

Instruction Manual



# Protein Extraction kit

Cat. No. C20000020 (AL-002-0050)

Version 1 / 14.10.2013

diagenode



# Content

Introduction	4
General remarks before starting	4
Required materials and reagents	5
Precautions	5
Protocol	6

## Introduction

Protein extraction from tissues is the first step for many biochemical and analytical techniques (PAGE, Western blotting, mass spectrometry, etc.) or protein purification. Efficient disruption and homogenization of animal tissues are required to ensure high yields of proteins. Diagenode's Bioruptor® uses state-of-the-art ultrasound technology to efficiently disrupt and homogenize tissues in just one step. Bioruptor® offers unique benefits for tissue disruption and homogenization:

- Fast and simple
- No contamination between samples
- Efficient
- Gentle processing
- Reproducible
- Temperature controlled
- Multiplexing capability
- 15 ml TPX tubes prefilled with Protein Extraction beads (Format: 50 rxns)

## General remarks before starting

- Conditions for protein extraction (e.g. use of fresh or frozen tissue, composition of extraction buffer etc.) must be adjusted according to the nature of the proteins of interest and the assays to be run. SDS might be added to the extraction buffer to maximize the yield of soluble proteins. SDS extracts can be used for SDS electrophoresis and Western blotting. It is recommended to reduce the SDS concentration for 2D electrophoresis, enzyme-linked immunosorbent assay and mass spectrometry.
- For functional studies (e.g. the study of protein-protein interactions), avoid using ionic detergents and high concentrations of salt.

*Extraction buffer: use RIPA buffer as a starting point for optimization:*

*50 mM Tris-HCl (pH 7.4)*

*150 mM NaCl*

*1% NP-40*

*0.25% Na-deoxycholate*

*Protease Inhibitor Mix*

*SDS 0.1 - 2% (optional)*

*It is always recommended to optimize the buffer composition depending on a specific research project*

- Always use Protease Inhibitor Mix during extraction procedure to block the possible protein degradation.
- Keep extracted proteins at -80°C.

## Other required materials not provided

- Bioruptor® Standard (Cat. No. B01010001) or Plus (Cat. No. B01020001)
- Bioruptor® Water Cooler (Cat. No. B02010003; 115V or B02010002; 230 V)
- Single Cycle Valve for Bioruptor® Plus (Diagenode, Cat. No. VB-100-0001)
- 1.5 ml (Cat. No. C30010010-50 or C30010010-1000) or 15 ml TPX tubes (Cat. No. C30010009) for sonication
- Tube holder for 1.5 ml tubes (Cat. No. B01200011)
- Tube holder pack for extraction kits (Cat. No. B01200015)
- Protease Inhibitor Mix (Cat. No. C12010012 or C12010011)
- Buffer for protein extraction from tissue (not supplied)
- Reagents for protein quantification (optional)

## Precautions

- This product is intended for laboratory research use only.  
**CAUTION:** Not for diagnostic use. The safety and efficacy of this product in diagnostic or other clinical uses has not been established.
- RNA extraction reagent is harmful and not intended for medical use. Handle with great care.
- When using, wear appropriate protective gear (gloves, goggles, etc.).
- If the product enters eye or adheres to skin, wash with large amounts of water for at least 15 minutes and consult a doctor.
- Handle this product in accordance with the manual.
- Diagenode is not responsible for problems caused if this product is not handled in accordance with the manual.

