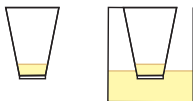


IncuCyte™ Chemotaxis Cell Migration Quick Start Guide

1 COAT INSERT
(optional)**Coat Membrane Surface**

Coat membrane with extracellular matrix.

2 HARVEST &
SEED CELLS**Cell Seeding**

Seed cells (1,000 or 5,000/well) into IncuCyte ClearView Plate.

3 TREAT CELLS
(optional)**Compound Addition (optional)**

Prepare 3x concentrations of treatment and add to wells. Mix by trituration.

4 ALLOW CELLS
TO SETTLE**Cell Settling**

Place plate on level surface and allow cells to settle.

5 ADD CHEMO-
ATTRACTANT**Reservoir Addition**

Add chemoattractant or controls to the reservoir plate. Place the insert into reservoir.