

T-PER™ Tissue Protein Extraction Reagent

78510

Pub. No. MAN0011386

Rev. B.0

Pub Part No. 2160848.5

Number	Description
78510	T-PER Tissue Protein Extraction Reagent , 500mL, contains a proprietary detergent in 25mM bicine, 150mM sodium chloride; pH 7.6

Storage: Store product at room temperature. Product shipped at ambient temperature.

Introduction

The Thermo Scientific T-PER Tissue Protein Extraction Reagent is for total protein extraction from tissue samples. The simple composition of this reagent provides versatility because components such as protease inhibitors, salts, reducing agents or chelating agents may be added to the reagent before proceeding with cell lysis. The prepared cell lysate may be used for reporter assays (e.g., luciferase, β -galactosidase, chloramphenical acetyltransferase), protein kinase assays (e.g., PKA, PKC, tyrosine kinase), immunoassays (e.g., Western blots, ELISAs, RIAs) and protein purifications.

T-PER Reagent is mild, requires minimal amount of detergent for cell lysis, and the detergent is dialyzable. Additionally the reagent is compatible with standard protein assays such as Thermo Scientific Pierce BCA Protein Assay (Product No. 23225) and Coomassie Plus (Bradford) Assay Kit (Product No. 23236).

Procedure for Tissue Lysis

1. If desired, add protease inhibitors (e.g., Thermo Scientific Halt Protease Inhibitor Cocktail, EDTA-Free, Product No. 87785) to the T-PER Reagent just before use.
2. Weigh tissue samples. Use a ratio of ~1g of tissue to 20mL T-PER Reagent. Use a smaller volume of reagent if a more concentrated protein extract is required.
3. Add the appropriate amount of T-PER Reagent to the tissue sample and homogenize.
4. Centrifuge the sample at $10,000 \times g$ for 5 minutes to pellet cell/tissue debris.
5. Collect supernatant and continue with downstream analysis or further purification.

Related Products

XP04200BOX	Novex™ Tris-Glycine protein gels (see thermofisher.com/proteingels for a complete listing)
NW04120BOX	Bolt™ Bis-Tris Plus protein gels (see thermofisher.com/proteingels for a complete listing)
90083	DNase I Solution (2,500U/mL)
87785	Halt™ Protease Inhibitor Cocktail, EDTA-Free (100X) , 1mL
87786	Halt Protease Inhibitor Cocktail , contains sufficient reagents to treat 100mL of sample
78501	M-PER™ Mammalian Protein Extraction Reagent , 250mL
78248	B-PER™ Bacterial Protein Extraction Reagent , 500mL
89826	NE-PER™ Nuclear and Cytoplasmic Extraction Reagents
78853	Y-PER™ Yeast Protein Extraction Reagent , 500mL
23236	Pierce™ Coomassie Plus (Bradford) Assay Kit
23225	Pierce BCA Protein Assay Kit

Product References

- Jepsen, K.H., *et al.* (2002). A syndrome of joint laxity and impaired tendon integrity in lumican and fibrodulon deficient mice. *J Biol Chem* **277**:35532-40.
- Lukashevich, I.S., *et al.* (2003). Arenavirus-mediated liver pathology: acute lymphocytic choriomeningitis virus infection of *Rhesus macaques* is characterized by high-level interleukin-6 expression and hepatocyte proliferation. *J Virol* **77**:1727-37.
- Sheng, J.G., *et al.* (2002). Disruption of corticocortical connections ameliorates amyloid burden in terminal fields in a transgenic model of Abeta amyloidosis. *J Neurosci* **22**(22):9794-9.

For research use only. Not for use in diagnostic procedures.

This product ("Product") is warranted to operate or perform substantially in conformance with published Product specifications in effect at the time of sale, as set forth in the Product documentation, specifications and/or accompanying package inserts ("Documentation") and to be free from defects in material and workmanship. Unless otherwise expressly authorized in writing, Products are supplied for research use only. No claim of suitability for use in applications regulated by FDA is made. The warranty provided herein is valid only when used by properly trained individuals. Unless otherwise stated in the Documentation, this warranty is limited to one year from date of shipment when the Product is subjected to normal, proper and intended usage. This warranty does not extend to anyone other than the original purchaser of the Product ("Buyer").

No other warranties, express or implied, are granted, including without limitation, implied warranties of merchantability, fitness for any particular purpose, or non infringement. Buyer's exclusive remedy for non-conforming Products during the warranty period is limited to replacement of or refund for the non-conforming Product(s).

There is no obligation to replace Products as the result of (i) accident, disaster or event of force majeure, (ii) misuse, fault or negligence of or by Buyer, (iii) use of the Products in a manner for which they were not designed, or (iv) improper storage and handling of the Products.

Current product instructions are available at thermofisher.com. For a faxed copy, call 800-874-3723 or contact your local distributor.

© 2017 Thermo Fisher Scientific Inc. All rights reserved. Unless otherwise indicated, all trademarks are property of Thermo Fisher Scientific Inc. and its subsidiaries. Printed in the USA.