



2x Cell Lysis Buffer for ELISA

Catalog Number EA-0001

(For Research Use Only)

2X Cell Lysis Buffer (10ml)

Prepare 1X Cell Lysis buffer by adding 10ml ddH₂O to 10ml 2X Cell Lysis Buffer.

Cell Lysate from cultured cells on a 96-well plate

1. Remove the cell medium and rinse cells once with ice-cold 1X PBS.
2. Freeze the cells and thaw the cells on ice.
3. Add 100ul of 1x Cell lysis buffer and incubate on ice for 10 minutes with gentle shaking.
4. Centrifuge the sample at 3,000 RPM for 5 minutes.
5. Transfer 90ul of supernatant to a well of ELISA plate.

Cell lysate from tissue

1. Weight tissue sample and add 1ml of 1X Cell lysis buffer to 100mg of tissue.
2. Homogenized tissues with grinder (PowerGen 125 or equivalent) on ice.
3. Sonicate lysates briefly on ice.
4. Centrifuge the sample at 10,000 RPM for 5 minutes to pellet the tissue debris.
5. Collect supernatant and measure the protein concentration of supernatant. The supernatant then can be used for assay. The supernatant also can be aliquoted, and frozen at -80°C. Avoid multiple freezes/thaws.
6. Dilute the cell lysate in the provided Diluent Buffer to the final 10ug/100ul per well.