

Incucyte[®]
Reagents, Consumables
and Software

Simplifying Progress



Incucyte® Reagents, Consumables and Software

Sartorius offers a range of reagents, consumables and software specifically designed around your application needs to enable long-term, kinetic live-cell imaging and analysis. Obtain specific and robust measurements

of cell health, movement and function with optimized, turnkey solutions allowing you to connect phenotype and function with pathological processes. Getting biologically relevant information has never been easier!

Find Your Solution to Meet Your Application Needs:

Cell Health

Cell Function

Cell Movement & Morphology

Assays for 3D Models

Key Advantages

- Derive meaningful data with sensitive, non-perturbing reagents for long term, live-cell studies.
- Unlock your productivity with lab-tested protocols and purpose-built, intuitive software.
- Generate data rich information within every sample with fluorescent reagent combinations.
- Support visualization and automation of cell movement studies with Incucyte[®] Consumables.

Incucyte® Reagents At-a-Glance

Application	Reagent Consumabl		Software Module		Compatible Instrument		
Cell Health				SX5	S3	S3 for Neuro	SX1
Proliferation:							
Label-free, Confluence					•	•	•
Label-free, Cell Counting						•	
Fluorescent Labeling, Cell Counting							•
Viability							•
Cell Cycle							
Apoptosis							•
Cytotoxicity							
Mitochondrial Membrane Potential						•	
ATP Metabolism							
Cell Function				SX5	S3	S3 for Neuro	SX1
mmune Cell Killing	•				-	•	
Antibody Internalization						•	•
Live-Cell Immunocytochemistry							
Phagocytosis						•	
NETosis							
Angiogenesis							
Neuronal Activity							
Cell Movement and Morphology				SX5	S3	S3 for Neuro	SX1
Chemotaxis Migration and Invasion		•	•				•
Scratch Wound Migration and Invasion			•			•	•
Immune Cell Activation and Proliferation							
Neurite Outgrowth			•			•	•
Assays for 3D Models				SX5	S3	S3 for Neuro	SX1
Spheroid Growth (Single and Multi-Spheroid)			•	•	•		
Spheroid Invasion (Single Spheroid)			•	•	•	•	•
Spheroid Immune Cell Killing					•	•	

Proliferation, Cell Counting, Viability and Cell Cycle

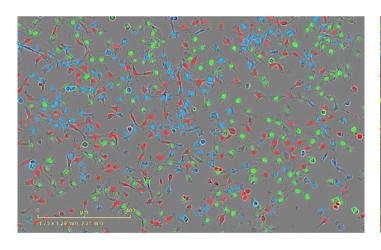
Incucyte's live-cell labeling reagents and purpose-built software enables long-term (>48 hr) analysis of cell proliferation, viability and cell cycle state.

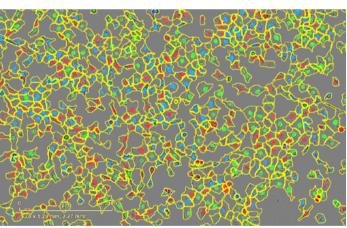
- Minimize experimental artifacts using non-perturbing, live-cell reagents—preservative free, sterile, or concentrated formulations to protect cell health.
- Quantify cell proliferation kinetically using label or labelfree strategies in adherent or non-adherent cell cultures.
- Investigate a range of cell models (mono-, co- or triculture) to answer relevant scientific questions.
- Maximize each sample with multiplexed readouts of proliferation, cell cycle or cell health that are easily validated by eye.

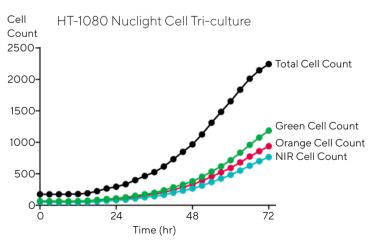
Application Spotlight: Incucyte® Live-Cell Proliferation Assays

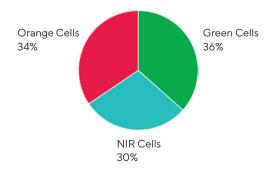
Perform long term, kinetic measurements of proliferation, with or without labels. Incucyte® Nuclight Reagents homogenously label a variety of cell types for non-

perturbing, continuous analysis of the same population of cells. The Incucyte® Cell-by-Cell Analysis Software Module enables label-free cell identification and counting.





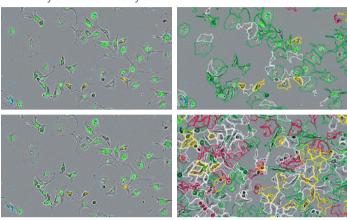


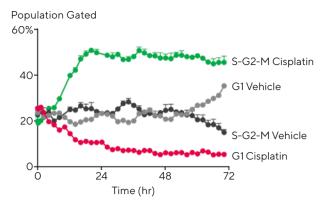


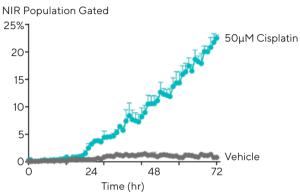
HT-1080 fibrosarcoma cells stably expressing Nuclight Green, Orange, or NIR were monitored for 72 hours. Representative images taken at 48 hours, with and without the label-free Cell-by-Cell Analysis mask, automatically identify the entire population of cells and quantify percentages of green, orange, or NIR expressing cells.

Application Spotlight: Incucyte® Cell Cycle Assay

Continuously quantify treatment effects on cell cycle progression of the same population of cells over multiple cell divisions using Incucyte® Cell Cycle Lentivirus Reagents. Gain deeper insight by multiplexing with Incucyte® Annexin V NIR Dye on the Incucyte® SX5.







