

Detailed T Cell Incucyte[®] Transendothelial Migration Chemotaxis Protocol

Materials

- Freshly isolated blood
- Lympholyte[®]-H Cell Separation Media (Cedarlane CL5010)
- Dynabeads[®] Human T-activator CD3/CD28 (Life Technologies 11131D)
- RPMI 1640 Medium (Life Technologies 11875-085)
- Dialyzed Fetal Bovine Serum (FBS; Omega Scientific FB-03)
- HUVEC Cells (LONZA C2517A)
- EGM-2 MV Bullet Kit (LONZA CC-3202)
- Collagen Type I Rat Tail (BD Biosciences 354236)
- Fibronectin (Sigma Aldrich F1141)
- Acetic Acid (0.02N) – used in Collagen coating step
- D-PBS -/- (w/o Ca²⁺, Mg²⁺, Life Technologies 10010) – used in Fibronectin coating step
- D-PBS +/- (with Ca²⁺, Mg²⁺, Life Technologies 14040) – used in monolayer wash step
- SDF-1a (R&D systems 350-NS-050)
- Incucyte[®] Clearview 96-well Plate (Sartorius Cat. No. 4582, 4599, or 4648)

Create HUVEC Monolayer (Day 1)

1. Prepare extracellular matrix coating of either 50 µg/mL Collagen diluted with 0.02N Acetic Acid or 5 µg/mL Fibronectin diluted with D-PBS (-/-).
2. Using reverse pipetting, add 150 µL of coating solution into the reservoir of an Incucyte[®] Clearview 96-well Plate. Gently place the insert into the reservoir and reverse pipette 20 µL of the fibronectin or collagen solution into the insert an Incucyte[®] Clearview Plate.
3. Incubate for 1 hour at ambient temperature.
4. During incubation, harvest and count HUVEC cells and prepare a cell seeding stock of 100,000 cells/mL in full growth medium (EGM-2).
5. Aspirate the coating from the reservoir plate and replace with 200 µL of D-PBS (-/-) and gently return the insert into the reservoir plate.
6. To the insert, directly add 60 µL D-PBS (-/-) to the inserts containing coating, then aspirate the entire volume.
7. Immediately seed 60 µL of the HUVEC seeding stock using a multi-channel pipette and reverse pipetting technique into every well of the insert plate (60 µL per well, 6,000 cells per well).

Calculation: 100,000 cells/mL x 0.06 mL = 6,000 cells per insert well.

8. Allow the HUVECs to settle at ambient temperature on a level surface for 15 minutes.
9. Place the Incucyte® Clearview Plate containing HUVEC cells at 37 °C and incubate for 24 hours.

Initiating TEM (Day 2)

1. After HUVEC monolayer has formed, gently wash the monolayer 2X with D-PBS (+/+), using partial washes
Note: It is important not to disrupt the HUVEC monolayer. It is recommended to gently remove about half of the growth medium then add 60 µL D-PBS for both washes. At the final wash step, remove as much of the medium/D-PBS as possible without disrupting the monolayer.
2. Prepare T cell seeding stock at 83,333 cells/mL in defined EBM-2 (EBM media + 2% FBS + hydrocortisone + ascorbic acid + heparin).
3. Using a manual multi-channel pipette and reverse pipetting technique, seed cells (60 µL per well, 5,000 cells per well) into every well of the insert plate of an Incucyte® Clearview Plate, being careful not to disrupt the HUVEC monolayer.
Calculation: 83,333 cells/mL x 0.06 mL = 5,000 cells per insert well.
4. Centrifuge the Incucyte® Clearview Plate for 3 minutes at 50 x g in order to quickly bring the T cells to the monolayer surface. Alternatively, if centrifugation is not possible, allow the T cells to settle on the endothelial monolayer at ambient temperature for 45–60 minutes.

5. Using a manual multi-channel pipette, add 200 µL of the chemoattractant and control medium to the appropriate wells of the second Incucyte® Clearview Reservoir Plate.
6. Carefully transfer the insert plate containing the cells into the pre-filled second reservoir plate containing assay medium ± chemoattractant.
7. Place the Incucyte® Clearview Plate into the Incucyte® Live-Cell Analysis System and allow the plate to warm to 37 °C for at least 15 minutes.

Note: After 15 minutes, wipe away any condensation that remains on the outside of the plate lid or bottom of the reservoir.

8. In the Incucyte® software, schedule 24 hour repeat scanning (10X) for every 30 minutes.

Note: This schedule is only for a scanning a single plate. Fewer scans will be required if scheduling multiple plates.

a. Objective: Ensure 10X objective is installed

b. Vessel Type: Select "Incucyte® Clearview Plate"

c. Channel Selection: Select "Phase"

d. Scan Mode: Select "Chemotaxis (Top/Bot)" scan type and desired Scan Pattern

Note: The Incucyte® Live-Cell Analysis System estimates a scan time of 20 minutes per plate (phase only); however, the actual scan time can take longer.

North America

Sartorius Corporation
300 West Morgan Road
Ann Arbor, Michigan 48108
USA
Phone +1 734 769 1600

Europe

Sartorius UK Ltd.
Longmead Business Centre
Blenheim Road
Epsom
Surrey, KT19 9QQ
United Kingdom
Phone +44 1763 227400

Asia Pacific

Sartorius Japan K.K.
4th Floor, Daiwa Shinagawa North Bldg.
1-8-11, Kita-Shinagawa 1-chome
Shinagawa-Ku
Tokyo 140-0001
Japan
Phone +81 3 6478 5202