



## Protocol

### IncuCyte® Label-Free Cell Proliferation Assay

For counting and confluence measurements of adherent or non-adherent cell lines

This protocol provides an overview of the IncuCyte® Cell Proliferation Assay methodology. It is compatible with the IncuCyte® Live-Cell Analysis System for kinetic, label-free analysis of cell confluence or cell counts using your choice of cells and treatments.

The highly flexible assay format can be combined with our range of IncuCyte® cell health and viability reagents for multiplexed measurements of cytotoxicity and apoptosis alongside proliferation in the same well.

#### Required materials

- Flat bottom tissue culture plate (e.g., Corning Cat. No. 3595)

#### Optional materials

- IncuCyte® Cell-By-Cell Analysis Software Module (Sartorius Cat. No. 9600-0031)—for label-free cell counting
- Poly-L-ornithine (Sigma Cat. No. P4957)—for non-adherent cells
- Fibronectin (Sigma Cat. No. A7906)—for non-adherent cells

#### General guidelines

- Following cell seeding, place plates at ambient temperature (15 minutes for adherent cell lines and 45 minutes for non-adherent cell lines) to ensure homogenous cell settling.
- Remove bubbles from all wells by gently squeezing a wash bottle containing 70–100% ethanol with the inner straw removed, to blow vapor over the surface of each well.
- After placing the plate in the IncuCyte Live-Cell Analysis System, allow the plate to warm to 37° C for 30 minutes prior to scanning.

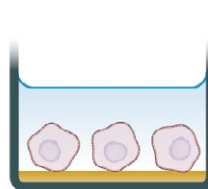
#### Quick guide

##### 1. Coat wells



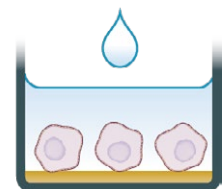
Coat wells of plate (50  $\mu$ L/well) with appropriate matrix. Optional for adherent cell lines.

##### 2. Plate cells



Seed cells (100  $\mu$ L/well, 1,000–10,000 for adherent and 5,000–50,000 for non-adherent) into a 96-well plate.

##### 3. Add treatments



Add desired treatments (100  $\mu$ L/well, 1x for adherent cell lines, 2x for non-adherent cell lines).

## Adherent cell line protocol

### Day 0

#### 1. Coat wells (optional)

- 1.1 Depending on cell line used, coat a 96-well flat bottom plate with relevant coating matrix according to manufacturer's recommendation.
- 1.2 Prior to cell seeding, prepare cell treatments at 2x final assay concentration in enough cell culture medium to achieve a volume of 100  $\mu$ L per well.

#### 2. Plate cells

- 2.1 Seed your choice of cells (100  $\mu$ L per well) at an appropriate density into a 96-well plate, such that by day 1 the cell confluence is approximately 10–20%. The seeding density will need to be optimized for the cell line used; however, we have found that 1,000 to 2,500 cells per well (10,000 to 25,000 cells/mL seeding stock) are reasonable starting points.
  - a. Monitor cell growth using the IncuCyte® Live-Cell Analysis System to capture phase contrast images every two hours and analyze using the integrated confluence algorithm.

### Day 1

#### 3. Add treatments

- 3.1 Prepare 1x concentrations of desired cell treatments in cell culture medium. The volumes may be varied; however, we recommend preparing enough volume of each desired treatment/dilution in order to achieve 100  $\mu$ L per well.
- 3.2 Remove the cell plate from the incubator and aspirate medium from wells.
- 3.3 Add treatments and controls to appropriate wells of the 96-well plate.
- 3.4 Place the cell plate into the IncuCyte Live-Cell Analysis System and allow the plate to warm to 37° C for 30 minutes prior to scanning.
  - a. Objective: 4x, 10x or 20x
  - b. Channel selection: Phase Contrast (+ "Green" or "Red" if fluorescent label or cell health reagents are used)
  - c. Scan type: Standard
  - d. Scan interval: Typically, every 1 to 2 hours, until your experiment is complete.

NOTE: Label free cell counting can be enabled on IncuCyte Live-Cell Analysis System with use of the IncuCyte Cell-by-Cell Analysis Software Module.

- a. Scan type: Standard/Adherent Cell-by-Cell
- b. Objective: 10x

For further details of this analysis module and its application see: [www.essenbioscience.com/cell-by-cell](http://www.essenbioscience.com/cell-by-cell)

## Non-adherent cell line protocol

### Day 1

#### 1. Seed cells and add prepared treatments

- 1.1 Coat a 96-well flat bottom plate with relevant coating matrix. We recommend coating with 50  $\mu$ L of either 0.01% poly-L-ornithine solution or 5  $\mu$ g/mL fibronectin diluted in 0.1% BSA. Coat plates for 1 hour at ambient temperature, remove solution from wells, then allow plates to dry for 30–60 minutes prior to cell addition.
- 1.2 Prior to cell seeding, prepare cell treatments at 2x final assay concentration in enough cell culture medium to achieve a volume of 100  $\mu$ L per well.