# Protocol

## I. Protein extraction from Tissues

- » *This protocol has been validated for up to 50 mg of tissue. Do not use more tissue per sample. For larger quantity cut the tissue and proceed to the disruption in separate tubes. When proceeding 20 - 50 mg of tissue 15 ml TPX tubes are recommended with a final volume of 1 - 2 ml. Less tissue could be sonicated in 1.5 ml TPX tubes with a final volume of 100 - 300 µl.*
  - » *Minimize the time of tissue collection to prevent protein degradation.*
  - » *Dissected tissues can be snap-frozen in liquid nitrogen and stored at -80°C until protein extraction*
1. Pre-cool Bioruptor® to 4°C using the Bioruptor® Water Cooler (Cat. No. B02010003; 115V or B02010002; 230 V).
  2. Add Protease Inhibitor Mix (200x) to the cold protein extraction buffer: 5 µl per 1 ml of extraction buffer. Scale accordingly.
  3. Add the required volume of a cold extraction buffer to the TPX tubes pre-filled with Protein Extraction Beads.
  4. Add tissue pieces to the TPX tubes. Make sure that the final volume is in the recommended range: 100 - 300 µl for 1.5 ml TPX microtubes and 1 - 2 ml for 15 ml TPX tubes.
  5. Vortex tubes briefly and proceed to sonication by using the Bioruptor® with the following settings:  
Power: H position (High)  
Sonication cycle: 30 sec ON/30 sec OFF  
Total sonication time: 5 - 15 cycles  
Temperature: 4°C
    - » *To guarantee homogeneity of sonication, the tube holder should be always completely filled with tubes.*
  6. Stop the Bioruptor® after each 5 cycles, vortex samples and check the sample visually for disruption.
    - » *Please note that the optimization might be required depending on the sample format (fresh or frozen tissue), tissue type and tissue amount. The shortest sonication time should be chosen to prevent protein damage. Incomplete disruption may occur with fibrous tissues (i.e. muscles).*
  7. Transfer the supernatant to a new tube and centrifuge samples at 14,000 rpm for 15 min at 4°C to remove any remaining insoluble material.
    - » *The Protein Extraction Beads might be washed once with extraction buffer for maximum recovery of total protein but this will lead to the sample dilution.*
  8. Transfer the supernatant containing soluble proteins to a new tube.
  9. Take an aliquot for the quantification and the further analysis if needed. Store proteins extracts in small aliquots at -80°C.
    - » *Different protein concentration assays exist including: absorbance at 280 nm, Lowry Assay, Bradford Assay, Bicinchoninic Assay (BCA) etc.. Many commercial kits for protein quantification are also available. Please note that measuring the protein concentration in an SDS extract requires that the assay is compatible with the detergent and reducing agent in the solution.*

## II. Protein extraction from Cultured Cells

- » *This protocol has been validated using RIPA buffer but it may be necessary to optimize the buffer*

composition depending on a specific research project.

- » We recommend using 100 µl of an appropriate lysis buffer per  $1 \times 10^6$  cells.
- » For Western blotting, cells might be lysed directly in 1x Laemmli buffer. After sonication, centrifuge extract at 14,000 rpm for 15 min. Transfer the supernatant to a new tube and boil for 3 min. The supernatant can be used in Western blot. Note that protein quantification by common methods is not compatible with Laemmli buffer.

1. Pre-cool Bioruptor® to 4°C using the Bioruptor® Water Cooler (Cat. No. B02010003; 115V or B02010002; 230 V).
2. Add Protease Inhibitor Mix (200x) to the ice-cold cell lysis buffer: 5 µl per 1 ml of extraction buffer. Scale accordingly.
3. **For monolayer cells:**  
Rinse the monolayer cells 3 times with cold PBS. For the final rinse, use a cell scraper and transfer the cell suspension to a TPX tube. Centrifuge cells at 1,500 rpm for 10 min at 4°C and aspirate as much supernatant as possible. Proceed to step 4.

**For suspension cells:**

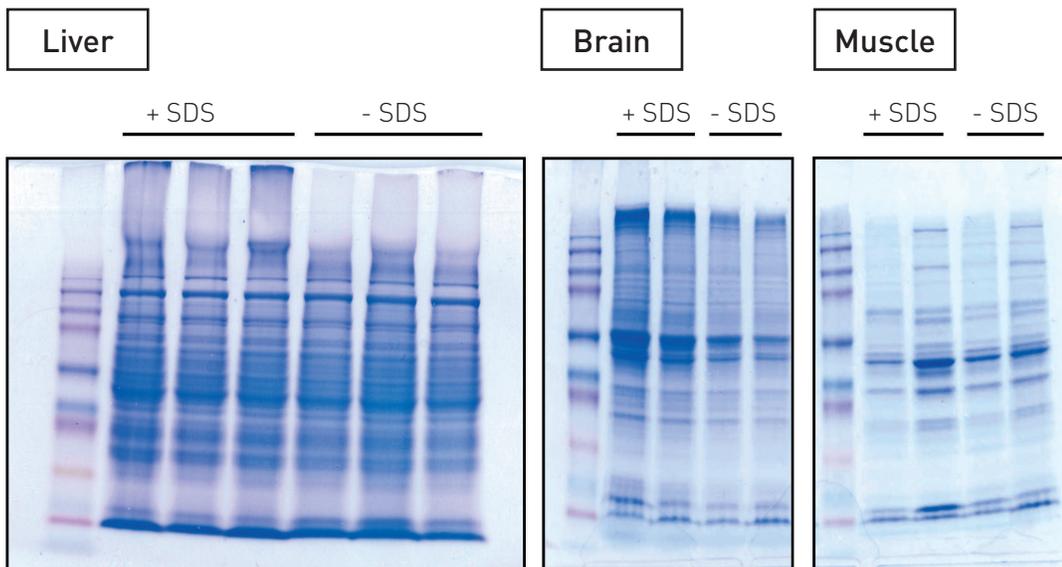
Centrifuge suspension at 1,500 rpm for 10 min at 4°C and aspirate the supernatant. Resuspend the pellet in cold PBS, transfer to a TPX tube and centrifuge at 1,500 rpm for 10 min at 4°C. Aspirate the supernatant. Repeat 2 more times. Proceed to the step 4.

