

Pierce™ Phosphatase Inhibitor Mini Tablets

Catalog Numbers A32957

Doc. Part No. 2162615 Pub. No. MAN0016070 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™ Phosphatase Inhibitor Mini Tablets protect phosphoproteins from degradation by endogenous phosphatases released during protein extraction and purification. Each tablet contains a mixture of four broad-specificity inhibitors that target serine, threonine, and tyrosine phosphatases (Table 1).

Table 1 Phosphatase inhibitors included in the tablets.

Inhibitor	Target Protease/Phosphatase
Sodium fluoride	Serine and threonine phosphatases
Sodium orthovanadate	Tyrosine phosphatases
Sodium pyrophosphate	Serine and threonine phosphatases
β-glycerophosphate	Serine and threonine phosphatases

Contents

Cat. No.	Product	Contents	Storage
A32957	Pierce™ Phosphatase Inhibitor Mini Tablets	20 tablets - each tablet sufficient for 10 mL of extract <i>(Replaces Pierce Phosphatase Inhibitor Tablets (Product No. 88667))</i>	Store at 4°C.

Additional information

- Each tablet can prepare 10 mL of 1X solution; however, samples containing particularly high levels of phosphatase activity may require a final concentration of 2-3X.
- The tablets are compatible with common protein assays and Thermo Scientific™ Pierce™ Cell Lysis Reagents.

Use inhibitor tablet

- Dispense tablet from vial and place into 10 mL of solution.
- Vortex to dissolve.

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.

Revision history: Pub. No. MAN0016070

Revision	Date	Description
A.0	12 Aug 2016	New manual

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.

Corporate entity: Life Technologies Corporation | Carlsbad, CA 92008 USA | Toll Free in USA 1 800 955 6288

©2016 Thermo Fisher Scientific Inc. All rights reserved. All trademarks are the property of Thermo Fisher Scientific and its subsidiaries unless otherwise specified.