

.....  
Now Includes Protocol for  
Bio-Plex Phospho-Histone H3  
Lysate Preparation!

## **Bio-Plex™ Cell Lysis Kit Product Insert**

For use with  
Bio-Plex phosphoprotein assays and  
Bio-Plex total target assays

For technical service, call your local Bio-Rad office,  
or in the US, call 1-800-4BIORAD (1-800-424-6723)

4110051 Rev D

For research use only.  
Not for diagnostic procedures.

**BIO-RAD**

## **Introduction**

The Bio-Plex cell lysis kit has been developed specifically to prepare cell culture and tissue lysate samples for analysis with Bio-Plex phosphoprotein and total target assays. The cell lysates can be tested for the presence of phosphorylated proteins using Bio-Plex phosphoprotein assays or for the abundance of target proteins using Bio-Plex total target assays. This cell lysis kit can also be used to prepare cell lysates for western blot analysis (request bulletin 3033).

## **Product Description**

The following components are provided with the Bio-Plex cell lysis kit:

Cell wash buffer

Cell lysis buffer

Cell lysis buffer, factor 1 (250x)

Cell lysis buffer, factor 2 (500x)

## Storage and Stability

The cell wash buffer and cell lysis buffer should be stored at 4°C. Factors 1 and 2 should be stored at -20°C and can be frozen and thawed up to 5 times. All components are guaranteed for 6 months from the date of purchase when stored as specified.

## Materials Required but Not Supplied

Phenylmethylsulfonyl fluoride (PMSF),  
Sigma catalog #P7626

Dimethyl sulfoxide (DMSO),  
Sigma catalog #D2650

Laemmli sample buffer,  
Bio-Rad catalog #161-0737

2-Mercaptoethanol,  
Bio-Rad catalog #161-0710

## Lysate Preparation

### Adherent and Suspension Cell Preparation

1. Rinse the samples with *cell wash buffer* as follows:

**Adherent Cells** — Stop the treatment reaction by aspirating the culture medium and quickly rinsing the cells with ice-cold cell wash buffer. The volume of cell wash buffer required is the same as the volume of aspirated cell culture medium. Keep the cells on ice.

**Suspension Cells** — Stop the treatment reaction by adding ice-cold wash buffer to the cells. The volume of cell wash buffer required is twice that of the culture medium. Centrifuge the cells at 1,000 rpm for 5 min at 4°C. Aspirate the supernatant.

**Tissue Samples** — Rinse the tissue sample with cell wash buffer once. Cut the tissue into 3 x 3 mm pieces and transfer them to a 2 ml tissue grinder.

2. Prepare 500 mM PMSF by dissolving 0.436 g PMSF in 5 ml DMSO. Store as 0.5 ml aliquots at  $-20^{\circ}\text{C}$ . Aliquots can be frozen and thawed up to 5 times.
3. Prepare an adequate volume of lysing solution (refer to the table on the left). For 10 ml of lysing solution, add 40  $\mu\text{l}$  of *factor 1* and 20  $\mu\text{l}$  of *factor 2* to 9.9 ml of *cell lysis buffer*. Vortex gently to mix and set aside on ice. Then add 40  $\mu\text{l}$  of 500 mM PMSF.

#### Lysing Solution Volume Guide

Culture vessel	Culture medium volume	Lysing solution volume	Notes
96-well plate	100 $\mu\text{l}$ /well	75 $\mu\text{l}$ /well	Grow cells to 80–85% confluence  Recommend leaving external wells empty due to edge effect
10 cm culture dish	10 ml	2–3 ml	Grow cells to 80–90% confluence

4. Lyse the samples:

#### Adherent and Suspension Cells

- a) Immediately add the lysing solution to the cells. The amount of lysing solution needed depends on the cell concentration in the culture vessel (see table on the left).
- b) Agitate the cells as follows:

**Culture Plate** — For suspension cells, place the plate on ice and pipet the contents of the wells up and down 5 times. For adherent cells, scrape the cells with a cell scraper. For both, agitate the plate on a microplate shaker at 300 rpm for 20 min at  $4^{\circ}\text{C}$ .

**Other Culture Vessel** — Transfer the cell lysate to a centrifuge tube and rotate for 20 min at  $4^{\circ}\text{C}$ .

HINT: Freeze-thawing the lysate once using dry ice or a  $-20^{\circ}\text{C}$  freezer may increase the extent of the lysis. Alternatively, briefly sonicate (eg., with a Sonifier 450 as follows: Duty cycle = 40, Output = 1, Pulse sonicating = 18 times).

- c) Centrifuge the samples at 4,500 g for 20 min at  $4^{\circ}\text{C}$ .

## **Tissue Samples**

- a) Immediately add 500  $\mu$ l of lysing solution to the tissue grinder and grind the tissue sample on ice using about 20 strokes.
  - b) Transfer the ground tissue to a clean microcentrifuge tube and freeze the sample at  $-70^{\circ}\text{C}$ .
  - c) Thaw the samples, then sonicate on ice as suggested above.
  - d) Centrifuge the samples at 4,500 g for 4 min.
5. Collect the supernatant without disturbing the pellet.

## **Suggested Protocol for Lysate Preparation of Histone H3 Assay:**

1. Follow steps 1-3 in lysate preparation.
4. Lyse the samples:
  - a) Immediately add the lysing solution to the cells. The amount of lysing solution needed depends on the cell concentration in the culture vessel (see table on the left).
  - b) Briefly sonicate (e.g., with a Sonifer 450 as follows: Duty cycle = 40, Output = 1, Pulse sonicating = two 10 minute pulses with a 1 minute break in between).
  - c) Agitate the cells. Transfer the cell lysate to a centrifuge tube and rotate for 20 min at  $4^{\circ}\text{C}$ .
  - d) Centrifuge the samples at 4,500 g for 20 min at  $4^{\circ}\text{C}$ .
5. Collect the supernatant without disturbing the pellet.

### **For Bio-Plex Phosphoprotein Assays and Bio-Plex Total Target Assays**

1. Determine the lysate protein concentration. The protein concentration should be 200–900 µg/ml. It may be necessary to test-lyse your samples with different volumes of lysing solution to obtain the specified protein concentration range.
2. Add an equal volume of *assay buffer* to the lysate.
3. If the lysate is not tested immediately, store at –20°C. The lysate is stable for up to 5 freeze-thaw cycles. For Bio-Plex Histone H3 assay, freeze lysate (overnight) at –20°C and thaw before testing.
4. For further assay instructions refer to the Bio-Plex phospho-protein detection instruction manual.

### **For Western Blot Analysis**

1. Determine the lysate protein concentration. The protein concentration should be 200–900 µg/ml. It may be necessary to test-lyse your samples with different volumes of lysing solution to obtain the specified protein concentration range.
2. If the lysate is not tested immediately, store at –20°C. The lysate is stable for up to 5 freeze-thaw cycles.
3. Prepare fresh sample loading buffer using a 1:20 dilution of 2-mercaptoethanol and Laemmli sample buffer. Alternatively, another sample loading buffer can be used.
4. Dilute 1 part sample with 2 parts sample loading buffer.
5. For further instructions, refer to Bio-Rad's Laemmli sample buffer manual.

## **Safety Considerations**

Eye protection and gloves are recommended while using this product. Consult the MSDS for additional information.