

Protocol

IncuCyte® Cell Proliferation Assay

For label-free proliferation 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® S3 Live-Cell Analysis System 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 3595)
- Poly-L-ornithine (Sigma P4957) – optional, for non-adherent cells
- Fibronectin (Sigma A7906) – optional, 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 S3 Live-Cell Analysis System, allow the plate to warm to 37 °C for 30 minutes prior to scanning.

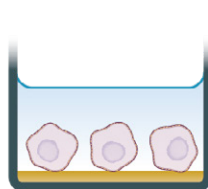
Quick Guide

1. Coat wells



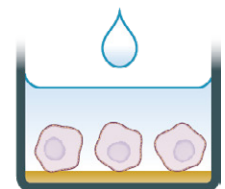
Coat wells of plate (50 µL/well) with appropriate matrix. Optional for adherent cell lines.

2. Plate cells



Seed cells (100 µL/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 µL/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 μ L per well.

2. Plate Cells

- 2.1 Seed your choice of cells (100 μ 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® S3 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 μ 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 S3 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 S3 Live-Cell Analysis System with use of the IncuCyte Cell-by-Cell Analysis Software Module (PN 9600 0031).

- 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

Non-Adherent Cell Line Protocol

For label-free proliferation measurements of adherent or non-adherent cell lines

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 μ L of either 0.01% poly-L-ornithine solution or 5 μ 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 μ L per well.

2. Plate Cells

- 2.1 Seed your choice of cells (100 μ 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 μ 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 S3 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 S3 Live-Cell Analysis System with use of the IncuCyte Cell-by-Cell Analysis Software Module (PN 9600 0031).

- 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

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.

Product	Cat No.	Amount
IncuCyte® NuLight Green BacMam 3.0 Reagent for nuclear labeling	4622	1 mL
IncuCyte® NuLight Green Lentivirus Reagent (EF-1 α , Puro) for nuclear labeling	4624	0.2 mL
IncuCyte® NuLight Red Lentivirus Reagent (EF-1 α , Puro) for nuclear labeling	4625	0.2 mL
IncuCyte® NuLight Green Lentivirus Reagent (EF-1 α , Bleo) for nuclear labeling	4626	0.2 mL
IncuCyte® NuLight Red Lentivirus Reagent (EF-1 α , Bleo) for nuclear labeling	4627	0.2 mL
IncuCyte® NuLight Green Lentivirus Reagent (EF- α , Puro) for nuclear labeling	4475	0.6 mL
IncuCyte® NuLight Red Lentivirus Reagent (EF-1 α , Puro) for nuclear labeling	4476	0.6 mL
IncuCyte® NuLight Green Lentivirus Reagent (EF-1 α Bleo) for nuclear labeling	4477	0.6 mL
IncuCyte® NuLight Red Lentivirus Reagent (EF-1 α , Bleo) for nuclear labeling	4478	0.6 mL
IncuCyte® Cytotox Red Reagent for counting dead cells	4632	5 μ L x 5
IncuCyte® Cytotox Green Reagent for counting dead cells	4633	5 μ 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 Green Reagent for apoptosis	4440	20 μ L
IncuCyte® Cell-by-Cell Analysis Software Module	9600-0031	1 Module

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

For additional product or technical information, please e-mail us at AskAScientist@sartorius.com visit our website at [essenbioscience.com](https://www.essenbioscience.com)

or call

+1-734-769-1600 (USA)

+44-1707-358688 (Europe)

+81-3-5579-6200 (Japan)