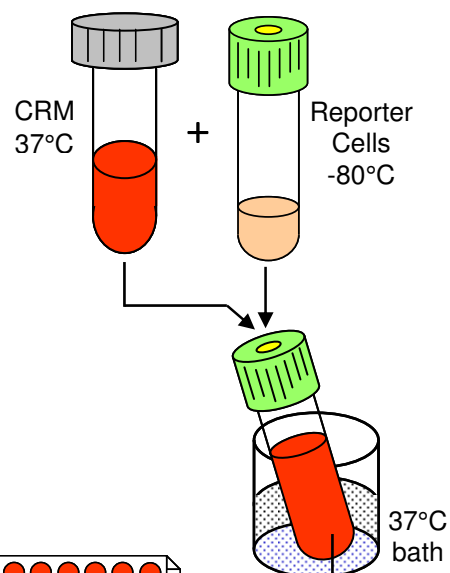
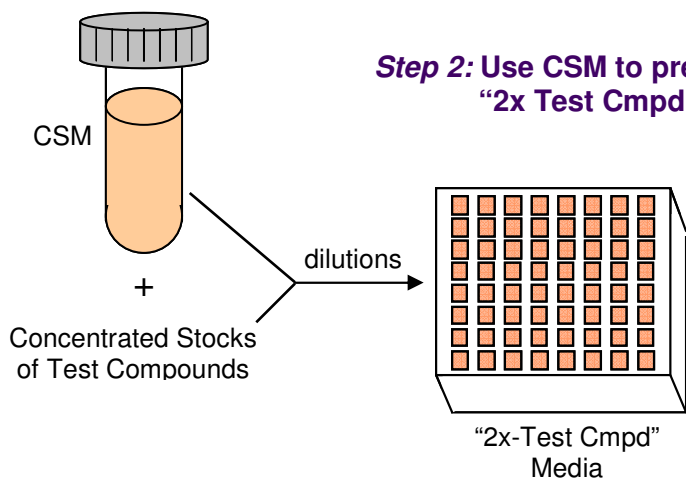




DAY 1

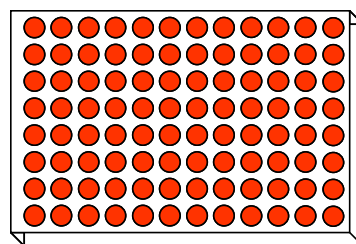
Step 1: Thaw CRM ⇒ 37°C
Thaw CSM ⇒ room temperature

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

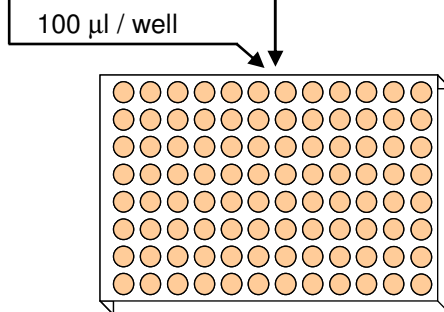


Steps 3 & 4: Rapid thaw of Reporter Cells
 ⇒ -80°C Cells + 37°C CRM
 ⇒ 37°C ≥ 3 min.

Step 5: Dispense Reporter Cells
 ⇒ 100 µl per well



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



Step 7: Incubate assay plate
 22 – 24 hr

Step 8: Transfer “Detection Substrate” & “Detection Buffer” to 4°C for overnight thaw

Incubate
 (37°C, ≥ 90% humidity, 5% CO₂)

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



INDIGO Biosciences, Inc.
 The Nuclear Receptor Company™

DAY 2

Step 9: Equilibrate "Detection Substrate" & "Detection Buffer" to room temp.

Step 10: Prepare Luminometer

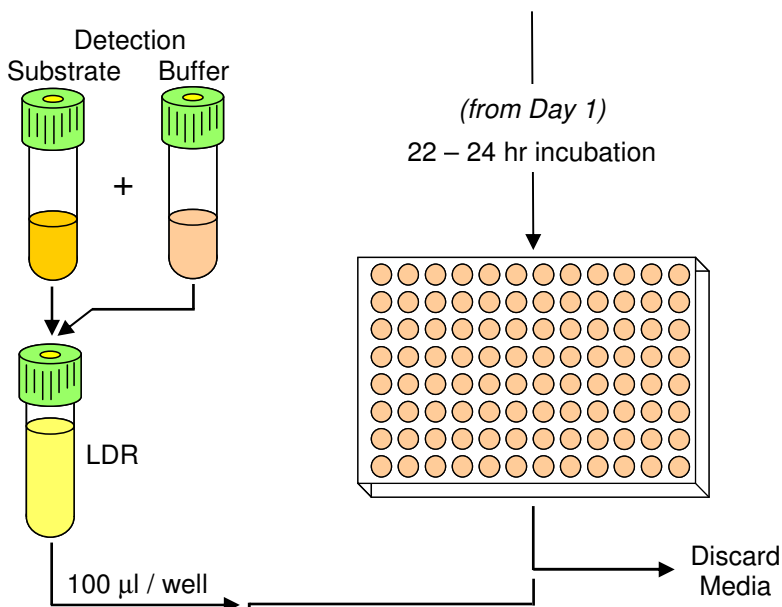
Step 11: Combine Detection reagents to generate "LDR"

Step 12: Discard media from assay plate

Step 13: Dispense LDR
 ⇒ 100 µl per well

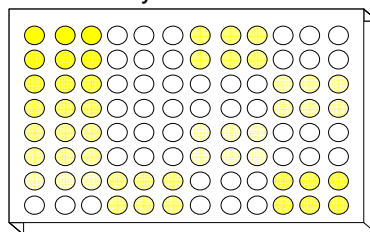
Step 14: Sit 15 minutes, room temperature

Steps 15: Read assay plate, 0.5 sec / well



15 minute rest

Quantify Luminescence



Data Analyses

