



# CryoKit Protocol


### Shipped from Essen BioScience

<p><b>Room Temperature Shipper (Media/Plates)</b></p> 	<p><b>Dry Ice Shipper (Cells/Media Supplements)</b></p> 
-----------------------------------------------------------------------------------------------------------------------------------------	-------------------------------------------------------------------------------------------------------------------------------------------

### At Customer Site

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

**Day 0**



NHDF thawed and seeded into 96-well plate

