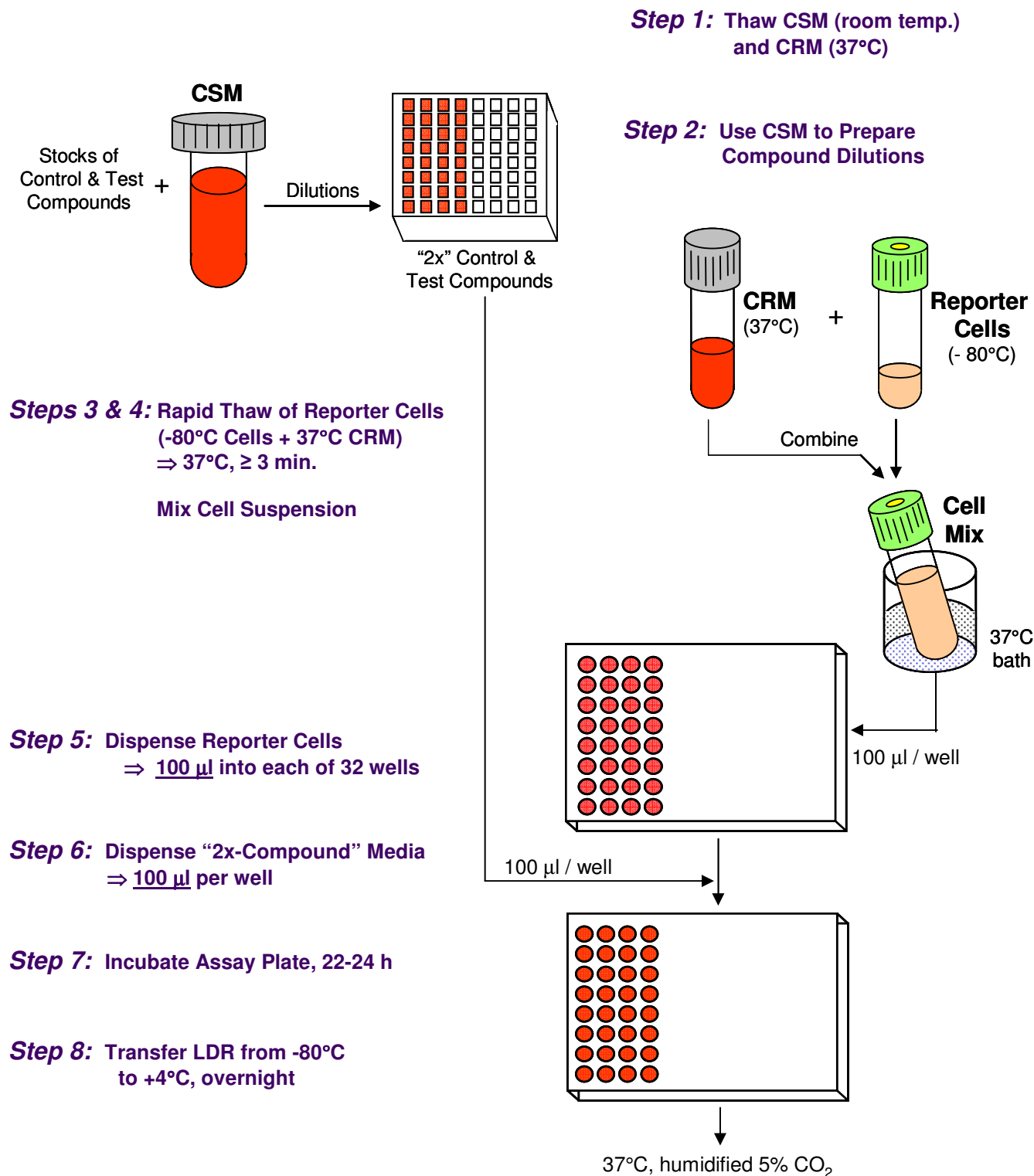




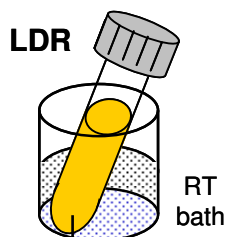
DAY 1





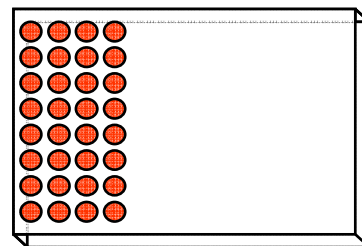
DAY 2

Step 9: Equilibrate LDR to room temp. ⇒ Mix



(from DAY 1)

22 - 24 h
 37°C, 5% CO₂



Step 10: Discard CSM

Step 11: Dispense LDR
 ⇒ 100 µl per well

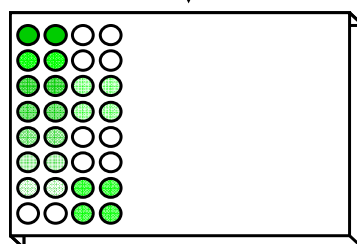
100 µl / well

Discard Media

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

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

Steps 13 & 14: Prepare Luminometer
 ⇒ Read Assay Plate



Data Analyses

