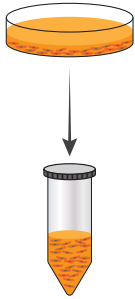
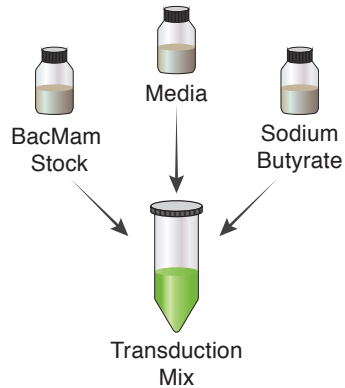


Day 1: Transduce and Plate Cells

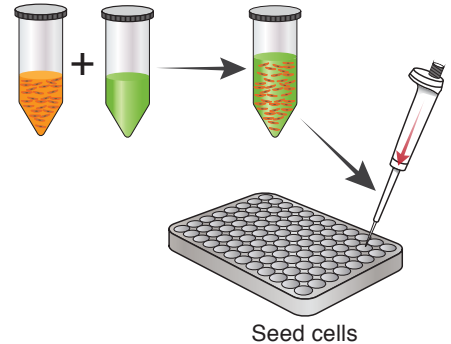
Step 1 Prepare Cells



Step 2 Prepare Viral Transduction Reaction

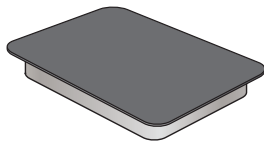
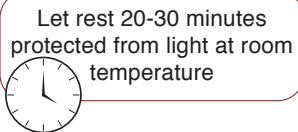
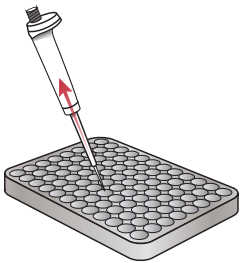


Step 3 Mix Cells and Transduction Mix



Day 2: Live Cell Assay

Step 1 Replace culture media with DPBS



Step 2 Measure Fluorescence

