

Protein Sample Preparation

Protocol

Bulletin 6198

Microbial Cultures

Use the MicroRotor™ cell lysis kit (bacteria) or the MicroRotor cell lysis kit (yeast). Alternatively, use the protocol detailed here, which relies on cell lysis with ultrasonic waves in combination with a solubilization in SDS under elevated temperature. This ensures deactivation and denaturation of proteases.

- 1 Centrifuge cells ($\sim 5 \times 10^7$) for 3 min at $5,000 \times g$ and resuspend the pellet in an equal volume of 37°C PBS and centrifuge again. Repeat two more times to remove all interfering material (extracellular proteases and growth media).
- 2 Add $200 \mu\text{l}$ of hot (95°C) SDS sample solubilization buffer to the pellet and vortex thoroughly.
- 3 Sonicate the sample solution 10 times for 1 sec each at $\sim 60 \text{ W}$ and $\sim 20 \text{ kHz}$. Incubate at 95°C for 5 min.
- 4 Cool the sample to 20°C and dilute with $\sim 250 \mu\text{l}$ $2\times$ SDS-PAGE sample buffer. Incubate for another 20 min at room temperature.
- 5 Centrifuge the sample solution at 20°C for 30 min at $14,000 \times g$ and harvest the supernatant.
- 6 Perform the protein assay. The protein concentration should be $\sim 5 \mu\text{g}/\mu\text{l}$.

Protein Fractions from Chromatography

When checking fraction purity or the enrichment of a particular protein after a chromatographic separation, you can observe the presence of high concentrations of salt, detergent, denaturants, and organic solvents. For example, in ion exchange chromatography, proteins are eluted by a salt gradient. But, the salt concentration of the corresponding fractions can be as high as 0.5 M , a concentration that interferes with SDS-PAGE.

Remove salt and other contaminants by one of the following approaches:

- Buffer exchange — use Bio-Spin® or Micro Bio-Spin™ columns, which are filled with size exclusion media equilibrated in Tris buffer. These columns accommodate a sample volume $50\text{--}100 \mu\text{l}$ and remove compounds $<6 \text{ kD}$ within 10 min. Mix the purified sample with $2\times$ SDS-PAGE sample buffer
- Precipitation — use the ReadyPrep™ 2-D cleanup kit (based on an acetone/TCA precipitation) for simultaneous removal of interfering substances and concentration of dilute samples ($<50 \text{ ng/ml}$)

TIPS

Reproducible lysis and protein solubilization of bacteria and yeast is challenging because the cells may release proteases and other enzymes into the growth medium. Wash the cultures thoroughly with isotonic buffers and take precautions to inactivate the proteolytic activity after cell lysis. Extensive disruption of microbial cells is required, usually with the help of a French press, bead impact instruments, or sonicator

REAGENTS

- SDS sample solubilization buffer
- SDS-PAGE sample buffer ($2\times$)
- Phosphate buffered saline (PBS)

EQUIPMENT

- Centrifuge
- Sonicator

PRODUCT LINKS

- MicroRotor cell lysis kit (bacteria)
- MicroRotor cell lysis kit (yeast)
- RC DC protein assay
- SDS-PAGE sample buffer ($2\times$)
- Bio-Spin Columns
- Micro Bio-Spin Columns
- ReadyPrep 2-D cleanup kit

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
