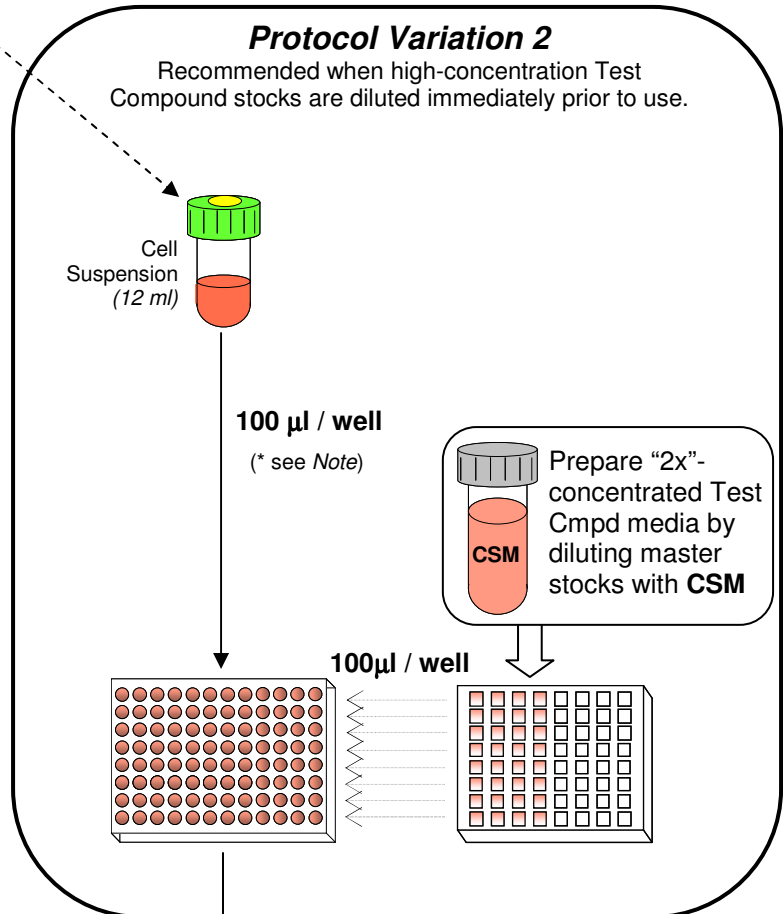
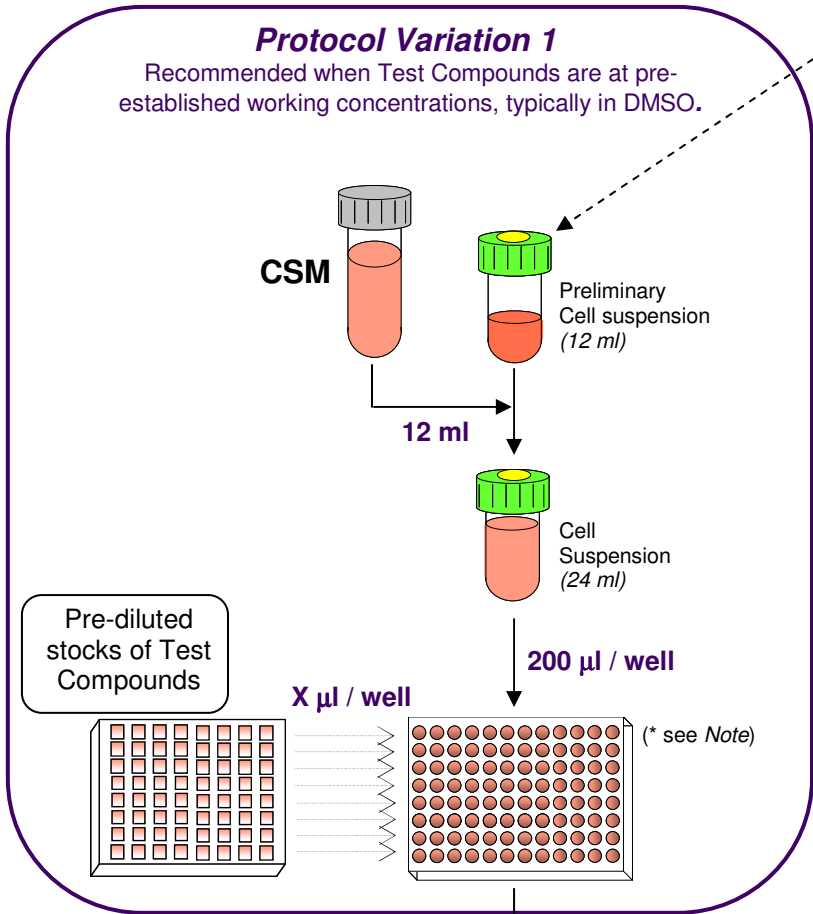
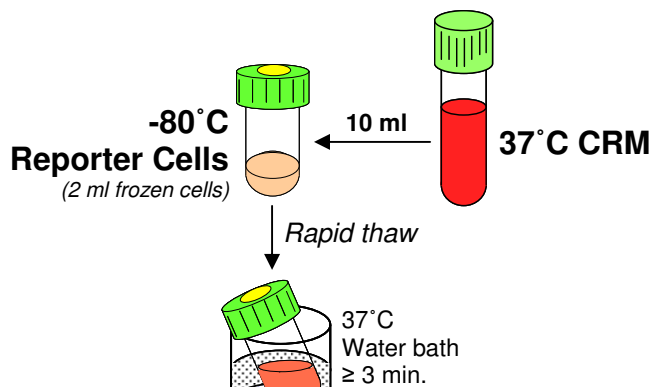


