

ab152163 –

Cell Lysis Buffer

Instructions for Use

For the preparation of non-denaturing cell lysate from adherent or suspension cells

This product is for research use only and is not intended for diagnostic use.

Table of Contents

1. Product Overview	3
2. Additional Materials Required	3
3. Directions for Usage	4

1. Product Overview

ab152163 is a 10X Cell Lysis Buffer to lyse adherent or suspension cells under non-denaturing conditions. It is recommended for preparing samples for use in western blotting and immunoprecipitation.

2. Additional Materials Required

- Sterile/distilled water
- PBS
- Phenylmethylsulfonyl fluoride (PMSF)

3. Directions for Usage

1. Thaw 10X Cell Lysis Buffer
2. Dilute in Sterile/DI water to achieve 1 X Cell Lysis Buffer
3. Add Phenylmethylsulfonyl fluoride (PMSF) to a final concentration of 1 mM.

Note: Make up fresh solution before each use.

4. Preparation of Non-denaturing Cell Lysate:

4.1. Adherent cells

- 4.1.1. Grow cells until they reach ~90% confluency.
- 4.1.2. Wash flasks twice with PBS.
- 4.1.3. Add appropriate vol. 1x Cell Lysis Buffer containing 1 mM PMSF to cells.
- 4.1.4. Transfer cell lysates into Eppendorf tubes and sonicate for 10–15 seconds.
- 4.1.5. Spin at 14,000 rpm, 4°C for 10 minutes.
- 4.1.6. Aliquot the supernatant into new Eppendorf tubes proceed with assay or store at -80°C.

4.2. **Suspension cells**

- 4.2.1. Harvest and count cells
- 4.2.2. Wash cells twice with PBS.
- 4.2.3. Add appropriate volume of 1x Cell Lysis Buffer containing 1 mM PMSF to cells.
- 4.2.4. Transfer cell lysates into Eppendorf tubes and sonicate for 10–15 seconds.
- 4.2.5. Spin at 14,000 rpm, 4°C for 10 minutes.
- 4.2.6. Aliquot the supernatant into new Eppendorf tubes and proceed with assay or store at -80°C.

UK, EU and ROW

Email: technical@abcam.com
Tel: +44 (0)1223 696000
www.abcam.com

US, Canada and Latin America

Email: us.technical@abcam.com
Tel: 888-77-ABCAM (22226)
www.abcam.com

China and Asia Pacific

Email: hk.technical@abcam.com
Tel: 108008523689 (中國聯通)
www.abcam.cn

Japan

Email: technical@abcam.co.jp
Tel: +81-(0)3-6231-0940
www.abcam.co.jp