4. Add ice-cold cell lysis buffer and resuspend the pellet. Incubate on ice for 10 min.
  - » *The viscosity may appear at this step*
5. Vortex tubes briefly and proceed to sonication by using the Bioruptor® with the following settings:  
*Power: H position (High)*  
*Sonication cycle: 30 sec ON/30 sec OFF*  
*Total sonication time: 5-10 cycles*  
*Temperature: 4°C*
  - » *To guarantee homogeneity of sonication, the tube holder should be always completely filled with tubes.*
6. Stop the Bioruptor® after 5 cycles, briefly vortex samples and visually check the samples: Samples should be in solution (viscosity should be reduced)
  - » *Please note that the optimization might be required depending on sample format (cell density, cell type etc.). The shortest sonication time should be chosen to prevent protein damage.*
7. Transfer the supernatant to a new tube and centrifuge samples at 14,000 rpm for 15 min at 4°C to remove any remaining insoluble material.
8. Take an aliquot for the quantification and the further analysis if needed. Store protein extracts at -80°C.
  - » *Different protein concentration assays exist including: absorbance at 280 nm, Lowry Assay, Bradford Assay, Bicinchoninic Assay (BCA) etc. Many commercial kits for protein quantification are also available. Please note that measuring the protein concentration in an SDS extract requires that the assay is compatible with the detergent and reducing agent in the solution.*



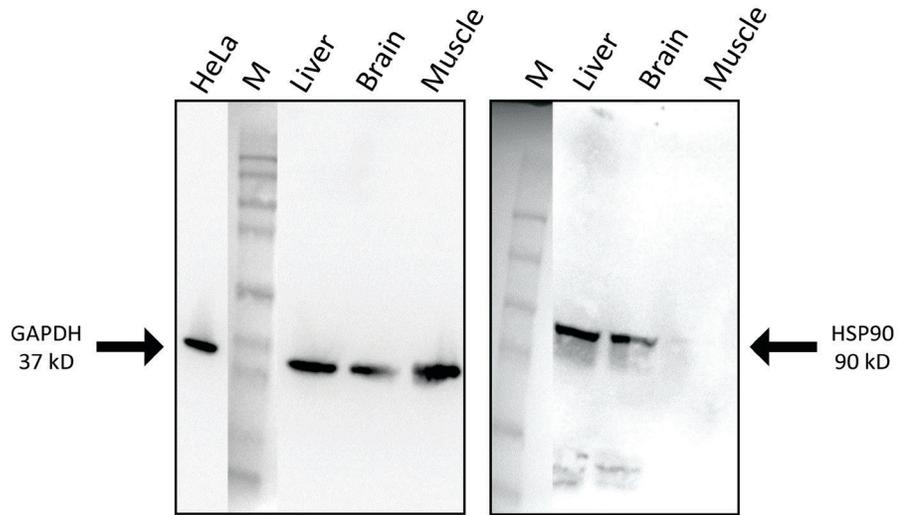
**Figure 1. Protein Extraction Beads are required for efficient tissue disruption using the Bioruptor® Plus**

Complete disruption is observed in the sample containing Diagenode's Protein Extraction Beads (left) after 5 cycles while non-disrupted tissue is still present in the sample without the Protein Extraction Beads (right).



**Figure 2. Total proteins effectively extracted from tissues using Bioruptor® Plus**

Various mouse tissues were disrupted in RIPA buffer supplemented with or without 2% SDS. Total proteins were separated by SDS-PAGE and stained with Coomassie Blue dye.



**Figure 3. Western blot analysis of GAPDH and HSP90 proteins in tissue and cultured cell extracts.**

Expected bands of 37 kD and 90 kD are observed for GAPDH (left panel) and HSP90 (right panel), respectively, in liver, brain and skeletal muscle. Note that HSP90 is expressed in muscle in an extremely low level (H. Quraishi and I. R. Brown, *J Neurosci Res.* 1996 Feb 1;43(3):335-45). Whole cell extract from HeLa cells is loaded as positive control. HeLa cells were lysed using the Bioruptor® Plus.







[orders@diagenode.com](mailto:orders@diagenode.com)  
[orders.na@diagenode.com](mailto:orders.na@diagenode.com)  
[www.diagenode.com](http://www.diagenode.com)