

# IncuCyte® Live-Cell Analysis System

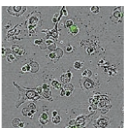
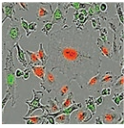
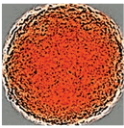
## Reagents, Cell Lines and Consumables

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

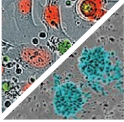
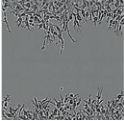
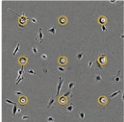
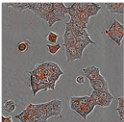
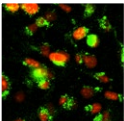
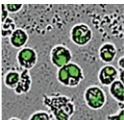
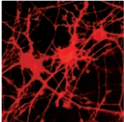
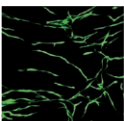
Please see other side for additional product listings.

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 – Cell Count		IncuCyte® NuLight Live-Cell Labeling Reagents	Quantity	Cat. No.
 <p>IncuCyte® NuLight live-cell labeling reagents and pre-labeled cell lines enable real-time counting of viable cells over hours, days and weeks. Calculate growth rates and doubling times, explore cocultures and multiplex with apoptosis and cytotoxicity readouts.</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	
	IncuCyte® NuLight Cell Lines		Quantity	Cat. No.
	A549 NuLight Green Cells	One vial: 1x10 <sup>6</sup> cells	4492	
	A549 NuLight Red Cells	One vial: 1x10 <sup>6</sup> cells	4491	
	HeLa NuLight Green Cells	One vial: 1x10 <sup>6</sup> cells	4490	
	HeLa NuLight Red Cells	One vial: 1x10 <sup>6</sup> cells	4489	
HT-1080 NuLight Green Cells	One vial: 1x10 <sup>6</sup> cells	4486		
HT-1080 NuLight Red Cells	One vial: 1x10 <sup>6</sup> cells	4485		
Jurkat NuLight Red Cells	One vial: 1.5x10 <sup>6</sup> cells	4613		
MCF7 NuLight Green Cells	One vial: 1x10 <sup>6</sup> cells	4528		
MCF7 NuLight Red Cells	One vial: 1x10 <sup>6</sup> cells	4524		
Neuro-2a NuLight Green Cells	One vial: 2x10 <sup>6</sup> cells	4511		
Neuro-2a NuLight Red Cells	One vial: 2x10 <sup>6</sup> cells	4512		
Spheroids		IncuCyte® Cytolight Live-Cell Labeling Reagents	Quantity	Cat. No.
 <p>IncuCyte® Cytolight live-cell labeling reagents enable real-time measurement of spheroid growth and shrinkage.</p>	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](http://shop.incucyte.com)

Learn more at [essenbioscience.com/consumables](http://essenbioscience.com/consumables)  
or contact us at [sales@essenbio.com](mailto:sales@essenbio.com)

## Other Live-Cell Assays

		<b>Product</b>	<b>Quantity</b>	<b>Cat. No.</b>
 <p>Directly measure tumor cell killing and proliferation in real time.</p>	<b>Immune Cell Killing, Clustering &amp; Proliferation</b>			
	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>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>	<b>Cell Migration and Invasion</b>			
	IncuCyte® ImageLock 96-well Plates	50 plates	4379	
IncuCyte® Cell Invasion Accessories	One kit	4444		
 <p>A novel, integrated solution for real-time visualization and automated analysis of chemotactic cell migration in a 96-well format.</p>	<b>Chemotaxis</b>			
	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>Perform rapid, kinetic, high throughput assays suitable for antibody screening or therapeutic profiling.</p>	<b>Antibody Internalization</b>			
	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>Reveal the dynamics of surface protein expression with a new approach to immunocytochemistry.</p>	<b>Immunocytochemistry</b>			
	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>Visualize and quantify the phagocytosis of cells or Bioparticles® in real time with pH sensitive pHrodo® indicators.</p>	<b>Phagocytosis</b>			
	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>Neuron specific live-cell labeling reagents, cell kits and pre-labeled cells. Measure neurite outgrowth and retraction in monoculture or coculture.</p>	<b>Neurite Dynamics – Monoculture &amp; Coculture</b>			
	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 <sup>6</sup> cells	4511	
	Neuro-2a NuLight Red Cells	One vial: 2x10 <sup>6</sup> cells	4512	
 <p>Measure vascular tube formation using our range of human coculture angiogenesis assay kits and components.</p>	<b>Angiogenesis</b>			
	<b>PrimeKit (GFP-labeled HUVEC and NHDF coculture model)</b>			
	IncuCyte® Angiogenesis 96-well PrimeKit	One kit	4452	
	IncuCyte® Angiogenesis PrimeKit VEGF/Suramin Supplement Kit	Two vials	4437	
	IncuCyte® Angiogenesis PrimeKit Optimized Assay Medium	One bottle: 125 mL	4541	
	HUVEC CytoLight Green Cells	One vial: 1.7x10 <sup>5</sup> cells	4453	
CytoLight Green Lentivirus (CMV Promoter, no selection)	One vial: 0.6mL	4513		