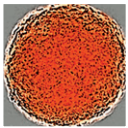
Reagents, Cell Lines and Consumables

Use IncuCyte® Reagents, Cell Lines, and Consumables with your live-cell analysis system for a turnkey solution to a broad range of cell health, movement and other applications.

Improve efficiency while maintaining cell health with IncuCyte® Reagents that are non-perturbing to cell proliferation and morphology.

Cell Health Assays

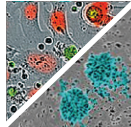
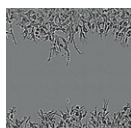
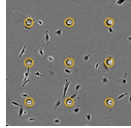
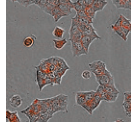
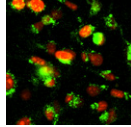
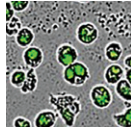
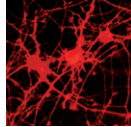
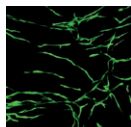
Please see other side for additional product listings.

Apoptosis and Cytotoxicity	Product	Quantity	Cat. No.	
 <p>Visualize and measure apoptotic cell death and cytotoxicity using the IncuCyte® range of no-wash, mix and read reagents.</p>	IncuCyte® Caspase-3/7 Green Apoptosis Reagent	One vial: 20 µL	4440	
	IncuCyte® Caspase-3/7 Red Apoptosis Reagent	One vial: 20 µL	4704	
	IncuCyte® Annexin V Red Reagent	One vial: 100 tests	4641	
	IncuCyte® Annexin V Green Reagent	One vial: 100 tests	4642	
	IncuCyte® Cytotox Red Reagent	Five vials: 5 µL	4632	
	IncuCyte® Cytotox Green Reagent	Five vials: 5 µL	4633	
Proliferation and Cell Cycle	IncuCyte® NuLight Live-Cell Labeling Reagents	Quantity	Cat. No.	
 <p>Count viable cells over hours, days, or weeks with IncuCyte® NuLight live cell labeling reagents and pre-labeled cell lines. Gain additional insight by multiplexing with apoptosis and cytotoxicity readouts or exploring co-cultures.</p>	NuLight Green Lentivirus (EF-1 Alpha Promoter, Puromycin selection)	One vial: 0.2 mL	4624	
	NuLight Red Lentivirus (EF-1 Alpha Promoter, Puromycin selection)	One vial: 0.2 mL	4625	
	NuLight Green Lentivirus (EF-1 Alpha Promoter, Bleomycin selection)	One vial: 0.2 mL	4626	
	NuLight Red Lentivirus (EF-1 Alpha Promoter, Bleomycin selection)	One vial: 0.2 mL	4627	
	NuLight Green Lentivirus (EF-1 Alpha Promoter, Puromycin selection)	One vial: 0.6 mL	4475	
	NuLight Red Lentivirus (EF-1 Alpha Promoter, Puromycin selection)	One vial: 0.6 mL	4476	
	NuLight Green Lentivirus (EF-1 Alpha Promoter, Bleomycin selection)	One vial: 0.6 mL	4477	
	NuLight Red Lentivirus (EF-1 Alpha Promoter, Bleomycin selection)	One vial: 0.6 mL	4478	
	NuLight Rapid Red Reagent	One vial: 50 µL	4717	
	<p>Further characterize growth or probe drug mechanisms of action in real-time with the IncuCyte® Cell Cycle live-cell labeling reagent.</p>	IncuCyte® NuLight Cell Lines	Quantity	Cat. No.
A549 NuLight Green Cells		One vial: 1x10 ⁶ cells	4492	
A549 NuLight Red Cells		One vial: 1x10 ⁶ cells	4491	
HeLa NuLight Green Cells		One vial: 1x10 ⁶ cells	4490	
HeLa NuLight Red Cells		One vial: 1x10 ⁶ cells	4489	
HT-1080 NuLight Green Cells		One vial: 1x10 ⁶ cells	4486	
HT-1080 NuLight Red Cells		One vial: 1x10 ⁶ cells	4485	
Jurkat NuLight Red Cells		One vial: 1.5x10 ⁶ cells	4613	
MCF7 NuLight Green Cells		One vial: 1x10 ⁶ cells	4528	
MCF7 NuLight Red Cells		One vial: 1x10 ⁶ cells	4524	
Neuro-2a NuLight Green Cells		One vial: 2x10 ⁶ cells	4511	
Neuro-2a NuLight Red Cells		One vial: 2x10 ⁶ cells	4512	
IncuCyte® Cell Cycle Live-Cell Labeling Reagent		IncuCyte® Cell Cycle Live-Cell Labeling Reagent	Quantity	Cat. No.
 <p>IncuCyte® CytoLight live-cell labeling reagents enable real-time measurement of spheroid growth and shrinkage.</p>		NEW IncuCyte® Cell Cycle Red/Green Lentivirus (EF-1 Alpha Promoter, Puromycin selection)	One vial: 0.2 mL	4779
Spheroids	IncuCyte® CytoLight Live-Cell Labeling Reagents	Quantity	Cat. No.	
	CytoLight Green Lentivirus (EF-1 Alpha Promoter, Puromycin selection)	One vial: 0.6mL	4481	
	CytoLight Red Lentivirus (EF-1 Alpha Promoter, Puromycin selection)	One vial: 0.6mL	4482	