	Product	Description	Cat. No.	Instrument Compatibility			
Software	Perform label-free cell counts and subsequent cell-by-cell classification based on shape, size or fluorescence intensity to quantify dynamic changes in cell subsets within heterogeneous living cell cultures.						
	Incucyte® Cell-by-Cell Analysis Software Module	1 module	9600-0031	SX5, S3, S3 for Neuroscience, SX1			
luclear Dye .abeling Reagents	Cell permeable DNA stains that specifically quantification of cell proliferation and viabili		and is ideally sui	ted for mix-and-read, live-cell			
	Incucyte® Nuclight Rapid Red Dye	One vial: 50 μL	4717	SX5 (configured with green/red optica module), S3, SX1			
	Incucyte® Nuclight Rapid NIR Dye	One vial: 50 μL	4804	SX5, S3 for Neuroscience			
*Nuclear Lentivirus Labeling Reagents	Lentivirus reagents provide homogenous ex function for live-cell quantification of cell pro			rescent protein without altering cell			
	Incucyte® Nuclight Green Lentivirus (puro)	0.2 mL	4624	SX5, S3, SX1			
		0.6 mL	4475	SX5, S3, SX1			
	Incucyte® Nuclight Red Lentivirus (puro)	0.2 mL	4625	SX5 (configured with green/red optica			
		0.6 mL	4476	module), S3, SX1			
	Incucyte® Nuclight Green Lentivirus	0.2 mL	4626	SX5, S3, SX1			
	(bleo)	0.6 mL	4477	SX5, S3, SX1			
	Incucyte® Nuclight Red Lentivirus	0.2 mL	4627	SX5 (configured with green/red opti			
	(bleo)	0.6 mL	4478	module), S3, SX1			
	Incucyte® Nuclight Orange Lentivirus (puro)	0.2 mL	4771	SX5, S3 for Neuroscience			
	Incucyte® Nuclight NIR Lentivirus (puro)	0.2 mL	4805	SX5, S3 for Neuroscience			
Cell Cycle entivirus Reagents	Fluorescent ubiquitination-based cell cycle fluorescent proteins to distinguish between						
	Incucyte® Cell Cycle Green/Red Lentivirus (puro)	0.2 mL	4779	SX5 (configured with green/red optica module), S3, SX1			
	Incucyte [®] Cell Cycle Green/Orange Lentivirus (puro)	0.2 mL	4809	SX5			

^{*}Pre-labeled Nuclight cell lines are also available for purchase. Please visit shop.incucyte.com for more information.

Apoptosis, Cytotoxicity, Mitochondrial Membrane Potential and ATP Metabolism

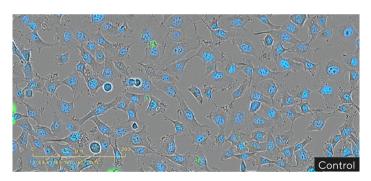
Incucyte's non-perturbing cell health reagents allow for kinetic readouts over multiple days for the evaluation of time-dependent and cell-specific responses to treatments.

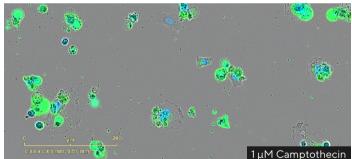
- Preserve cell health using non-perturbing, live-cell reagents to track true, artifact-free biological responses.
- Reduce loss of precious or compromised cells with mix-
- and-read or novel genetically encoded reagents and optimized protocols.
- Perform multi-parametric analysis using cell health reagent combinations within a single well to generate data rich information
- Validate results and confirm changes in cell morphology with HD phase images at every time point.

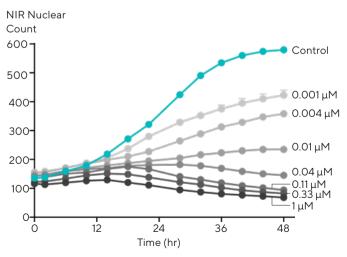
Application Spotlight: Incucyte® Apoptosis Assay

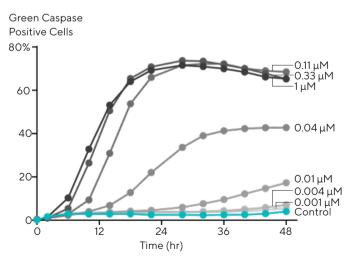
Evaluate cell death with two unique assays for apoptotic pathway analysis using specifically formulated reagents. Utilize Incucyte® Caspase-3/7 Reagents allow direct detection of caspase-3/7 activation or Incucyte® Annexin

V Reagents to measure exposed phosphatidylserine in apoptotic cells. Enhance your insight with multiplexed measurements of nuclear counts of cell death.









HT-1080 fibrosarcoma cells stably expressing Nuclight NIR Lentivirus (pseudo-colored blue) were treated with decreasing concentrations of Camptothecin in the presence of Incucyte Caspase-3/7 Green Dye (pseudo-colored green). Representative images validate kinetic data of both cell viability and apoptotic death.

	Product	Description	Cat. No.	Instrument Compatibility		
Software	Perform label-free cell counts and subseque quantify dynamic changes in cell subsets with	· · · · · · · · · · · · · · · · · · ·		The state of the s		
	Incucyte® Cell-by-Cell Analysis Software Module	1 module	9600-0031	SX5, S3, S3 for Neuroscience, SX1		
Caspase Activity Reagents	Inert, non-fluorescent (DEVD) substrates the caspase-3/7 to release a DNA-binding fluore		ell membrane wh	nere they can be cleaved by activated		
	Incucyte® Caspase-3/7 Green Dye	One vial: 20 μL (100-200 tests)	4440	SX5, S3, SX1		
	Incucyte® Caspase-3/7 Red Dye	One vial: 20 μL (100-200 tests)	4704	SX5 (configured with green/red optica module), S3, SX1		
	Incucyte® Caspase-3/7 Dye for Metabolism	One vial: 20 µL (100-200 tests)	4776	SX5 (configured with SX5 Metabolism Optical Module)		
Apoptosis Plasma Membrane Integrity	Membrane impermeable, highly-selective phextracellular surface of cells undergoing apo		PS) cyanine fluor	rescent dyes label PS exposure on the		
Reagents	Incucyte® Annexin V Green Dye	One vial: 100-200 tests	4642	SX5, S3, SX1		
	Incucyte® Annexin V Red Dye	One vial: 100-200 tests	4641	SX5 (configured with green/red optica module), S3, SX1		
	Incucyte® Annexin V Orange Dye	One vial: 100-200 tests	4759	SX5, S3 for Neuroscience		
	Incucyte® Annexin V NIR Dye	One vial: 100-200 tests	4768	SX5, S3 for Neuroscience		
Cytotoxicity Reagents	Highly sensitive cyanine-based dyes that do not enter living cells. When the cell membrane is compromised, it enters the cell, binds to DNA, and becomes fluorescent. Dye does not enter cells with intact cell membranes.					
	Incucyte® Cytotox Green Dye	Five vials: 5 μL (100 tests each)	4633	SX5, S3, SX1		
	Incucyte® Cytotox Red Dye	Five vials: 5 μL (100 tests each)	4632	SX5 (configured with green/red optica module), S3, SX1		
	Incucyte® Cytotox NIR Dye	One vial: 100 uL (500-100 tests)	4846	SX5, S3 for Neuroscience		
New!	Fluorescent dye which diffuses across the intermembrane space and accumulates in proportion to the mitochondrial					
Mitochondria	membrane potential. Shifts in fluorescent intensity denote mitochondrial membrane potential state. Requires Incucyte® Cell-by-Cell Analysis Software Module for analysis.					
Membrane Potential (MMP)	New! Incucyte® MMP Orange Reagent Kit:	One kit	4775	SX5, S3 for Neuroscience		
	MMP Orange Dye	One vial: 30 µL (200 tests)				
	• FCCP	One vial: 10 µL (16 tests)	_			
	Oligomycin A	One vial: 10 µL (16 tests)	_			
New! ATP Metabolism	Direct, kinetic measurement of ATP to analyz	e changes of cance	r cell metabolisr	m in in advanced cell models.		
Software	Enables analysis of ATP dynamics be capturing cell morphology in each well of a 96- or 384-		es while qualitat	tively monitoring associated changes in		
	New! Incucyte® ATP Analysis Software Module	1 module	9600-0033	SX5 (configured with SX5 Metabolism Optical Module)		
Metabolism Reagent	Genetically-encoded fluorescent ATP indica	ator for direct meas	surements of cy	tosolic ATP in living cells.		
	New! Incucyte® CytoATP Lentivirus	One vial: 0.2mL	4772	SX5 (configured with SX5 Metabolism Optical Module)		

