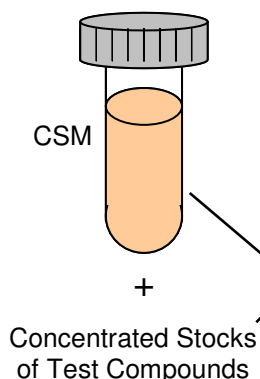




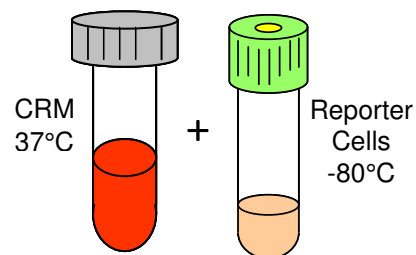
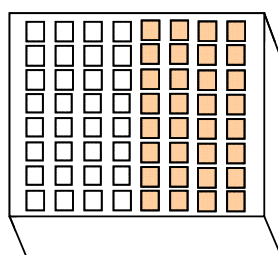
DAY 1

Step 1: Thaw CRM ⇒ 37°C
Thaw CSM ⇒ room temperature

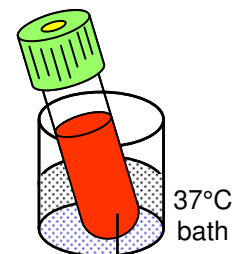


**Step 2: Use CSM to prepare
 “2x Test Cmpd” media**

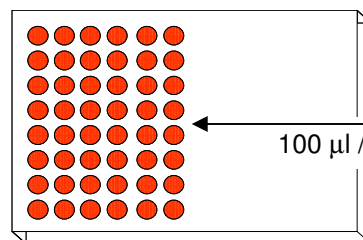
dilutions



**Steps 3 & 4: Rapid thaw of
 Reporter Cells**
 ⇒ -80°C Cells + 37°C CRM
 ⇒ 37°C ≥ 3 min.

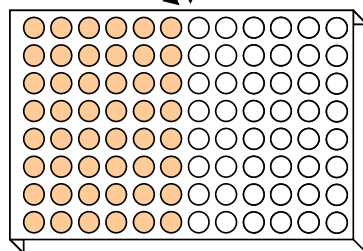


Step 5: Dispense Reporter Cells
 ⇒ 100 µl per well



**Step 6: Dispense “2x-Test Cmpd”
 media ⇒ 100 µl per well**

100 µl / well



Step 7: Incubate assay plate
 22 – 24 hr

**Step 8: Transfer “Detection
 Solutions I & II” to 4°C
 for overnight thaw**

Incubate
 37°C, humidified 5% CO₂

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



INDIGO Biosciences, Inc.
The Nuclear Receptor Company™

DAY 2

Step 9: Equilibrate “Detection Solutions I & II” to room temperature

Step 10: Prepare luminometer

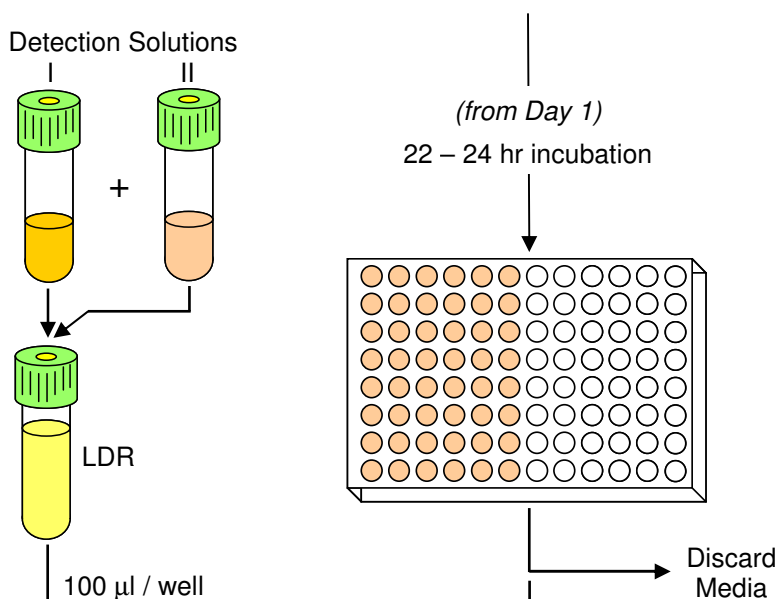
Step 11: Combine “Detection Solutions I & II” to generate “LDR”

Step 12: Discard media from assay plate

Step 13: Dispense LDR
⇒ 100 µl per well

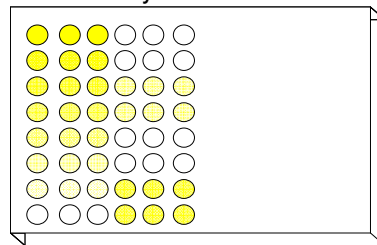
Step 14: Sit 15 minutes, room temperature

Steps 15: Read assay plate, 0.5 sec / well



15 minute rest

Quantify Luminescence



Data Analyses