Ordering reagents is now much easier!
 Visit our online store: shop.incucyte.com

Learn more at essenbioscience.com/consumables
 or contact us at sales@essenbio.com

Other Live-Cell Assays

		Product	Quantity	Cat. No.
 <p>Immune Cell Killing, Clustering & Proliferation</p> <p>Directly measure tumor cell killing and proliferation in real time.</p>		IncuCyte® Caspase-3/7 Green Apoptosis Reagent	One vial: 20 µL	4440
		IncuCyte® Caspase-3/7 Red Apoptosis Reagent	One vial: 20 µL	4704
		IncuCyte® Annexin V Red Reagent	One vial: 100 tests	4641
		IncuCyte® Annexin V Green Reagent	One vial: 100 tests	4642
		IncuCyte® Cytolight Rapid Green Reagent	One vial: 100–1,000 tests	4705
		IncuCyte® Cytolight Rapid Red Reagent	Five vials: 100–1,000 tests	4706
 <p>Cell Migration and Invasion</p> <p>ImageLock™ plates enable precise, scan-to-scan repeat imaging of the same field of view. CoolBoxes and accessories enable fine temperature control of biomatrices.</p>		ImageLock™ 96-well Plates	50 plates	4379
		IncuCyte® Cell Invasion Accessories	One kit	4444
 <p>Chemotaxis</p> <p>A novel, integrated solution for real-time visualization and automated analysis of chemotactic cell migration in a 96-well format.</p>		IncuCyte® ClearView 96-Well Cell Migration Plate	1 plate	4582
		IncuCyte® ClearView 96-Well Cell Migration Plates	10 plates	4648
		IncuCyte® ClearView 96-Well Cell Migration Plates	50 plates	4599
		IncuCyte® ClearView Reservoir Plates	10 plates	4600
		IncuCyte® ClearView Reservoir Plates	100 plates	4601
 <p>Antibody Internalization</p> <p>Perform rapid, kinetic, high throughput assays suitable for antibody screening or therapeutic profiling.</p>		IncuCyte® Human FabFluor-pH Red Antibody Labeling Reagent	One vial: 50 µg	4722
		IncuCyte® Mouse IgG1 FabFluor-pH Red Antibody Labeling Reagent	One vial: 50 µg	4723
		IncuCyte® Rat FabFluor-pH Red Antibody Labeling Reagent	One vial: 50 µg	4737
		IncuCyte® Mouse IgG2a FabFluor-pH Red Antibody Labeling Reagent	One vial: 50 µg	4750
		IncuCyte® Mouse IgG2b FabFluor-pH Red Antibody Labeling Reagent	One vial: 50 µg	4751
 <p>Immunocytochemistry</p> <p>Reveal the dynamics of surface protein expression with a new approach to immunocytochemistry.</p>		IncuCyte® Mouse IgG2a FabFluor-488 Antibody Labeling Reagent	One vial: 50 µg	4743
		IncuCyte® Mouse IgG2b FabFluor-488 Antibody Labeling Reagent	One vial: 50 µg	4744
		IncuCyte® Mouse IgG1 FabFluor-488 Antibody Labeling Reagent	One vial: 50 µg	4745
 <p>Phagocytosis</p> <p>Visualize and quantify the phagocytosis of cells or Bioparticles® in real time with pH sensitive pHrodo® indicators.</p>		IncuCyte® pHrodo® Red Cell Labeling Kit	One Kit	4649
		IncuCyte® pHrodo® Red E. coli Bioparticles®	One vial: 2 mg	4615
		IncuCyte® pHrodo® Green E. coli Bioparticles®	One vial: 2 mg	4616
		IncuCyte® pHrodo® Red Zymosan Bioparticles®	One vial: 1 mg	4617
		IncuCyte® pHrodo® Green Zymosan Bioparticles®	One vial: 1 mg	4618
		IncuCyte® pHrodo® Red S. aureus Bioparticles®	One vial: 2 mg	4619
		IncuCyte® pHrodo® Green S. aureus Bioparticles®	One vial: 2 mg	4620
 <p>Neurite Dynamics: Monoculture & Coculture</p> <p>Neuron specific live-cell labeling reagents, cell kits and pre-labeled cells. Measure neurite outgrowth and retraction in monoculture or coculture.</p>		IncuCyte® NeuroLight Red Lentivirus - Synapsin Promotor	One vial: 0.45mL	4584
		IncuCyte® NeuroPrime Cell Kit	One kit	4585
		Neuro-2a NuLight Green Cells	One vial: 2x10 ⁶ cells	4511
		Neuro-2a NuLight Red Cells	One vial: 2x10 ⁶ cells	4512
 <p>Angiogenesis</p> <p>Measure vascular tube formation using our range of human coculture angiogenesis assay kits and components.</p>		PrimeKit (GFP-labeled HUVEC and NHDF coculture model)		
		IncuCyte® Angiogenesis 96-well PrimeKit	One kit	4452
		IncuCyte® Angiogenesis PrimeKit VEGF/Suramin Supplement Kit	Two vials	4437
		IncuCyte® Angiogenesis PrimeKit bFGF/Suramin Supplement Kit	Two vials	4438
		IncuCyte® Angiogenesis PrimeKit Optimized Assay Medium	One bottle: 125 mL	4541
		HUVEC Cytolight Green Cells	One vial: 1.7x10 ⁵ cells	4453
Cytolight Green Lentivirus (CMV Promoter, no selection)	One vial: 0.6mL	4513		