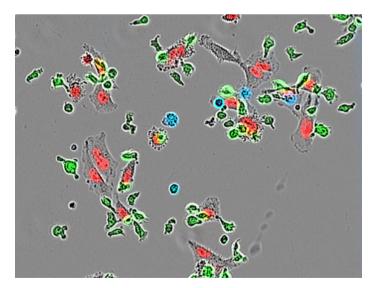
Immune Cell Activation, Killing, and NETosis

Incucyte® Immune Cell Assays are an integrated solution for real-time visualization and automated analysis of a range of immune cell functions from T-cell activation and killing to programmed neutrophil cell death.

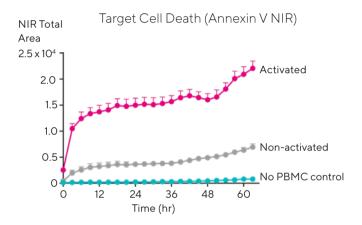
- Derive meaningful data with sensitive, non-perturbing reagents and HD phase images.
- Make multiplexed measurements of death, viability and proliferation in 2D or 3D in vitro assay models.
- Visualize and quantify dynamic cell interactions overtime in complex co-cultures.

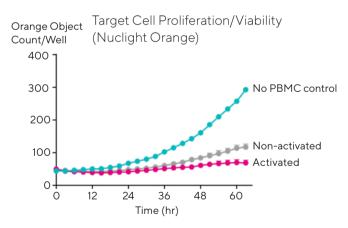
Application Spotlight: Incucyte® Immune Cell Killing

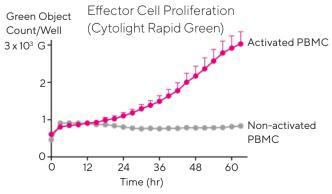
Capture, visualize and automatically quantify dynamic biological changes and cellular interactions of immune cell-mediated killing of tumor cells. With Incucyte cell health and proliferation reagents you can make multiplexed measurements of tumor cell death (Annexin V NIR Dye), tumor cell proliferation/viability (Incucyte Nuclight Orange Lentivirus) and immune cell health (Incucyte Cytolight Rapid Green Dye) in the same population of cells overtime.



MDA-MB-231 adenocarcinoma cells transduced with Incucyte Nuclight Orange co-cultured with Incucyte Cytolight Rapid Green labeled activated or non-activated PBMCs in the presence of Incucyte Annexin V NIR. Quantification of NIR (pseudo-colored blue) fluorescence area indicates target cell death and object count of orange (pseudo-colored red) fluorescence, target cell proliferation/viability (pseudo-colored green). Effector cell proliferation was quantified based on green object count over the course of the experiment.







