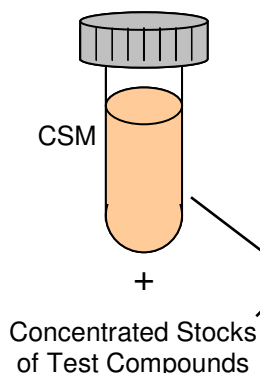




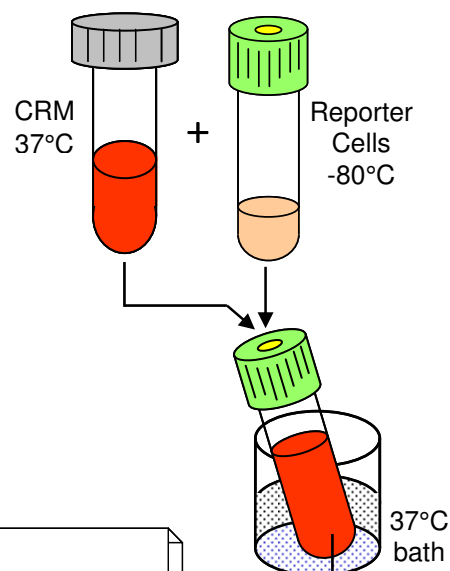
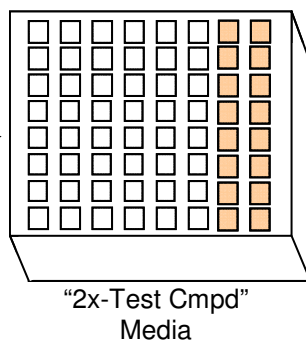
DAY 1

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

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



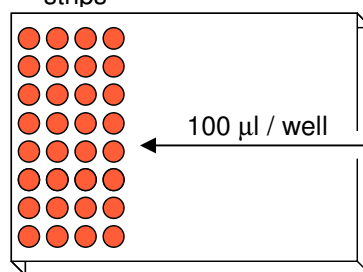
dilutions



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

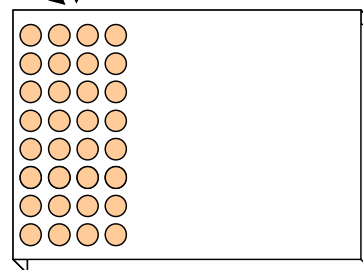
Step 5: Dispense Reporter Cells
 ⇒ 100 µl per well

Four 8-well strips



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



DAY 2

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

Step 10: Prepare Luminometer

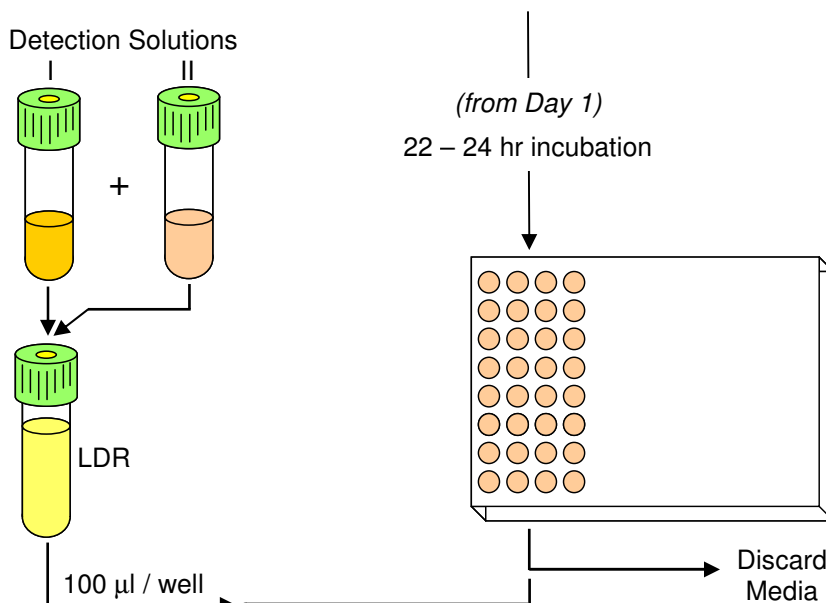
Step 11: Combine “Detection Solutions I & II” to generate “LDR”

Step 12: Discard media from assay plate

Step 13: Dispense LDR
 ⇒ 100 µl per well

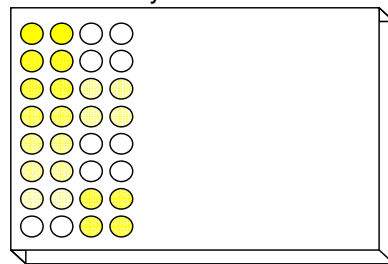
Step 14: Sit 15 minutes, room temperature

Steps 15: Read assay plate, 0.5 sec / well



15 minute rest

Quantify Luminescence



Data Analyses