#### 2. Plate cells

- 2.1 Seed your choice of cells (100  $\mu$ L per well) at an appropriate density into a 96-well plate. The seeding density will need to be optimized for the cell line used; however, we have found that 5,000 to 50,000 cells per well (50,000 – 500,000 cells/mL seeding stock) are reasonable starting points.

NOTE: If studying immune cell clustering and proliferation, prepare cell activation treatments at 5x final concentration, and immediately add 50  $\mu$ L per well containing cells. It is advised that some control wells containing only vehicle are included in the plate.

#### 3. Add treatments

- 3.1 Immediately after cell seeding, add treatments and controls to appropriate wells of the 96-well plate containing cells. Triturate wells to appropriately mix the treatment to ensure cell exposure at 1x.
- 3.2 Place the cell plate into the IncuCyte Live-Cell Analysis System and allow the plate to warm to 37° C for 30 minutes prior to scanning.
  - a. Objective: 4x (recommended 1 image per well or whole well) or 10x
  - b. Channel selection: Phase Contrast (+ "Green" or "Red" if fluorescent label or cell health reagents are used)
  - c. Scan type: Standard
  - d. Scan interval: Typically, every 1 to 2 hours, until your experiment is complete.

NOTE: Label-free cell counting can be enabled on IncuCyte Live-Cell Analysis System with use of the IncuCyte Cell-by-Cell Analysis Software Module.

- a. Scan type: Standard/Non-Adherent Cell-by-Cell
- b. Objective: 20x

For further details of this analysis module and its application see: [www.essenbioscience.com/cell-by-cell](http://www.essenbioscience.com/cell-by-cell)

## Related products and applications:

A comprehensive range of fluorescent nuclear labeling and cell health reagents are available for use with the IncuCyte Live-Cell Analysis System to enable multiplexed measurements of cytotoxicity and proliferation alongside apoptosis.

### Compatible with the IncuCyte® Live-Cell Analysis System Cat. No. 4647

Product	Cat. No.	Amount
IncuCyte® NuLight Green Lentivirus Reagent (EF-1 $\alpha$ , Puro) for nuclear labeling	4624	0.2 mL
IncuCyte® NuLight Red Lentivirus Reagent (EF-1 $\alpha$ , Puro) for nuclear labeling	4625	0.2 mL
IncuCyte® NuLight Green Lentivirus Reagent (EF-1 $\alpha$ , Bleo) for nuclear labeling	4626	0.2 mL
IncuCyte® NuLight Red Lentivirus Reagent (EF-1 $\alpha$ , Bleo) for nuclear labeling	4627	0.2 mL
IncuCyte® NuLight Green Lentivirus Reagent (EF-1 $\alpha$ , Puro) for nuclear labeling	4475	0.6 mL
IncuCyte® NuLight Red Lentivirus Reagent (EF-1 $\alpha$ , Puro) for nuclear labeling	4476	0.6 mL
IncuCyte® NuLight Green Lentivirus Reagent (EF-1 $\alpha$ , Bleo) for nuclear labeling	4477	0.6 mL
IncuCyte® NuLight Red Lentivirus Reagent (EF-1 $\alpha$ , Bleo) for nuclear labeling	4478	0.6 mL
IncuCyte® Cytotox Red Reagent for counting dead cells	4632	5 $\mu$ L x 5
IncuCyte® Cytotox Green Reagent for counting dead cells	4633	5 $\mu$ L x 5
IncuCyte® Annexin V Red Reagent for apoptosis	4641	100 tests
IncuCyte® Annexin V Green Reagent for apoptosis	4642	100 tests
IncuCyte® Caspase-3/7 Red Reagent for apoptosis	4704	20 $\mu$ L
IncuCyte® Caspase-3/7 Green Reagent for apoptosis	4440	20 $\mu$ L

### Compatible with the IncuCyte® Live-Cell Analysis System for Neuroscience Cat. No. 4763

Product	Cat. No.	Amount
IncuCyte® NuLight Orange Lentivirus Reagent (EF-1 $\alpha$ , Puro) for nuclear labeling	4771	0.2 mL
IncuCyte® Annexin V Orange Reagent for apoptosis	4760	100 tests
IncuCyte® Annexin V NIR Reagent for apoptosis	4768	100 tests

A complete suite of cell health applications is available to fit your experimental needs. Find more information at [www.sartorius.com/incucyte](http://www.sartorius.com/incucyte)

For additional product or technical information, please email us at [AskAScientist@sartorius.com](mailto:AskAScientist@sartorius.com)

Essen Bioscience,  
a Sartorius Company  
300 West Morgan Road  
Ann Arbor, Michigan, 48108  
USA

[www.sartorius.com/incucyte](http://www.sartorius.com/incucyte)  
Email: [AskAScientist@sartorius.com](mailto:AskAScientist@sartorius.com)

USA +1.734.769.1600  
UK +44.1707.358688  
Japan +81.3.5826.4795