

IncuCyte® Live-Cell Analysis System

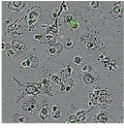
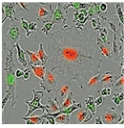
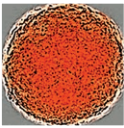
Reagents, Cell Lines and Consumables

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

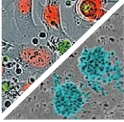
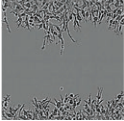
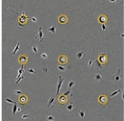
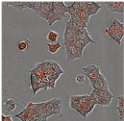
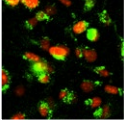
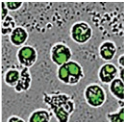
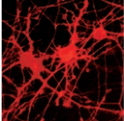
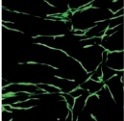
Please see other side for additional product listings.

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>		IncuCyte® NuLight Green BacMam 3.0 Reagent for Nuclear Labeling	One vial: 1 mL	4622
		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
		NEW NuLight Rapid Red Reagent	One vial: 50 µL	4717
	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	
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

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

Other Live-Cell Assays

 <p>Directly measure tumor cell killing and proliferation in real time.</p>	Immune Cell Killing, Clustering & Proliferation	Product	Quantity	Cat. No.
	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>	Cell Migration and Invasion	Product	Quantity	Cat. No.
	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>	Chemotaxis	Product	Quantity	Cat. No.
	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>	NEW Antibody Internalization	Product	Quantity	Cat. No.
	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>	NEW Immunocytochemistry	Product	Quantity	Cat. No.
	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>	Phagocytosis	Product	Quantity	Cat. No.
	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>	Neurite Dynamics – Monoculture & Coculture	Product	Quantity	Cat. No.
	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>Measure vascular tube formation using our range of human coculture angiogenesis assay kits and components.</p>	Angiogenesis	PrimeKit (GFP-labeled HUVEC and NHDF coculture model)	Quantity	Cat. No.
	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	