	Product	Description	Cat. No.	Instrument Compatibility		
Software	Perform label-free cell counts and subseque and cell surface protein expression for evalu subsets of cells in complex immune-cell killi	ation of immune cel				
	Incucyte® Cell-by-Cell Analysis Software Module	1 module	9600-0031	SX5, S3, S3 for Neuroscience, SX1		
Nuclear Lentivirus Labeling Reagents	Lentivirus reagents provide homogenous ex function for live-cell quantification of cell pr			rescent protein without altering cell		
	Incucyte® Nuclight Green Lentivirus	0.2 mL	4624	SX5, S3, SX1		
	(puro)	0.6 mL	4475	SX5, S3, SX1		
	Incucyte® Nuclight Red Lentivirus (puro)	0.2 mL 0.6 mL	4625 4476	SX5 (configured with green/red optica module), S3, SX1		
	Incucyte® Nuclight Green Lentivirus	0.2 mL	4626	SX5, S3, SX1		
	(bleo)	0.6 mL	4477	SX5, S3, SX1		
	Incucyte® Nuclight Red Lentivirus	0.2 mL	4627	_ SX5 (configured with green/red optica		
	(bleo)	0.6 mL	4478	module), S3, SX1		
	Incucyte® Nuclight Orange Lentivirus (puro)	0.2 mL	4771	SX5, S3 for Neuroscience		
	Incucyte® Nuclight NIR Lentivirus (puro)	0.2 mL	4805	SX5, S3 for Neuroscience		
Cytoplasmic Dye Labeling Reagents	Live-cell cytoplasmic labeling dyes that free into a cell membrane-impermeant form, pro effector cells.					
	Incucyte® Cytolight Rapid Green Dye	One vial: 15 µg	4705	SX5, S3, SX1		
	Incucyte® Cytolight Rapid Red Dye	Five vials: 50 μg	4706	SX5 (configured with green/red optica module), S3, SX1		
	Incucyte® Cytolight Rapid Orange Dye	One vial: 1 mg	4839	SX5, S3 for Neuroscience		
Caspase Activity Reagents	Inert, non-fluorescent (DEVD) substrates that freely cross the cell membrane where they can be cleaved by activated caspase-3/7 to release a DNA-binding fluorescent label. Recommended for quantifying apoptosis in adherent target cells					
	Incucyte® Caspase-3/7 Green Dye	One vial: 20 μL (100-200 tests)	4440	SX5, S3, SX1		
	Incucyte® Caspase-3/7 Red Dye	One vial: 20 μL (100-200 tests)	4704	SX5 (configured with green/red optica module), S3, SX1		
Apoptosis Plasma Membrane Integrity	Membrane impermeable, highly-selective pextracellular surface of cells undergoing apo					
Reagents	Incucyte® Annexin V Green Dye	One vial: 100-200 tests	4642	SX5, S3, SX1		
	Incucyte® Annexin V Red Dye	One vial: 100-200 tests	4641	SX5 (configured with green/red optica module), S3, SX1		
	Incucyte® Annexin V Orange Dye	One vial: 100-200 tests	4759	SX5, S3 for Neuroscience		
	Incucyte® Annexin V NIR Dye	One vial: 100-200 tests	4768	SX5, S3 for Neuroscience		
Cytotoxicity Reagents	Highly sensitive cyanine-based dyes enter the binding to DNA. Dye does not enter cells with NETosis as extracellular DNA is released and	th intact cell membr	anes. Allows for	rapid visualization and quantification of		
	Incucyte® Cytotox Green Dye	Five vials: 5 μL (100 tests each)	4633	SX5, S3, SX1		
	Incucyte® Cytotox Red Dye	Five vials: 5 μL (100 tests each)	4632	SX5 (configured with green/red optica module), S3, SX1		
	Incucyte® Cytotox NIR Dye	One vial: 100 uL (500-100 tests)	4846	SX5, S3 for Neuroscience		

 $^{{}^\}star \text{Pre-labeled Nuclight cell lines are also available for purchase. Please visit \textbf{shop.incucyte.com} for more information.}$

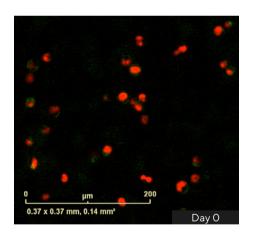
Antibody Internalization and Immunocytochemistry

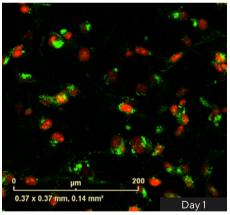
Incucyte's antibody labeling reagents are novel fluorescently labeled Fabs that can be mixed with Fc-containing antibodies and applied directly to living cells for long-term monitoring of spatial and temporal protein dynamics.

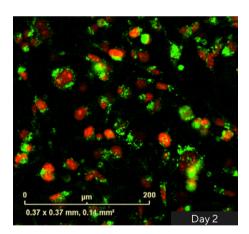
- Increase productivity with rapid single-step labeling paired with mix-and-read protocols for efficient testing of antibody panels.
- Associate changes in surface protein expression or antibody internalization with cell function and morphology over time.
- Combine sensitive, kinetic fluorescent measurement of protein dynamics with images and movies for visual confirmation of biology in every well.

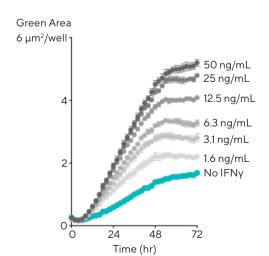
Application Spotlight: Monitoring Dynamic Cell Surface Protein Expression

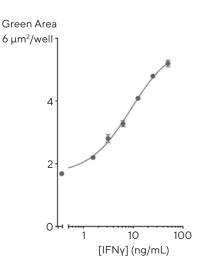
Quantify cell surface protein expression and distribution in live-cells to study long-term protein dynamics and their relationship to function and morphology using Incucyte® Fabfluor-488 Antibody Labeling Reagents.







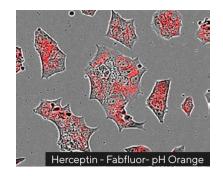


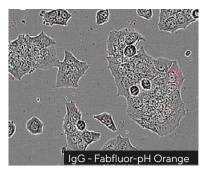


Incucyte Fabfluor-488 was conjugated to anti-PD-L1 Ab (BioLegend) and added to Nuclight Red MDA-MB-231 breast cancer cells in the absence and presence of IFN-γ (+ Incucyte® Opti-Green background suppressor). Quantification of the green fluorescent area shows that IFN-γ induces a time- and concentration- dependent increase in PD-L1 expression.

Application Spotlight: Antibody Internalization

Efficiently evaluate the full-time course of antibody internalization for real-time analysis of internalization rates under physiological conditions using Incucyte® Fabfluor-pH Antibody Labeling Reagents.



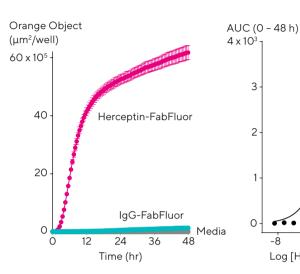


Log [Herceptin] (g/mL)

3

2

HD phase and orange fluorescence images (10X) show HER-2 positive BT-474 cells treated with Incucyte® Fabfluor-pH Orange labeled Herceptin display orange (pseudo-colored red), cytosolic fluorescence while cells treated with an isotype control display no cellular fluorescence. Time-course data shows a rapid increase in orange object area over time in cells treated with labeled Herceptin, but not with IgG1 isotype control.



