



Signosis
Innovative Plate Assay Solutions

Firefly Luciferase Lysis Buffer 5x

(Catalog# LS-001)

(For Research Use Only)

Product description

Luciferase Cell Culture lysis buffer 5x is used with the Signosis Firefly Luciferase Substrate (Cat# LUC015 and LUC100), developed for reporter quantitation in mammalian cells. This lysis buffer provides rapid lysis of a variety of cell types, including adherent and nonadherent mammalian cells, plant cells and disuse homogenates.

Material provided

One bottle of 5x Luciferase Cell Culture lysis buffer (10 ml)

Handling upon arrival

Store 5x Luciferase Cell Culture lysis buffer at -20°C

Protocol

1. Add 4 volume of water to 1 volume of Luciferase Cell Culture Lysis buffer 5x. Equilibrate 1x cell culture lysis buffer to room temperature.
2. Carefully remove the growth medium from cells to be assayed. Rinse cells with PBS, being careful to not dislodge attached cells.
3. Add enough 1x cell culture lysis buffer to cover the cells. (eg. 400 μ l/60mm culture dish, 900 μ l/100mm culture dish or 20 μ l per well of 96 well plate).
4. Rock culture dish for 10 minutes to ensure complete coverage the cells with lysis buffer. Scrape attached cells from the dish. Transfer cells and all liquid to a microcentrifuge tube. Place the tube on ice.
Note: If using 96- well plates and a plate-reading luminometer, the cell lysate may be left in the well.
5. Vortex the microcentrifuge tube 10-15 seconds, then centrifuge at 12,000 g for 15 seconds at room temperature.
6. The cell lysates maybe assayed directly or stored at -70°C

Materials required but not provided

- PBS buffer (Mg²⁺ and Ca²⁺ free)