

Pierce™ Protease Inhibitor XL Capsules, EDTA-Free

Catalog Numbers A37989

Doc. Part No. 2162709 Pub. No. MAN0017533 Rev. A.0

WARNING! Read the Safety Data Sheets (SDSs) and follow the handling instructions. Wear appropriate protective eyewear, clothing, and gloves. Safety Data Sheets (SDSs) are available from thermofisher.com/support.

Product description

The Thermo Scientific™ Pierce™ Protease Inhibitor XL Capsules, EDTA-Free protect proteins from degradation by endogenous proteases released during protein extraction and purification procedures. This large-volume format is provided in a capsule where the user twists the capsule to release the particles into solution. Each capsule is sufficient for 500 mL of solution and is a broad-spectrum formulation (Table 1) that inhibits against the major classes of proteases for better protection than phenylmethylsulfonyl fluoride (PMSF) alone (Figure 1).

Table 1 Protease inhibitors included in the tablets.

Inhibitor	Target Protease
AEBSF	Serine proteases
Bestatin	Aminopeptidases
E-64	Cysteine proteases
Pepstatin A	Aspartic acid proteases

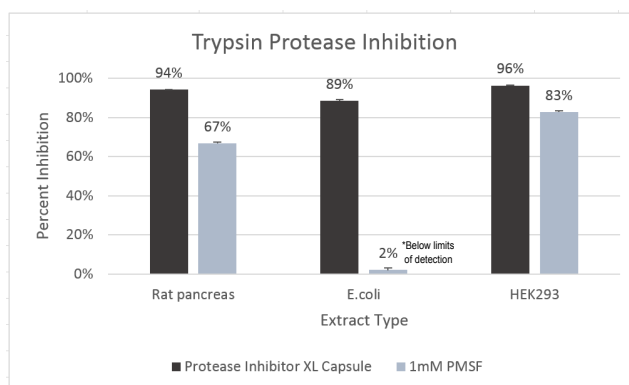


Fig. 1 Performance comparison between Pierce™ Protease Inhibitor XL Capsules, EDTA-Free and PMSF.

Contents

Cat. No.	Product	Contents	Storage
A37989	Pierce™ Protease Inhibitor XL Capsules, EDTA-Free	10 capsules - each tablet sufficient for 500 mL of extract	Store at 4°C.

Additional information

- Tablets are compatible with common protein assays and Thermo Scientific™ Pierce™ Cell Lysis Reagents.
- The outside capsule is not intended to be dissolved. Make sure to gently open the capsule to release the contents for use. However, if the whole capsule is accidentally put into solution, the capsule will dissolve, but at a much slower rate than the contents alone.
- The capsules are not recommended for volumes less than 500 mL.

Use inhibitor capsule

- Gently twist the capsule to release inhibitor particles into 500 mL of solution.
- Place solution on a magnetic stir plate with a stir bar to dissolve.

Troubleshooting

Observation	Possible cause	Recommended action
Capsule not protecting against a protease family.	Inhibitor for a specific family was not a cocktail component.	Consider using another Thermo Scientific™ protease inhibitor product, such as Halt™ Protease Inhibitor Cocktail, or add individual inhibitors for the specific protease to the samples.
	A specific inhibitor was not at the required concentration.	Consider using another Thermo Scientific™ protease inhibitor product, such as Halt™ Protease Inhibitor Cocktail, or add individual inhibitors for the specific protease to the samples.

Limited product warranty

Life Technologies Corporation and/or its affiliate(s) warrant their products as set forth in the Life Technologies' General Terms and Conditions of Sale found on Life Technologies' website at www.thermofisher.com/us/en/home/global/terms-and-conditions.html. If you have any questions, please contact Life Technologies at www.thermofisher.com/support.



Manufacturer: Pierce Biotechnology, Inc. | Thermo Fisher Scientific | 3747 N. Meridian Road | Rockford, Illinois 61101 USA

The information in this guide is subject to change without notice.

DISCLAIMER: TO THE EXTENT ALLOWED BY LAW, LIFE TECHNOLOGIES AND/OR ITS AFFILIATE(S) WILL NOT BE LIABLE FOR SPECIAL, INCIDENTAL, INDIRECT, PUNITIVE, MULTIPLE, OR CONSEQUENTIAL DAMAGES IN CONNECTION WITH OR ARISING FROM THIS DOCUMENT, INCLUDING YOUR USE OF IT.

Important Licensing Information: These products may be covered by one or more Limited Use Label Licenses. By use of these products, you accept the terms and conditions of all applicable Limited Use Label Licenses.

©2018 Thermo Fisher Scientific Inc. All rights reserved. All trademarks are the property of Thermo Fisher Scientific and its subsidiaries unless otherwise specified.