	Product	Description	Cat. No.	Instrument Compatibility		
Software	Perform label-free cell counts and subseque quantify dynamic changes in cell subsets wit	•		The state of the s		
	Incucyte [®] Cell-by-Cell Analysis Software Module	1 module	9600-0031	SX5, S3, S3 for Neuroscience, SX1		
Fabfluor-pH Antibody Labeling	Novel pH-sensitive Fc-targeting antibody fra antibody internalization.	agment, fluorescent	tly label your ant	tibody of choice for real-time analysis of		
Reagents	Incucyte® Human Fabfluor-pH Orange Antibody Labeling Dye	One vial: 50 μg	4812	SX5, S3 for Neuroscience		
	Incucyte® Human Fabfluor-pH Red Antibody Labeling Dye	One vial: 50 μg	4722	SX5 (configured with green/red optica module), S3, SX1		
	Incucyte® Mouse IgG1 Fabfluor-pH Red Antibody Labeling Dye	One vial: 50 μg	4723	SX5 (configured with green/red optica module), S3, SX1		
	Incucyte® Mouse IgG2a Fabfluor-pH Red Antibody Labeling Dye	One vial: 50 μg	4750	SX5 (configured with green/red optica module), S3, SX1		
	Incucyte® Mouse IgG2b Fabfluor-pH Red Antibody Labeling Dye	One vial: 50 μg	4751	SX5 (configured with green/red optica module), S3, SX1		
	Incucyte® Rat Fabfluor-pH Red Antibody Labeling Dye	One vial: 50 μg	4737	SX5 (configured with green/red optica module), S3, SX1		
Fabflour Live-Cell	Novel fluorescently tagged Fc-targeting Fab fragments label your antibody of choice for cell surface protein expression.					
Immunocytochemistry Labeling reagents	Incucyte® Mouse IgG2a Fabfluor-488 Antibody Labeling Dye	One vial: 50 ug	4743	SX5, S3, SX1		
	Incucyte® Mouse IgG2b Fabfluor-488 Antibody Labeling Dye	One vial: 50 ug	4744	SX5, S3, SX1		
	Incucyte® Mouse IgG1 Fabfluor-488 Antibody Labeling Dye	One vial: 50 ug	4745	SX5, S3, SX1		

Phagocytosis

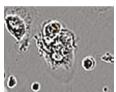
pHrodo® Reagents for Incucyte® are formulated for longterm and sensitive analysis of phagocytosis and efferocytosis of the same population of cells overtime.

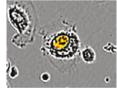
- Monitor and measure phagocytosis and efferocytosis using highly specific and sensitive labeling reagents, amenable to both adherent and non-adherent cell models.
- Study the model of your choice overtime with simple mix-and-read 96- or 384-well protocols.
- Confirm phagocytosis signals using HD phase-contrast and fluorescent images.

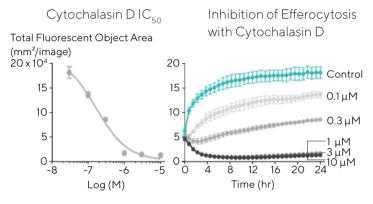
Application Spotlight: Efferocytosis Assay

Observe and measure efferocytosis over the entire assay time course using your choice of target cells labeled with pHrodo® Orange Cell Labeling Dye for Incucyte® (pseudo-colored orange, acquired with the Incucyte® S3 for Neuroscience).









Time-lapse visualization of iPSC-derived microglia (Axol BioSciences) engulfing pHrodo Orange labeled apoptotic Neuro-2A cells. Images verify the entry of an apoptotic target cell into the cytoplasm of the microglia. Inhibition of this engulfment by Cytochalasin D was quantified based on orange fluorescence object area over the course of the experiment.

	Product	Description	Cat. No.	Instrument Compatibility
pHrodo® Cell Labeling Kits	Fluorescent dyes label whole cells with a plue mediated cellular phagocytosis.	H-sensitive fluoroph	ore for real-tim	e analysis of efferocytosis and antibody
	pHrodo® Orange Cell Labeling Kit for Incucyte®	One kit	4766	SX5, S3 for Neuroscience
	pHrodo® Red Cell Labeling Kit for Incucyte®	One kit	4649	SX5 (configured with green/red optical module), S3, SX1
pHrodo® Bioparticles	Sterile fluorogenic reagents ideally suited t	o a simple mix-and-	read, real-time	live cell quantification of phagocytosis.
	pHrodo® Red E. coli Bioparticles® for Incucyte®	One vial: 2 mg	4615	SX5 (configured with green/red optical module), S3, SX1
	pHrodo®® Green E. coli Bioparticles® for Incucyte®	One vial: 2 mg	4616	SX5, S3, SX1
	pHrodo® Red Zymosan Bioparticles® for Incucyte®	One vial: 1 mg	4617	SX5 (configured with green/red optical module), S3, SX1
	pHrodo® Green Zymosan Bioparticles® for Incucyte®	One vial: 1 mg	4618	SX5, S3, SX1
	pHrodo® Red S. aureus Bioparticles® for Incucyte®	One vial: 2 mg	4751	SX5 (configured with green/red optical module), S3, SX1
	pHrodo® Green S. aureus Bioparticles® for Incucyte®	One vial: 2 mg	4737	SX5, S3, SX1

Angiogenesis

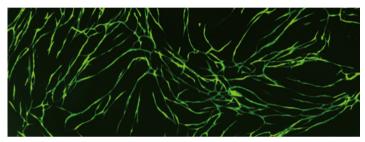
The Incucyte® Angiogenesis Primekit and associated software are a fully integrated solution to measure vascular tube formation in a relevant co-culture model over time.

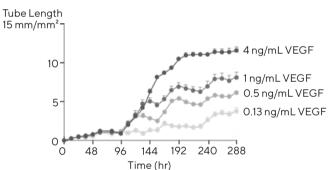
- Study all phases of the angiogenic process with an optimized, ready-to-use 96-well cryopreserved kit.
- Increase insight with automated, kinetic data analysis and visualization of tube formation and disruption in real-time
- Characterize the angiogenic effects of vascular growth factors on tube formation over the complete time course of the assay.

Application Spotlight: Angiogenesis Assay

Assess complex network development and disruption quantitatively and qualitatively with the Incucyte® Angiogenesis Primekit and Incucyte® Angiogenesis Analysis Software Module to study therapeutic interventions of vascular formation.

Concentration dependent vascular tube formation in Primekit Cytolight Green HUVEC and human dermal fibroblast co-culture model. (Top) Example image of VEGF (4 ng/ml) induced vascular structures. (Bottom) Concentration dependent vascular tube formation induced by Vasoactive Endothelial Growth Factor (VEGF) in Primekit Cytolight Green HUVEC and human dermal fibroblast co-culture model.





