

Protein Sample Preparation

Protocol

Bulletin 6195

Human Cells

Use the MicroRotor™ cell lysis kit (mammalian) or the protocol detailed here, which uses sonication and radioimmunoprecipitation assay (RIPA) buffer, for cell lysis and protein extraction.

Suspension Cultured Cells



- 1 Pellet the cells by centrifugation at $2,000 \times g$ for 5 min at 4°C .
- 2 Discard the supernatant and wash pelleted cells in cold PBS. Repeat steps 1 and 2 twice.
- 3 Add RIPA buffer to the pelleted cells and suspend the pellet with a pipet.

Monolayer Cultured Cells



- 1 Carefully remove (decant) culture medium from cells. Wash cells twice in cold PBS.
- 2 Add RIPA buffer to the cells and keep on ice for 5 min. Swirl the plate occasionally to spread the buffer around the plate.
- 3 Use a cell scraper to collect the lysate and transfer to a microcentrifuge tube.

- 4 Place the cell suspension on ice, incubate 5 min, and sonicate at appropriate intervals. Check lysis efficacy by light microscopy.
- 5 Centrifuge cell debris at $\sim 14,000 \times g$ for 15 min at 4°C and transfer supernatant to a new vial.
- 6 Perform a protein assay of the supernatant. A protein concentration of $3\text{--}5 \mu\text{g}/\mu\text{l}$ is best for PAGE.
- 7 Add 2x SDS-PAGE sample buffer to the protein solution to yield a 1x sample buffer concentration.

REAGENTS

- Phosphate buffered saline (PBS)
- RIPA solubilization buffer (use 1 ml RIPA buffer with 3×10^7 cells; store and use RIPA buffer at 4°C)
- SDS-PAGE sample buffer (2x)

EQUIPMENT

- Centrifuge
- Sonicator

PRODUCT LINKS

- MicroRotor cell lysis kit (mammalian)
- RC DC protein assay
- SDS-PAGE sample buffer (2x)

BIO-RAD

This is an excerpt from Bio-Rad's comprehensive Electrophoresis Guide (Bulletin 6640).



**Bio-Rad
Laboratories, Inc.**

Life Science
Group

Web site www.bio-rad.com **USA** 800 424 6723 **Australia** 61 2 9914 2800 **Austria** 01 877 89 01 **Belgium** 09 385 55 11 **Brazil** 55 11 5044 5699
Canada 905 364 3435 **China** 86 21 6169 8500 **Czech Republic** 420 241 430 532 **Denmark** 44 52 10 00 **Finland** 09 804 22 00
France 01 47 95 69 65 **Germany** 089 31 884 0 **Greece** 30 210 9532 220 **Hong Kong** 852 2789 3300 **Hungary** 36 1 459 6100 **India** 91 124 4029300
Israel 03 963 6050 **Italy** 39 02 216091 **Japan** 03 6361 7000 **Korea** 82 2 3473 4460 **Mexico** 52 555 488 7670 **The Netherlands** 0318 540666
New Zealand 64 9 415 2280 **Norway** 23 38 41 30 **Poland** 48 22 331 99 99 **Portugal** 351 21 472 7700 **Russia** 7 495 721 14 04
Singapore 65 6415 3188 **South Africa** 27 861 246 723 **Spain** 34 91 590 5200 **Sweden** 08 555 12700 **Switzerland** 061 717 95 55
Taiwan 886 2 2578 7189 **Thailand** 800 88 22 88 **United Kingdom** 020 8328 2000
