

IncuCyte® S3 Multi-Spheroid Assay

This protocol describes a solution for creating multi-spheroids using 96-well flat bottom plates coated with Matrigel®. 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 #3595)
- Matrigel®, protein concentration ≥ 8 mg/mL (Corning Cat #356234)
- Wet ice
- Serum-free cell culture media for Matrigel® dilutions
- Complete culture media for cell culture and assay
- Manual multi-channel pipette
- IncuCyte® S3 Spheroid Software Module, version 2018A (Essen Cat #9600-0019)

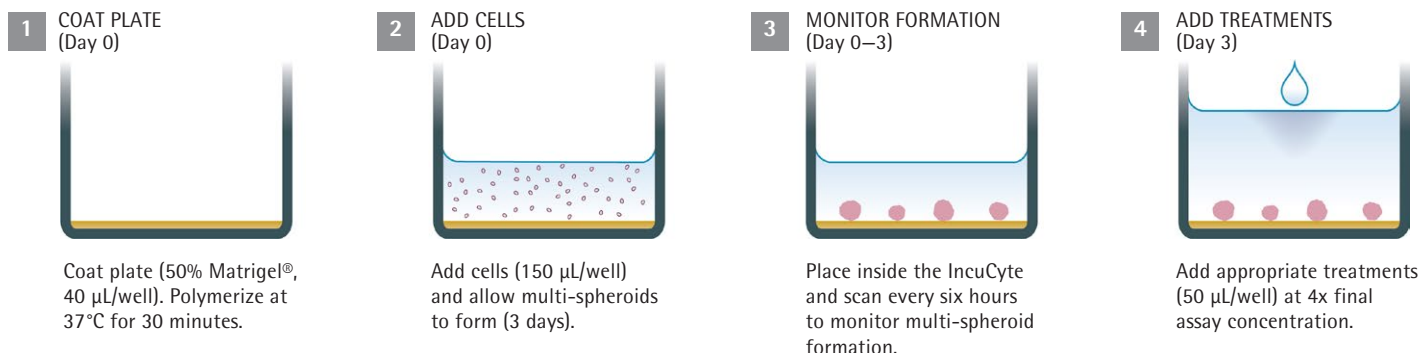
Recommended materials

- IncuCyte® Cool Accessories (Cat # 1500-0078)
 - CoolBox 96F System (Includes x1 Block with gelpack and Coolsink 96F)
- IncuCyte® NuLight Red or Green Lentivirus Reagent (Essen Bioscience, Cat #4476 or #4475)
- IncuCyte® CytoLight Red or Green Lentivirus Reagent (Essen Bioscience, Cat #4481 or #4482)

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 manufacture's guidelines for thawing and storing of 100% Matrigel®. Thaw Corning® Matrigel® overnight by submerging the vial in ice cold water 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.

Quick Guide



Protocol

IMPORTANT:

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

Day 0

1 Coat plate with Matrigel®

- In a cell culture hood, chill plates (10 – 15 minutes) on a pre-chilled Coolsink 96F within a Coolbox 96F box.
- 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®.

 - To coat a single 96-well plate, add 2.5 mL of cold serum-free culture media to a pre-chilled polypropylene tube.
 - 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.
- Pour prepared solution into a pre-chilled sterile reagent reservoir (keep on ice).
- 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.
 - 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®.

2 Seed cells

2.1 Seed cells of interest (150 µL per well) 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 a 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.

2.2 Place plate in a 37°C incubator for 30 minutes prior to scanning.

Day 0–3

3 Monitor multi-spheroid formation

- Place the cell plate into the IncuCyte® live-cell analysis System and schedule 24 hour repeat scanning:
 - Objective: 10x (96-well corning) 1 image per well
 - Channel selection: Phase Contrast + Brightfield and “Red” or “Green” if required
 - Scan type: Spheroid, Spheroid Type: Multi
 - Scan interval: Every 6 hours

Day 3

4 Add treatments

- 4.1 Once multi-spheroids have reached desired size, remove the plate from the IncuCyte and carefully add appropriate treatments at 4x final assay concentration (50 μ L per well).
- 4.2. Continue to monitor multi-spheroid growth (e.g. every 6 hours for 7 days).

5 Re-feed cultures (optional)

- 5.1 Re-feed cultures every 3 – 4 days. Remove plate from IncuCyte. Carefully remove 100 μ L of media per well and replace with 100 μ L of media containing test agents (1x final assay concentration).
- 5.2 Return plate to the IncuCyte and continue to monitor multi-spheroid growth.

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