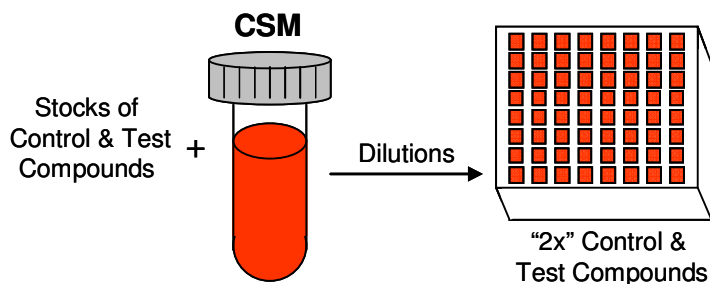


Nuclear Receptor Reporter Assay System
 384-well Format Assays
 ~ Protocol Quick Guide ~



INDIGO Biosciences, Inc.
 The Nuclear Receptor Company™

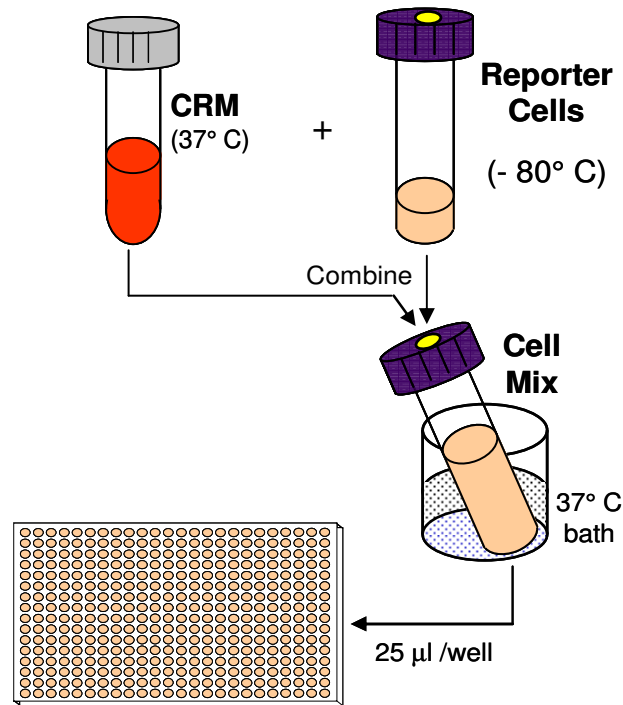
DAY 1



Step 1: Thaw CSM (room temp.) and CRM (37°C)

Step 2: Use CSM to Prepare Compound Dilutions

Steps 3 & 4: Rapid Thaw of Reporter Cells (-80°C Cells + 37°C CRM) ⇒ 37°C, ≥ 3 min.
 Mix Cell Suspension

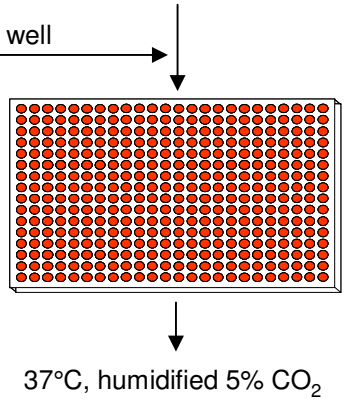


Step 5: Dispense Reporter Cells ⇒ 25 µl per well

Step 6: Dispense "2x-Compound" Media ⇒ 25 µl per well

Step 7: Incubate Assay Plate, 22-24 h

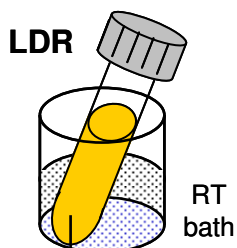
Step 8: Transfer LDR from -80°C to +4°C, overnight





DAY 2

Step 9: Equilibrate LDR to room Temp. ⇒ Mix



Step 10: Discard CSM

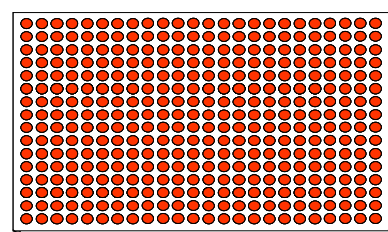
Step 11: Dispense LDR ⇒ 25 µl per well

Step 12: Shake Assay Plate; 5-10 min.

Steps 13 & 14: Prepare Luminometer ⇒ Read Assay Plate

(from DAY 1)

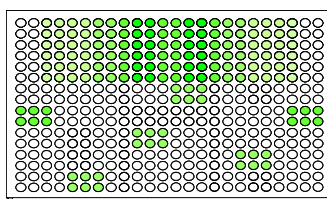
22 - 24 h
37°C, 5% CO₂



Discard Media

25 µl / well

5 - 10 min., r.t.
(Shake)



Data Analyses

