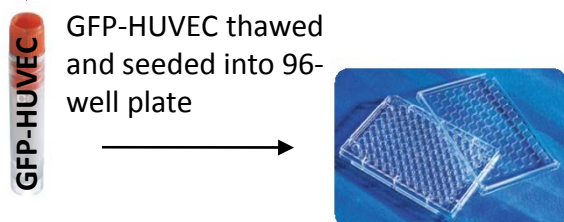
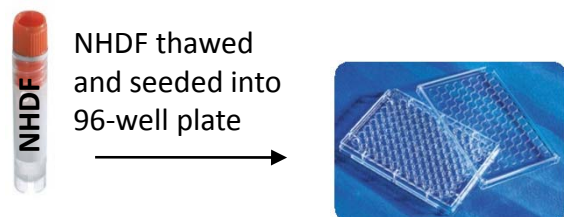


LiveKit Protocol

At Essen BioScience

Day 0 (Friday) – Preparation



Day 3 (Monday) – Shipping

Live-cell plate and media shipped to customer



Day 4-14 – Kit Validation

Partial plate including VEGF and suramin response performed at Essen BioScience using cells from same vial

At Customer Site

Day 4 (Tues) – Upon arrival unpack plate, feed with growth factor ± test agent, load plate in IncuCyte, store assay media at 4 °C



Day 5, 7, 10, 12 – Assay media/test agent replacement, monitor progress, process tube formation dynamically using integrated IncuCyte algorithms

Day 14 – Terminate Experiment

