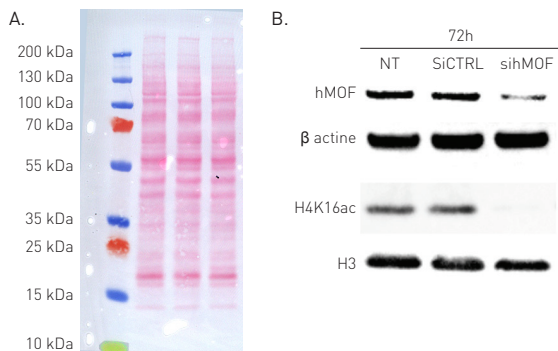


# BIORUPTOR® PROTOCOL

## WESTERN BLOT - PROCEDURE FOR SIMULTANEOUS EXTRACTION OF CYTOPLASMIC, NUCLEAR AND CHROMATIN PROTEINS WITH BIORUPTOR®

The regular protocol for the extraction of histone requires an acid extraction, making impossible the detection of any other cytoplasmic and nuclear proteins from the same extract. Using Bioruptor®, Slimane AIT-SI-ALI and his "epigenetic and cell fate" team from the University Paris Descartes developed a protocol allowing the simultaneous extraction of histone and other proteins.



**Figure:** Protein staining of cell lysates (containing both chromatin and soluble proteins) obtained with the protocol developed by Lauriane Fritsch from the laboratory of Slimane AIT-SI-ALI. A) Ponceau S Staining Western Blot of Hep G2 cells lysates obtained using the Bioruptor®. The rectangle indicates the location of histones. B) Western Blot of Hep G2 cells transfected and cultivated for 72h with a siRNA against hMOF, a histone H4 K16-specific acetyl transferase. Both soluble (hMOF and B-actine) and chromatin (Histone 3 and 4) proteins are obtained on the same cell extract.

### 1. Material required and buffer preparation

- Bioruptor® Plus (Diagenode, Cat. No. B01020001) with 15 ml tube holder and 15 mL sonication probes (Diagenode, Cat. No. B01200013) or 1.5 ml tube holder
- 15mL TPX tubes (Diagenode, Cat. No. C30010009) or 1.5 ml TPX tubes (Diagenode, Cat. No. C30010010)
- PBS
- Mnase (Sigma-Aldrich, Cat. No. n3755). Reconstitute in ultra pure water at 0.5U/μl.
- 0.1 M CaCl<sub>2</sub>
- RIPA buffer: commercially available or homemade buffer
  - 50mM tris pH 7,5
  - 150 mM NaCl
  - 1% NP40
  - 0,5% Na-deoxycholate
  - 0,1% SDS

### 2. Procedure

- Wash cells with PBS and pellet cells by centrifugation in 15 ml TPX tubes or 1,5 ml TPX tubes (depending on the volume: link to the Diagenode tube guide).
- Add 3x the volume of ice-cold RIPA buffer.
- Incubate 10 minutes on ice.
- Add reconstitute Mnase to a final concentration of 0.0025U/μl.
- Add CaCl<sub>2</sub> to a final concentration of 1 mM.
- Incubate 15 minutes at 37°C.
- Pre-cooled Bioruptor® at 4°C with the water cooling system (Diagenode, Cat. No. B02010002 or B02010003) or use cold distilled water.
- Install samples in Bioruptor® and sonicate at High Power for 8 cycles (sonication cycle: 15 sec ON, 60 sec OFF).
- Centrifuge at 13,200 rpm for 10 minutes at 4°C.
- Collect supernatant and discard pellet.
- Determine protein concentration.
- Store aliquots at -20°C or proceed with the western blot analysis. The use of 10 to 30ug is recommended.

### 3. Selected publications

1. Mozzetta C, Pontis J, Fritsch L, Robin P, Portoso M, Proux C, et al. The histone H3 lysine 9 methyltransferases G9a and GLP regulate polycomb repressive complex 2-mediated gene silencing. *Mol Cell*. 2014 Jan 23;53(2):277–89.
2. Guasconi V, Pritchard L-L, Fritsch L, Mesner LD, Francastel C, Harel-Bellan A, et al. Preferential association of irreversibly silenced E2F-target genes with pericentromeric heterochromatin in differentiated muscle cells. *Epigenetics*. 2010 Nov;5(8):704–9.
3. Fritsch L, Robin P, Mathieu JRR, Souidi M, Hinaux H, Rougeulle C, et al. A subset of the histone H3 lysine 9 methyltransferases Suv39h1, G9a, GLP, and SETDB1 participate in a multimeric complex. *Mol Cell*. 2010 Jan 15;37(1):46–56.



**Using Bioruptor® in a creative way?**

Share your protocol with us at [info@diagenode.com](mailto:info@diagenode.com)

