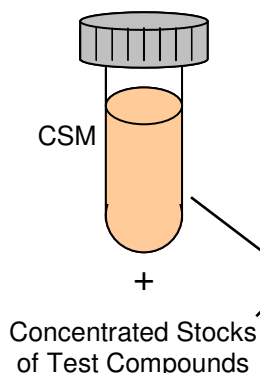




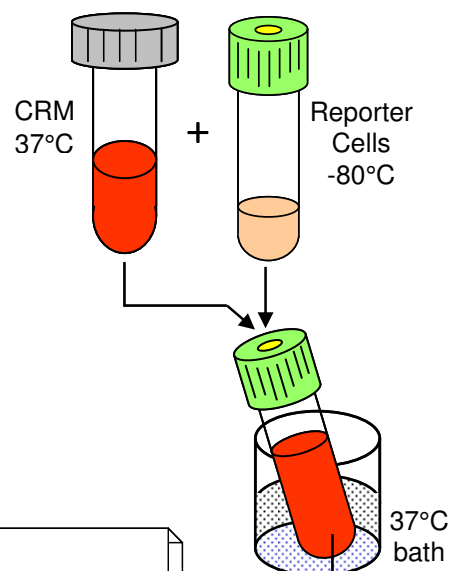
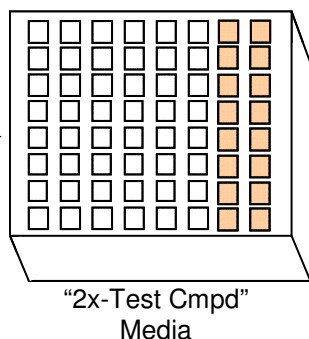
**DAY 1**

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

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

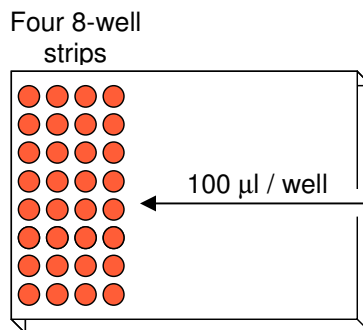


dilutions



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

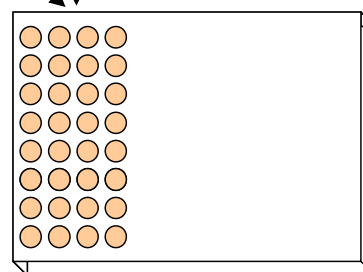
**Step 5: Dispense Reporter Cells**  
 ⇒ 100 µl per well



**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 Substrate” & “Detection Buffer” to 4°C for overnight thaw**

Incubate

(37°C, ≥ 90% humidity, 5% CO<sub>2</sub>)



**DAY 2**

**Step 9:** Equilibrate “Detection Substrate” & “Detection Buffer” to room temp.

**Step 10:** Prepare Luminometer

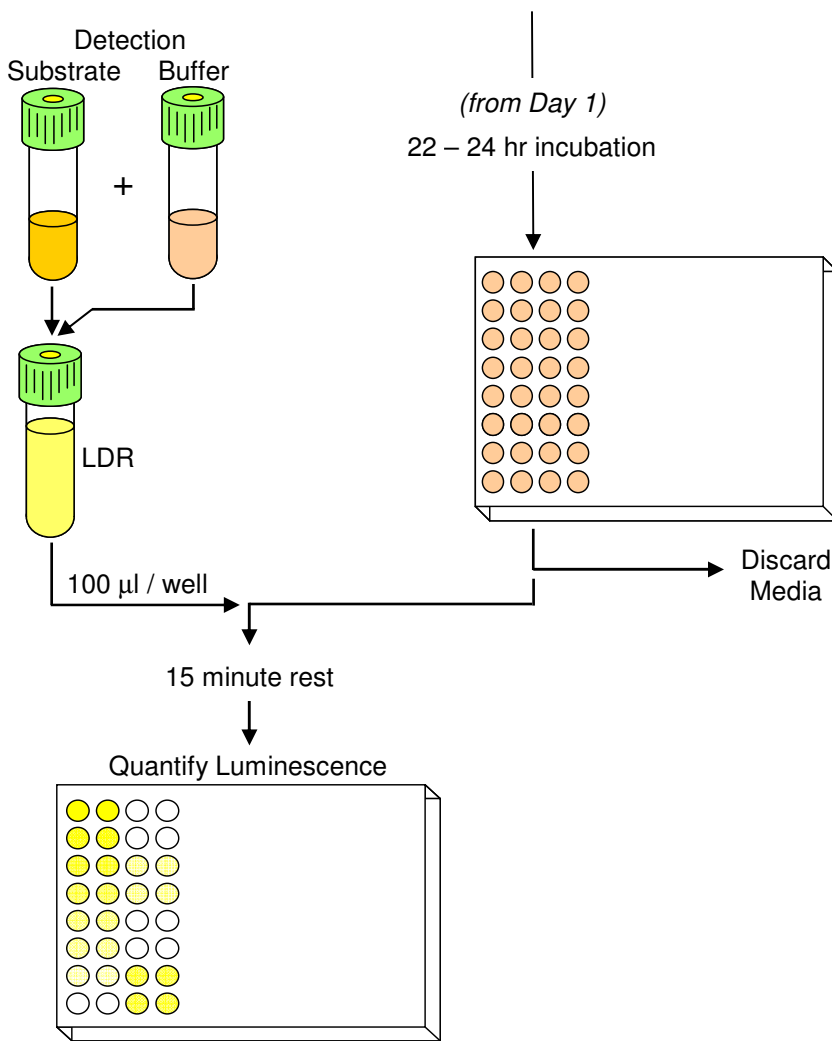
**Step 11:** Combine Detection reagents to generate “LDR”

**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



Data Analyses

