

ab286872 – EZLys Tissue Protein Extraction Reagent

For extraction of total protein from tissue samples.
For research use only - not intended for diagnostic use.

For overview, typical data and additional information please visit:

<http://www.abcam.com/ab286872>

Storage and Stability

Store buffer at 4°C, protected from light. Read the entire protocol before performing the assay.

Materials Supplied

Item	Storage Condition
Tissue Protein Extraction Reagent	+4 °C

Materials Required, Not Supplied

These materials are not included in the kit, but will be required to successfully utilize this assay:

- DNase I, protease inhibitor cocktail

Tissue Protein Extraction Reagent Protocol

1. Weigh the tissue sample. For each gram of tissue, use approximately 1015 ml of cold EZLys Tissue Protein Extraction Reagent and 100150 U of DNase I. For more concentrated protein extract, use 4 ml of Tissue Protein Extraction Reagent per gram of tissue.
Δ Note: In order to avoid protein degradation, we recommend that you add protease inhibitor cocktail.
2. Homogenize using Dounce homogenizer. Centrifuge at 10,000 X g at 4°C for 10 min to remove the debris.
3. Collect the supernatant. Determine the protein concentration using Bradford. Continue with downstream applications. For future use, aliquot the tissue protein extract, snap-freeze in liquid nitrogen and store at 70°C.

Technical Support

Copyright © 2021 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)