

Incucyte® Multi-Spheroid Assay

For the Quantification of Multi-Spheroid Growth and Health on a Layer of Matrigel®

This protocol describes a solution for creating multi-spheroids using 96-well flat bottom plates coated with Matrigel®, and subsequent determination of cell viability, apoptosis, or cytotoxicity using Incucyte® Cell Health

Reagents. The method utilizes the Incucyte® Live-Cell Analysis System for image-based brightfield and fluorescence measurements of multi-spheroid size (area), number and health.

Required Materials

- 96-well flat bottom TC-treated microplate (Corning Cat. No. 3595)
- Matrigel®, protein concentration ≥ 8 mg/mL (Corning Cat. No. 356234)
- Wet ice
- Serum-free cell culture media for Matrigel® dilutions
- Complete culture media for cell culture and assay
- Manual multi-channel pipette
- Incucyte® Spheroid Analysis Software Module, version 2018A or newer (Cat. No. 9600-0019)

Optional Materials

- Incucyte® Cool Accessories (Cat. No. 4444)
 - CoolBox 96F System (Includes x2 Block with gelpack and CoolSink 96F)
- Incucyte® Cytotox Red or Green Dye (Cat. No. 4632 or 4633)
- Incucyte® Annexin V Red or Green Dye (Cat. No. 4641 or 4642)
- Incucyte® Caspase 3/7 Green Dye (Cat. No. 4440)
- Incucyte® Nuclight Green, Red, Orange or NIR Lentivirus (Cat. No. 4475, 4476, 4771 or 4805)
- Incucyte® Cytolight Red or Green Lentivirus (Cat. No. 4481 or 4482)

Note: Transfect cells with Nuclight or Cytolight Lentivirus prior to performing multi-spheroid experiments by following the protocols supplied with the reagents.

General Guidelines

- All materials (e.g., culture-ware, reagents) that will come in contact with Matrigel® must be kept cold (on ice, stored at +4° C).
- Follow manufactures guidelines for thawing and storing of 100% Matrigel®. Thaw Corning® Matrigel® overnight by submerging the vial in ice cold water placed in the rear of a refrigerator (+4° C). Do not allow Matrigel® to warm to room temperature at any time as this will induce polymerization.
- We recommend sourcing a batch of Matrigel® with a concentration of ≥ 8 mg/mL.
- Following Matrigel® coating, cell seeding and after treatment addition, 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.

Protocol

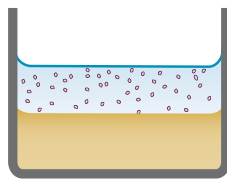
Quick Guide

1. Coat plate (Day 0)



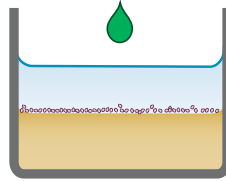
Coat plate (50% Matrigel®, 40 μ L/well). Polymerize at 37° C for 30 minutes.

2. Add cells (Day 0)



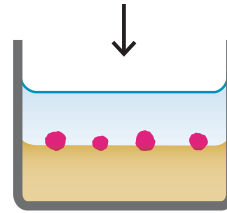
Add cells in media (100 or 150 μ L/well) with or without cell health reagent, respectively.

3. Add reagent (Day 0, optional)



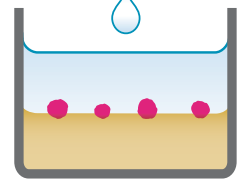
Add cell health reagent (50 μ L/well) at 3X final assay concentration.

4. Monitor formation (Day 0-3)



Place inside the Incucyte® Live-Cell Analysis System and scan every six hours to monitor multi-spheroid formation.

4. Add treatments (Day 3)



Add treatments and continue to monitor growth in Incucyte® Live-Cell Analysis System.

Important:

1. In advance of multi-spheroid experiments, it is important to have stored the Cool Pack accessories at the correct temperatures for at least 4 hours:
 - a. CoolBox x1 (block with gelpack: -20° C),
 - b. CoolSink 96F x1 (4° C)
2. Keep all culture-ware and reagents coming in contact with Matrigel® on ice during the entire process.
3. Store pipette tips used for dispensing diluted Matrigel® at +4° C.

Day 0

Coat Plate with Matrigel®

- 1.1 In a cell culture hood, chill plates (10–15 minutes) on a pre-chilled CoolSink 96F within a CoolBox 96F box.
- 1.2 In a cold polypropylene tube, dilute 100% Matrigel® 1:1 in cold serum-free culture media (keep all Matrigel® solutions on ice).

Note: To prevent incomplete gel formation, for coating we recommend using ≥ 4 mg/mL Matrigel®. As a guideline, a total volume of 5 mL diluted Matrigel® will adequately coat a single 96-well plate.
2. a. To coat a single 96-well plate, add 2.5 mL of cold serum-free culture media to a pre-chilled polypropylene tube.
2. b. Using a cold serological pipette, slowly pipette 2.5 mL of 100% Matrigel® into the serum-free media and, taking care to avoid bubbles, slowly mix by pipetting the solution up and down.
- 1.3 Pour prepared solution into a pre-chilled sterile reagent reservoir (keep on ice).