	Product	Description	Cat. No.	Instrument Compatibility				
Software	Assess complex network development and of vascular formation.	Assess complex network development and disruption quantitatively and qualitatively to study therapeutic interventions of vascular formation.						
	Incucyte® Angiogenesis Analysis Software Module	1 module	9600-0011	SX5, S3, SX1				
Angiogenesis Reagents	Cryopreserved kit of lentivirally-infected H (NHDF) for the study of angiogenic networ	, ,	FP cultured wit	h normal human dermal fibroblasts				
	Incucyte® Angiogenesis Primekit	One kit	4452	SX5, S3, SX1				
	Incucyte® Angiogenesis Primekit VEGF/Suramin Kit	Two vials	4437	SX5, S3, SX1				
	Incucyte® Angiogenesis Primekit Assay Media Kit	One bottle: 125 mL	4541	SX5, S3, SX1				
	Incucyte® Cytolight Green HUVEC	One vial: 1.7 x 10 ⁶ cells	4453	SX5, S3, SX1				
	Incucyte® Cytolight Green Lentivirus (no selection)	One vial: 0.6 mL	4513	SX5, S3, SX1				

Neurite Dynamics and Neuronal Activity

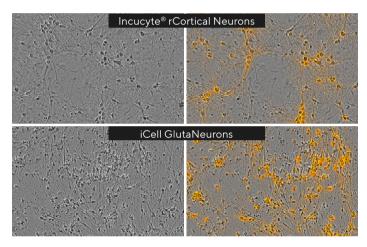
Incucyte's novel, live-cell labeling reagents and purposebuilt software quantify long-term changes in neurite dynamics and neuronal activity that enable continuous analysis of sensitive neuronal cell models.

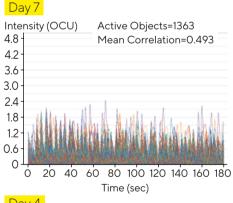
- Conduct long term studies of neuronal function with novel, non-perturbing fluorescent reagents.
- Capture transient events in your choice of cell model with non-invasive, repeated measurements of the same neuronal culture in physiologically relevant conditions.
- Analyze relevant morphological and functional metrics using intuitive, purpose-built Incucyte® software.

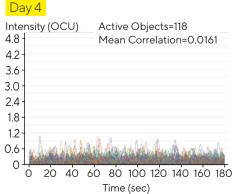
Application Spotlight: Neuronal Activity Assay

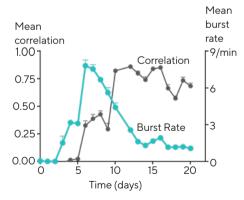
Access complex, neuronal activity and connectivity measurements from thousands of cells chronically to gain unprecedented functional insight into neuronal cell models using our novel Incucyte® Neuroburst Orange Lentivirus and Incucyte® Neuronal Activity Analysis Software Module.

Incucyte® rCortical Neurons and iCell GlutaNeurons (Cellular Dynamics International) express the Incucyte Neuroburst Orange Lentivirus, without perturbing the health and morphology of the cells.

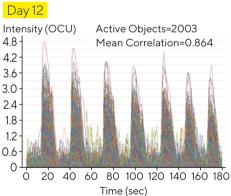


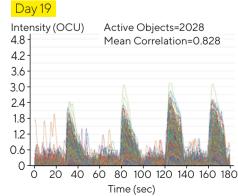






Kinetic quantification (center graph) of longitudinal, dynamic changes in neuronal activity mean burst rate and mean correlation) of iCellGluta Neurons expressing Neuroburst Orange Lentivirus, showing changes over time during neuronal network maturation. Active object traces (corner traces) provide detailed insight into the dynamic changes in neuronal activity and connectivity for every acquired timepoint.

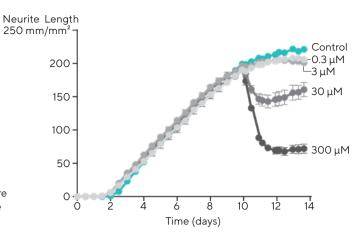


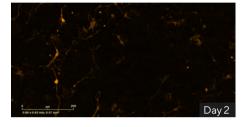


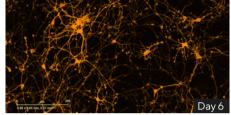
Application Spotlight: Neurite Dynamics Assay

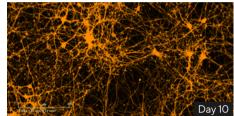
Generate kinetic, image-based and automated measurements using Incucyte® Neurotrack Analysis Software Module for continuous analysis of neurite outgrowth and stability – inside your incubator.

Incucyte® rCortical Neurons transduced with Incucyte® Neurolight Orange Lentivirus labeled cells cultured in the presence of Incucyte® rAstrocytes were treated with glutamate at day 10. Time course analysis of orange fluorescence neurite length reveals concentration-dependent treatment effects.









	Product	Description	Cat. No.	Instrument Compatibility			
Neuronal Activity	Record activity from over a thousand cells t	o study changes in ne	euronal network	activity and connectivity.			
Software	Purpose-built acquisition and analysis softw	vare for the detection	of calcium osc	illations in 96-well plates.			
	Incucyte® Neuronal Activity Analysis Software Module	1 module	9600-0032	SX5, S3 for Neuroscience			
Neuronal Activity	Fluorescently detect changes in activity usi	ng a novel genetically	/-encoded fluo	rescent calcium indicator.			
Reagents	Incucyte® Neuroburst Orange Lentivirus	One vial: 2 mL	4736	SX5, S3 for Neuroscience			
	Incucyte® Neuroactive Orange Kit	One kit	4761	SX5, S3 for Neuroscience			
Neurite Dynamics	Characterize neurite dynamics over time in mono- or culture models while assessing cell viability (refer to page 7, Annexin V Reagents)						
Software	Enables label-free or fluorescent analysis or well plates.	f neurite outgrowth, n	naturation and	disruption in each well of a 96- or 384-			
	Incucyte [®] Neurotrack Analysis Software Module	1 module	9600-0010	SX5, S3, SX1			
Neurite Labeling Reagents		Lentivirus reagents driven off a synapsin promoter provide homogenous expression of a fluorescent protein in target cells without altering cell function for live-cell quantification of neurite outgrowth.					
	Incucyte® Neurolight Orange Lentivirus	Two vials: 0.45 mL each	4808	SX5, S3 for Neuroscience			
	Incucyte® Neurolight Red Lentivirus	Two vials: 0.45 mL each	4807	SX5 (configured with green/red optical module), S3, SX1			
	Incucyte® Neuroprime Orange Kit	One kit	4760	SX5, S3 for Neuroscience			
	Incucyte® Neuroprime Red Kit	One kit	4585	SX5 (configured with green/red optical module), S3, SX1			
Neuronal Cells	Read-to-use cryopreserved cells from from the cortex of Sprague Dawley Rats at day 18 of gestation.						
	Incucyte® rCortical Neurons	One vial: 2 x 10° cells	4753	Assay dependent			
		2 x 10° cells					

