

IncuCyte[®] Cytotoxicity Assay

For the fluorescent quantification of cell death

This protocol provides an overview of the IncuCyte[®] Cytotoxicity Assay methodology which uses the mix-and-read IncuCyte[®] Green or Red Reagent to detect cell death in real time. The protocol is compatible with the IncuCyte[®] live-cell analysis system and your choice of cells (e.g., tumor, immune, neuronal) and treatments.

Furthermore, this protocol can be used with cells labeled using the IncuCyte[®] NuLight nuclear labeling reagents to provide multiplexed measurements of proliferation alongside cell death in the same well.

Required materials

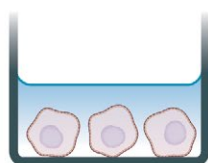
- IncuCyte[®] Red Cytotoxicity Reagent (EssenBioscience Cat #4632)
or
- IncuCyte[®] Green Cytotoxicity Reagent (EssenBioscience Cat #4633)
- Poly-L-ornithine (Sigma P4957)
 - optional, for non-adherent cells
- Fibronectin (Sigma F1141)
 - optional, for non-adherent cells
- Flat bottom tissue culture plate (e.g., Corning 3595, TPP 92096 for neuronal cell health)

General Guidelines

- We recommend medium with low levels of riboflavin to reduce the green fluorescence background. EBM, F12-K, and Eagles MEM have low riboflavin (<0.2 mg/L). DMEM and RPMI have high riboflavin (>0.2 mg/L).
- 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
- If monitoring cytotoxicity in primary neuronal cultures, we recommend use of the IncuCyte[®] Cytotox Red reagent to eliminate risk of green channel excitation issues in these sensitive cell types.

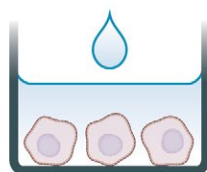
Adherent Cell Line Protocol

1 SEED CELLS



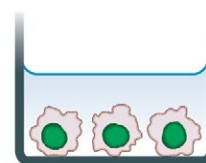
Seed cells (100 µL/well, 1,000 – 5,000) into a 96-well plate and incubate overnight.

2 PREPARE CYTOTOXICITY REAGENT AND TREAT CELLS



Prepare the desired treatments at 1x in medium containing IncuCyte Cytotoxicity reagent. Aspirate media from wells and add treatment (100 µL/well).

3 LIVE CELL FLUORESCENT ANALYSIS



Capture images every 2-3 hours (20x or 10x) in the IncuCyte[®] System. Analyze using integrated software.

Day 0:

1 Seed effector cells

1.1. Seed your choice of cells (100 µL per well) at an appropriate density into a 96-well plate, such that by day 1 the cell confluence is approximately 30%. The seeding density will need to be optimized for the cell line used; however, we have found that 1,000 to 5,000 cells per well (10,000 – 50,000 cells/mL seeding stock) are reasonable starting points.

NOTE: For non-proliferating cell lines (e.g., rate forebrain neurons) we recommend seeding at 15×10^3 to 30×10^3 cells per well, and culturing for 14 days for the neural network to establish, prior to evaluating cytotoxicity.

- a. Monitor cell growth using the IncuCyte system to capture phase contrast images every 2 hours and analyze using the integrated confluence algorithm.

Day 1:

2 Cytotoxicity reagent preparation and cell treatment addition

- 2.1. Dilute cytotoxicity reagent in desired medium formulation.
NOTE: All test agents will be diluted in this reagent-containing medium, so make up a volume that will accommodate all treatment conditions. The volumes/dilutions added to cells may be varied; however, a volume of 100 µL per well is generally sufficient for the duration of the assay.
- 2.2. Remove the cell plate from the incubator and aspirate off growth medium.
- 2.3. Add treatments and controls to appropriate wells of the 96-well plate.

3 Live-cell imaging of cytotoxicity

- 3.1. 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.
- Objective: 10x or 20x
 - Channel selection: Phase Contrast and Fluorescence
 - Scan type: Standard (2-4 images per well)
 - Scan interval: Typically, every 2 hours, until your experiment is complete.
- NOTE:** For neuronal cultures we recommend scanning every 6 to 12 hours to minimize risk of phototoxicity.

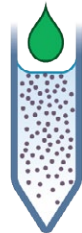
Non-Adherent Cell Line Protocol

1 COAT PLATE



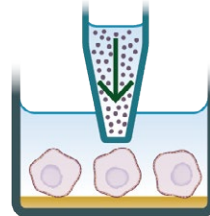
Coat plate with 0.01% poly-L-ornithine solution or 5 µL/mL fibronectin diluted in 0.1% BSA.

2 PREPARE INCUCYTE CYTOTOXICITY REAGENT AND TREATMENT



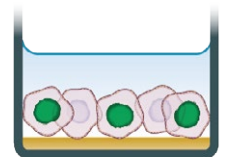
Dilute cytotoxicity reagent in medium at 1x and prepare cell treatments.

3 SEED CELLS AND ADD TREATMENT



Seed cells (100 µL/well, 5,000 – 25,000 cells) into the coated 96-well plate. Immediately add cytotoxicity ± treatments and triturate.

4 LIVE CELL FLUORESCENT ANALYSIS



Capture images every 2-3 hours (20x or 10x) in the IncuCyte® System.

Day 1:

1 Coat Plate

- 1.1. Coat a 96-well flat bottom plate with appropriate coating matrix. We recommend coating with 50 µL of either 0.01% poly-L-ornithine solution or 5 µg/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.

2 Prepare cytotoxicity reagent and treatments

- 2.1. Prior to cell seeding, dilute cytotoxicity reagent to a final concentration of 250 nM (1:4000 dilution) in desired medium formulation.
NOTE: All test agents will be diluted in this reagent-containing medium, so make up a volume that will accommodate all treatment conditions. The volumes/dilutions added to cells may be varied; however, a volume of 100 µL per well is generally sufficient for the duration of the assay.
- 2.2. Prepare cell treatments at 2x final assay concentration in enough cell culture medium the cytotoxicity reagent to achieve a volume of 100 µL per well.

3 Seed cells and add prepared treatments

- 3.1. Seed your choice of cells (100 µL per well) at an appropriate density into a 96-well plate in medium containing the cytotoxicity reagent. The seeding density will need to be optimized for the cell line used; however, we have found that 5,000 to 25,000 cells per well (50,000 – 250,000 cells/mL seeding stock) are reasonable starting points.
- 3.2. Immediately 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.

4 Live-Cell Imaging of cytotoxicity

- 4.1. 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.
- Objective: 4x (recommended 1 image per well or whole well) or 10x
 - Channel selection: Phase Contrast and Fluorescence
 - Scan type: Standard
 - Scan interval: Typically, every 2 hours, until your experiment is complete.

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 apoptosis and proliferation alongside cytotoxicity.

Product	Cat No.	Amount
IncuCyte® NuLight Red BacMam 3.0 Reagent for nuclear labeling	4621	1 mL
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-1 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

A complete suite of cell health applications is available to fit your experimental needs. Find more information at essenbioscience.com/applications

For additional product or technical information, please e-mail us at AskAScientist@essenbio.com
Visit our website at essenbioscience.com
or call
+1 734-769-1600 (USA),
+44 1707 358688 (Europe)
+81-3-5579-6200 (Japan)