- 1.4 Using pre-chilled pipette tips and reverse pipetting technique, coat the pre-chilled 96-well plate by carefully adding 40 μL of diluted Matrigel[®] into the center of each well.
 - a. While the plate is cold and Matrigel[®] is still liquid, gently rock the plate once within the CoolBox to ensure even coating of each well.

Note: To avoid cell penetration to the base of the plate, coat wells with a minimum of 40 μL . Use of reverse pipetting technique is important to minimize bubbles.
- 1.5 Remove any bubbles using a wash bottle containing 70–100% ethanol, with the inner straw removed, to blow vapor over the surface of each well.
- 1.6 Place the plate in a 37° C incubator for 30 minutes to polymerize the Matrigel[®].

Seed Cells

- 2.1 Seed cells of interest (100 μL per well if using cell health reagent, 150 μL if not using cell health reagent) at an appropriate density on top of polymerized Matrigel[®] base such that by day 3, multi-spheroids have formed with the desired size (e.g., 30–80 μm in diameter).

Note: Seeding density will need to be optimized for each cell type used. As an example and guide, we recommend seeding A549, MCF-7 and MDA-MB-231 at 1000–2000 cells per well or SKOV-3 at 2000–4000 cells per well.

Add Cell Health Reagent

Note: Annexin V Dye requires solubilization in assay media before use. Centrifuge briefly to collect solid in bottom of vial and add 100 μL assay media to achieve a 100% stock concentration.

- 3.1. Prepare Cell Health Reagent at 3X required concentration. This concentration may require optimization for specific cell lines however as a guide we recommend the use of Cytotox Dyes at 250 nM final assay concentration, Annexin V Dyes at 1% final assay concentration, and Caspase 3/7 Dye (green only) at 2.5 μM final assay concentration.

- a. Incucyte[®] Cytotox Dye: Dilute the stock solution 1:1333 in complete medium to make a 750 nM (3X final assay concentration) working solution.
 - b. Incucyte[®] Annexin V Dye: Dilute the stock solution 1:33 in complete medium to make a 3% (3X final assay concentration) working solution.
 - c. Incucyte[®] Caspase 3/7 Green Dye: Dilute the stock solution 1:666.7 in complete medium to make a 7.5 μM (3X final assay concentration) working solution.
- 3.2. Add the Cell Health Reagent solution(s) on top of the cells (50 μL total per well).
 - 3.3. Place plate in a 37° C incubator for 30 minutes prior to scanning.

Day 0–3

Monitor Multi-Spheroid Formation

- 4.1. Place the cell plate into the Incucyte[®] Live-Cell Analysis System and schedule 24 hour repeat scanning:
 - a. Objective: 10X (96-well corning) 1 image per well
 - b. Channel selection; Phase Contrast + Brightfield and Fluorescence depending on reagent used
 - c. Scan type: Spheroid, Spheroid Type: Multi
 - d. Scan interval: Every 6 hours

Day 3

Add Treatments

- 5.1. Once multi-spheroids have reached desired size, remove the plate from the Incucyte[®] Live-Cell Analysis System and carefully add appropriate treatments at:
 - a. No cell health reagent utilized: 4X final assay concentration (50 μL per well)
 - b. With cell health reagent utilized: 11X final assay concentration (15 μL per well).
- 5.2. Continue to monitor multi-spheroid growth (e.g., every 6 hours for 7 days).

Re-feed cultures (optional and not recommended when using Annexin, Cytotox, or Caspase 3/7 Dyes).

Analysis Guidelines

- Result: Size, number and viability | health measurements
- Suggested Metric: Brightfield Object Area (Total, Average), Object Count
- Secondary Metrics: Fluorescent metrics within a Brightfield Object Boundary
 - For cell health reagents, use Mean Intensity within Brightfield Object Boundary.
 - For fluorescently labelled cell lines, use Integrated Intensity within Brightfield Object Boundary.
- Spectral Unmixing: To analyze green reagent response in Nuclight Red-labeled cells, remove 12% red from the green channel. To analyze red reagent response in Nuclight or Cytolight Green-labeled cells, there is no need for spectral unmixing.

Find more information at www.sartorius.com/incucyte

Sales and Service Contacts

For further contacts, visit
www.sartorius.com

Essen BioScience, A Sartorius Company

www.sartorius.com/incucyte

E-Mail: AskAScientist@sartorius.com

Specifications subject to change without notice.

© 2020. All rights reserved. Incucyte, Essen BioScience, and all names of Essen BioScience products are registered trademarks and the property of Essen BioScience unless otherwise specified. Essen BioScience is a Sartorius Company. Publication No.: 8000-0592-C00

Status: 09 | 2020

North America

Essen BioScience Inc.
300 West Morgan Road
Ann Arbor, Michigan, 48108
USA
Telephone +1 734 769 1600
E-Mail: orders.US07@sartorius.com

Europe

Essen BioScience Ltd.
Units 2 & 3 The Quadrant
Newark Close
Royston Hertfordshire
SG8 5HL
United Kingdom
Telephone +44 1763 227400
E-Mail:
euorders.UK03@sartorius.com

APAC

Essen BioScience K.K.
4th Floor Daiwa Shinagawa North
Bldg.
1-8-11 Kita-Shinagawa
Shinagawa-ku, Tokyo
140-0001
Japan
Telephone: +81 3 6478 5202
E-Mail: orders.US07@sartorius.com