

StemKit Protocol

Shipped from Essen BioScience Room Temperature Shipper Dry Ice Shipper (Media/Plate) (Cells/Media Supplements) Assay **ECFC** Growth Media Growth Supplements CytoLight Green

At Customer Site

Upon arrival, store cells in liquid nitrogen, media at 4 °C and media supplements at -20 °C.

Day -2

Thaw ADSC and seed into T-225 flask

Day -1

- Thaw ECFC CytoLight Green and seed into Collagen coated T-75
- Harvest ADSCs and seed into 96-well plate

Day 0 (Assay initiation)

- Harvest ECFCs and seed over ADSC monolayer
- Add growth factor ± treatment agents
- Image in IncuCyte[™] FLR or ZOOM every 2-6 hours for 4-5 days
- Kinetic analysis using integrated IncuCyte™ algorithms
 - Tube Length, Tube Area, Branch Points



Day 4

Terminate experiment or investigate disruption of established networks