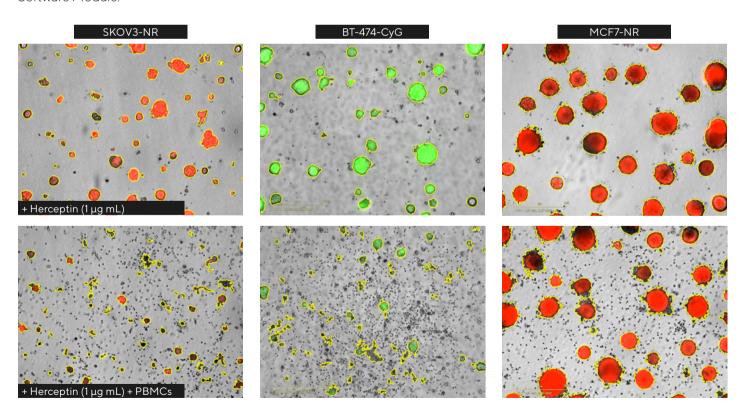
Spheroid Growth, Invasion, and Immune Cell Killing

The Incucyte® Spheroid Analysis Software Module is designed to acquire and analyze label or label-free spheroids to automatically monitor and quantify the formation, growth, shrinkage and invasive properties of advanced cell models in real time inside your tissue culture incubator.

- Flexible acquisition mode to enable studies of both single and multi-spheroid assays
- Enable long-term imaging with enhanced depth of focus Brightfield (DF® Brightfield) image acquisition
- Analyze relevant spheroid metrics using intuitive Incucyte® processing definitions

Application Spotlight: 3D Immune Cell Killing

Quantify and visualize immune cell –mediated killing of solid tumors in real time to using non-perturbing reagents and purpose-built, integrated Incucyte® Spheroid Analysis Software Module.



Tumor cells either stably expressing Incucyte Nuclight Red Lentivirus (SKOV3-NR, MCF7-NR) or Incucyte Cytolight Green Lentivirus (BT-474-CyG) were seeded on a bed of Matrigel® in flat bottom 96-well plates. Multi-spheroids were allowed to form (3 d) prior to addition of freshly isolated PBMCs (E:T, 5:1) and Herceptin. Incucyte Brightfield and fluorescence images (7 d; SKOV3-NR, MCF-NR or 10 d; BT-474-CyG) compare the effect of Herceptin on spheroid proliferation in absence (top panel) and presence (bottom panel) of PBMCs (Brightfield outline mask shown in yellow).

	Product	Description	Cat. No.	Instrument Compatibility			
Spheroid Software	Analyze growth, viability and invasion of si spheroids in flat bottom plates to detect c			i-well formats or measure multiple			
	Incucyte® Spheroid Analysis Software Module	1 module	9600-0031	SX5, S3, S3 for Neuroscience, SX1			
Nuclear Lentivirus Labeling Reagents	Lentivirus reagents provide homogenous function for live-cell quantification of cell			prescent protein without altering cell			
	Incucyte® Nuclight Green Lentivirus	0.2 mL	4624	SX5, S3, SX1			
	(puro)	0.6 mL	4475	SX5, S3, SX1			
	Incucyte® Nuclight Red Lentivirus	0.2 mL	4625	SX5 (configured with green/red optical			
	(puro)	0.6 mL	4476	module), S3, SX1			
	Incucyte® Nuclight Green Lentivirus	0.2 mL	4626	SX5, S3, SX1			
	(bleo)	0.6 mL	4477	SX5, S3, SX1			
	Incucyte® Nuclight Red Lentivirus	0.2 mL	4627	SX5 (configured with green/red optical			
	(bleo)	0.6 mL	4478	module), S3, SX1			
	Incucyte® Nuclight Orange Lentivirus (puro)	0.2 mL	4771	SX5, S3 for Neuroscience			
	Incucyte® Nuclight NIR Lentivirus Reagent (puro)	0.2 mL	4805	SX5, S3 for Neuroscience			
Apoptosis Plasma Membrane Integrity	Membrane impermeable, highly-selective phosphatidylserine (PS) cyanine fluorescent dyes label PS exposed on the extracellular surface of cells undergoing apoptosis.						
Reagents	Incucyte® Annexin V Green Dye	One vial: 100 – 200 tests	4642	SX5, S3, SX1			
	Incucyte® Annexin V Red Dye	One vial: 100 – 200 tests	4641	SX5 (configured with green/red optical module), S3, SX1			
	Incucyte® Annexin V Orange Dye	One vial: 100 – 200 tests	4759	SX5, S3 for Neuroscience			
	Incucyte® Annexin V NIR Dye	One vial: 100 – 200 tests	4768	SX5, S3 for Neuroscience			
Cytoplasmic Dye	Live-cell cytoplasmic labeling dyes that freely pass through cell membranes and into cells, where they are transformed						
_abeling Reagents	into a cell membrane-impermeant form, p	roviding spatial conte	xt for cell-to-ce	ll interactions.			
	Incucyte® Cytolight Rapid Green Dye	One vial: 15 µg	4705	SX5, S3, SX1			
	Incucyte® Cytolight Rapid Red Dye	Five vials: 50 μg	4706	SX5 (configured with green/red optical module), S3, SX1			
Cytoplasmic Lentivirus Labeling	Lentivirus reagents provide homogenous quantification of spheroid growth and shri		scent protein wi	thout altering cell function for live-cell			
Reagents	Incucyte® Cytolight Green Lentivirus (puro)	0.6 mL	4481	SX5, S3, SX1			
	Incucyte Cytolight Red Lentivirus (puro)	0.6 mL	4482	SX5 (configured with green/red optical module), S3, SX1			

 $^{{}^*\}text{Pre-labeled Nuclight cell lines are also available for purchase. Please visit \textbf{shop.incucyte.com} for more information.}$

Invasion and Migration Assays

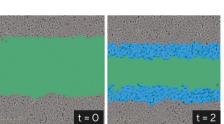
Incucyte® Scratch Wound and Chemotaxis Assays allow you to continuously monitor and analyze migration and invasion using purpose built consumables and automated software.

