

ab289882 – Nuclear/Cytosol Fractionation Kit

For isolation of nuclear/cytosol fractions without ultracentrifugation.

For research use only - not intended for diagnostic use.

For overview, typical data and additional information please visit:

<http://www.abcam.com/ab289882>

Storage and Stability

The entire kit can be stored at -20°C. After opening the kit, store all Buffers, Protease Inhibitor Cocktail I and DTT II at -20°C.

Materials Supplied

Item	Quantity	Storage Condition
Extraction Buffer VI	20 mL	-20°C
Lysis Buffer X	1.2 mL	-20°C
Extraction Buffer V	10 mL	-20°C
DTT II	100 µl	-20°C
Protease Inhibitor Cocktail I	1 vial	-20°C

Reagent Preparation

- Always keep all buffers on ice during the experiment.
- All centrifugation procedures should be performed at 4°C.

Protease Inhibitor Cocktail I: Add 250 µL DMSO to dissolve the 500X Protease Inhibitor Cocktail I before use.

Extraction Buffer V: Before starting the procedure, prepare enough Extraction Buffer V and Extraction Buffer VI for your experiment: Add 2 µL Protease Inhibitor Cocktail I and 1 µL DTT II to each of 1 mL of Extraction Buffer V and each of 1 mL of Extraction Buffer VI, individually.

ΔNote: If you plan on using the Histone Acetyltransferase Activity Assay Kit, ab65352, OMIT the DTT II as this will interfere with the Histone Acetyltransferase Activity Assay.

Assay Protocol

- This protocol is described for fractionation of up to 2 x 10⁶ cells. The procedure is also applicable for large-scale preparations (e.g., up to 10⁹ cells) by scaling up the volume.
- 1. Collect cells by centrifugation at 600 x g for 5 min at 4°C.
- 2. Add 0.2 mL Extraction Buffer VI Mix containing DTT II and Protease Inhibitors (prepared as in Reagent Preparation section). If using tissue samples, cut the tissue (100-200 mg) into small pieces, add ice cold PBS (1-2 mL), and homogenize in a tissue homogenizer. Pellet the cells by centrifugation at 500 x g for 2- 3 min and remove the supernatant. Add 0.2 mL of the Extraction Buffer VI.
- 3. Vortex vigorously on the highest setting for 15 sec to fully resuspend the cell pellet. Incubate the tube on ice for 10 min.
- 4. Add 11 µL of ice-cold Lysis X to the tube. Vortex 5 sec on the highest setting. Incubate on ice for 1 min.
- 5. Vortex 5 sec on the highest setting. Centrifuge the tube for 5 min at maximal speed in a microcentrifuge (16,000 x g).
- 6. Immediately transfer the supernatant (Cytoplasmic extract) fraction to a clean prechilled tube. Place the tube on ice.
- 7. Resuspend the pellet (contains nuclei) in 100 µL of ice-cold Extraction Buffer V (prepared as in Reagent Preparation section).
- 8. Vortex on the highest setting for 15 sec. Return the sample to ice.
- 9. Repeat Step 8 for every 10 min for a total 40 min.

10. Centrifuge the tube at full speed (16,000 x g) in a microcentrifuge for 10 min.
11. Immediately transfer the supernatant (Nuclear extract) to a clean pre-chilled tube. Place on ice. Store extract at -80°C for future use.

ΔNote: Nuclear extract prepared using the above procedure contains proteins in a concentration ~1 mg/mL. If higher concentration is desired, the nuclei can be resuspended in less volume of Extraction Buffer V (such as 20 µL) in Step 7.

Technical Support

Copyright © 2025 Abcam. All Rights Reserved. The Abcam logo is a registered trademark. All information / detail is correct at time of going to print.
For all technical or commercial enquiries please go to:

www.abcam.com/contactus

www.abcam.cn/contactus (China)

www.abcam.co.jp/contactus (Japan)