Nuclear Receptor Reporter Assay System
96-well Format Assays
~ Protocol Quick Guide ~

DAY 1

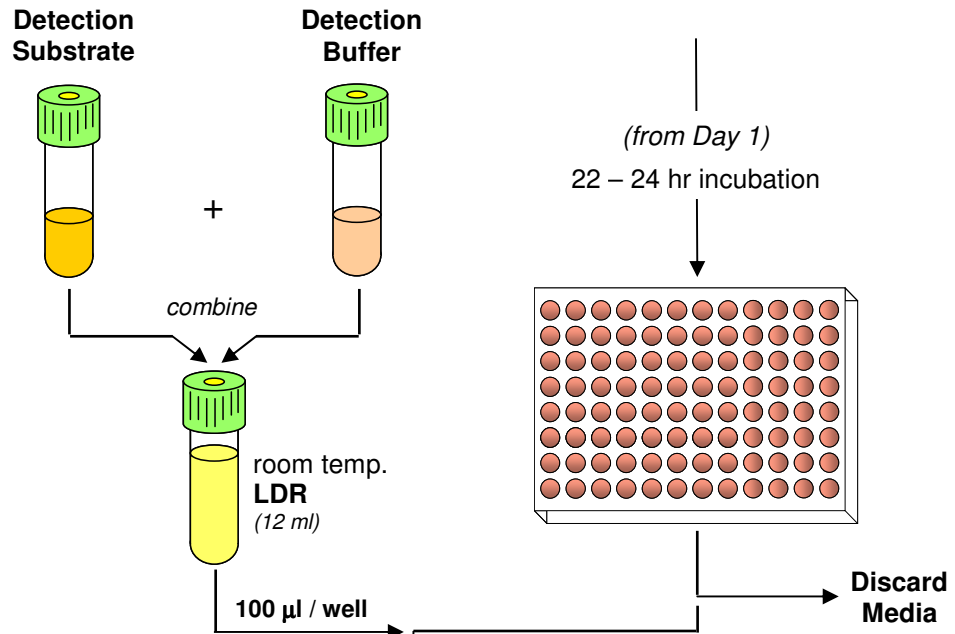
* *Note:* For microscopic examination of cells in assay wells, dispense the excess volume of cell suspension into a clear 96-well plate. Treat +/- test compounds, as desired.



Incubate 22 – 24 hrs
 37°C / humidified 5% CO₂

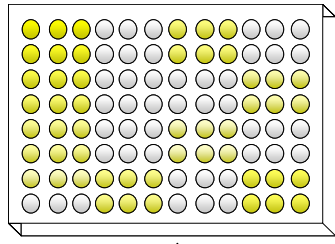


DAY 2



15 minute rest at room temp.

Quantify Luminescence
 Read time ≤ 0.5 sec / well



Data Analyses