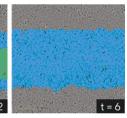
- Unlock your productivity with optimized protocols, peripherals and automated analysis
- Confirm cell movement and morphology with images and movies – all from inside your incubator
- Analyze relevant spheroid metrics using intuitive Incucyte® processing definitions

Application Spotlight: Scratch Wound Migration and Invasion Assay

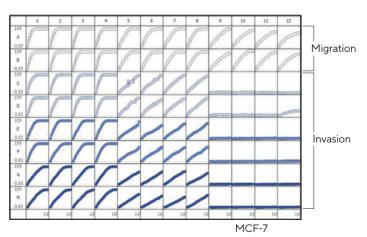
Make label-free, quantitative measurements of migration and invasion, while visualizing morphological changes due treatment effects using the integrated Incucyte® Scratch Wound Analysis Software Module.



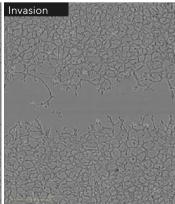




Incucyte® 96-well Woundmaker Tool creates 96 precise, uniform cell-free zones with the touch of a button in cell monolayers cultured in our Incucyte Imagelock plates. Wound closure is visualized and analyzed in real-time with the Incucyte® Scratch Wound Analysis Software Module.



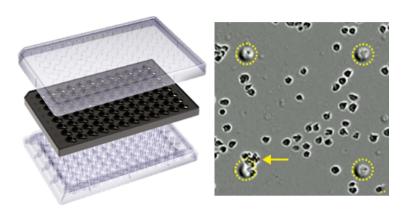




Microplate graph showing progression of differential cell migration and invasion of three different cell types using the Relative Wound Density metric. Representative images of HT-1080 fibrosarcoma cells highlights differences in morphology of cells during migration versus invasion.

Application Spotlight: Incucyte® Chemotaxis Assay

Maximize your productivity and insight into dynamic cell movement with automated, label-free imaging and analysis using optically clear Incucyte® Clearview 96-Well Chemotaxis Plates and Incucyte® Chemotaxis Analysis Software Module.



Total T-cell area normalized to initial value 1.2 —Control —3.125 nM —6.25 nM —12.5 nM —50 nM

20

Time (h)

-200 nM -100 nM

30

Chemotaxis of Primary T-Cells to SDF-1a

Incucyte Clearview 96-well Chemotaxis Plates provide an optically clear surface for label-free imaging and analysis of chemotactic cell migration or invasion. Cells are added to the upper chamber and chemoattractant to the lower reservoir plate. Chemotactic transmembrane migration is automatically quantified as the cells migrate through laser etched pores (yellow circles) toward chemoattractant. Example data of concentration-dependent SDF-1 α mediated chemotaxis of CD3/CD28 activated human T-cells.

0.4

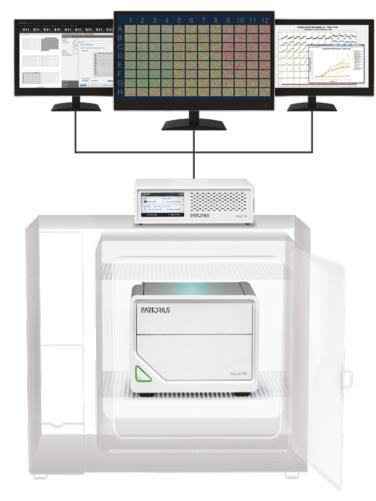
0.0

10

	Product	Description	Cat. No.	Instrument Compatibility			
Scratch Wound	Analyze 96-well Incucyte® Scratch Wound Assay images to analyze cell migration and invasion with or without labels.						
Cell Migration and Invasion	Incucyte® Scratch Wound Analysis Software Module	1 module	9600-0012	SX5, S3, SX1			
	Incucyte® 96-well Woundmaker Tool	1 wounder	4563	SX5, S3, SX1			
	Incucyte® Cell Migration Bundle	1 Scratch Wound Analysis Software Module 196-well Woundmaker Tool	4493	SX5, S3, SX1			
	Incucyte® Cell Migration/Invasion Bundle	1 Scratch Wound Analysis Software Module 196-well Woundmaker Tool 1 Biocision CoolBox Kit	4474	SX5, S3, SX1			
	Incucyte® Imagelock 96-well Plate	Pack of 10 plates	4806	SX5, S3, SX1			
		Pack of 50 plates	4379	SX5, S3, SX1			
Chemotaxis Cell	Track and quantify label-free and fluorescently labeled chemotaxis cell migration and invasion in microplate format.						
Migration and Invasion	Incucyte® Chemotaxis Analysis Software Module	1 module	9600-0015	SX5, S3, SX1			
	Incucyte® Clearview 96-Well Chemotaxis	1 plate	4582	SX5, S3, SX1			
	Plates	Pack of 10 plates	4648	SX5, S3, SX1			

From our workhorse, the Incucyte[®] S3, to our economical Incucyte® SX1, to the newest member, the frontier-breaking Incucyte® SX5, we deliver real-time insights to simplify progress for labs of any size.

Analyze your cells for days, weeks or even months as they sit stationary in the stable environment of your tissue culture incubator. With Incucyte's user-friendly interface and robust instrument portfolio, any cell biologist can gain dynamic insights into the health, morphology, movement and function of their cell models



Support an entire research team via concurrent, remote network access. Make real-time decisions from the comfort of your office or home.

Sales and Service Contacts

For further contacts, visit www.sartorius.com/incucyte



facebook.com/essenbioscience/



linkedin.com/company/essen-bioscience/



twitter.com/EssenBioScience

North America

Phone: +17347691600, ext. 1 E-Mail: orders.US07@sartorius.com

Europe

Phone: +44 (0) 1763 227400

E-Mail: euorders.UK03@sartorius.com

Japan

Phone: +81364785202

E-Mail: orders.US07@sartorius.com

China

Phone: +86 21 6878 2300

Rest of APAC and other countries around the world:

Phone: +65 6872 3966