INSTRUCTIONS



1988.4

Halt Protease Inhibitor Single-Use Cocktail EDTA-Free

78425

Number Description

78425

Halt Protease Inhibitor Single-Use Cocktail EDTA-Free (100X), $100\mu L \times 24$ microtubes, sufficient to treat 240mL of sample

Storage: Upon receipt store at 4°C. Product shipped on ice pack.

Introduction

The Thermo ScientificTM HaltTM Protease Inhibitor Single-Use Cocktail EDTA-Free protects proteins from degradation by endogenous proteases that may be released during protein extraction and purification. The ready-to-use cocktail contains several potent inhibitors (see Table 1) in DMSO that inhibit the major classes of proteases including aspartic acid, cysteine and serine proteases. The refrigerated, single-use format maintains the potency of inhibitors by eliminating freeze-thaw cycles. Also, each microtube conveniently contains 100µL that can treat up to 10mL of lysate.

Important Product Information

- Halt Protease Inhibitor Single-Use Cocktail is a broad-spectrum protease inhibitor and does not contain phosphatase inhibitors. The protease inhibitors used in the formulation at their 1X working concentrations are listed in Table 1.
- The protease inhibitor cocktail is supplied as a 100X stock in DMSO and is generally effective when used at a final concentration of 1X; however, some sample types contain particularly high levels of proteases and optimization of the cocktail's working concentration may be required.
- This cocktail is typically a liquid at 4°C but occasionally might be frozen. Frozen product does affect the function or performance of the cocktail. If frozen, equilibrate the tube to room temperature or hold the tube between thumb and forefinger until the cocktail becomes a liquid.

Table 1. Protease inhibitors included in the Halt Cocktail.		
Inhibitor	<u>1X</u>	
AEBSF	1mM	
Aprotinin	800nM	
Bestatin	50µM	
E64	15μM	
Leupeptin	20μΜ	
Pepstatin A	10μΜ	



Procedure

Note: Wear proper personal protective equipment when working with this DMSO-containing protease inhibitor cocktail.

 Remove one strip of cocktail from the foil pouch. Use scissors to carefully cut off one tube from the strip. Return unused tubes to 4°C immediately after removing the required tubes. Do not allow unused cocktail to equilibrate to room temperature.

Note: To preserve the integrity of the foil seals, use scissors to cut the strip. Do not tear the tubes from the strip as this will compromise the seals of neighboring tubes.

- 2. Equilibrate the tube to room temperature.
- 3. Vortex the tube before use to ensure a homogeneous suspension. Puncture the foil seal with a pipette tip.
- 4. Add Halt Protease Inhibitor Single-Use Cocktail EDTA-Free, at 10μL/mL directly to the lysis buffer or cell extract to produce a 1X final concentration.

Note: Up to 10mL of lysis buffer or cell extract can be treated with one tube of cocktail.

Troubleshooting

Problem	Possible Cause	Solution
Poor inhibition of protease activity	Sample contains high levels of protease	Add sufficient cocktail to produce a 2X or 3X final concentration
Cocktail does not provide sufficient protection against a particular protease family	Inhibitor for that family is not a component of the cocktail, or the cocktail does not contain the needed concentration	Add individual inhibitor to the sample for the specific protease family

Related Thermo Scientific Products

78430	Halt Protease Inhibitor Single-Use Cocktail (100X), $24 \times 100 \mu L$
87785	Halt Protease Inhibitor Cocktail, EDTA-Free (100X), 1mL
87786	Halt Protease Inhibitor Cocktail, 1mL
88660	Pierce Protease Inhibitor Tablets, 30 tablets
88661	Pierce Protease Inhibitor Tablets, EDTA-free, 30 tablets
78441	Halt Protease and Phosphatase Inhibitor Cocktail, EDTA-Free (100X), 1mL
88664	Pierce Protease and Phosphatase Inhibitor Tablets, EDTA-free, 20 tablets
78420	Halt Phosphatase Inhibitor Cocktail (100X), 1mL
88662	Pierce Phosphatase Inhibitor Tablets, 20 tablets

Products are 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"). 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 Buyer. Any model or sample furnished to Buyer is merely illustrative of the general type and quality of goods and does not represent that any Product will conform to such model or sample.

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 REPAIR, REPLACEMENT OF OR REFUND FOR THE NON-CONFORMING PRODUCT(S) AT SELLER'S SOLE OPTION. THERE IS NO OBLIGATION TO REPAIR, REPLACE OR REFUND FOR 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.

Unless otherwise expressly stated on the Product or in the documentation accompanying the Product, the Product is intended for research only and is not to be used for any other purpose, including without limitation, unauthorized commercial uses, in vitro diagnostic uses, ex vivo or in vivo therapeutic uses, or any type of consumption by or application to humans or animals.

Current product instructions are available at <u>www.thermoscientific.com/pierce</u>. For a faxed copy, call 800-874-3723 or contact your local distributor. © 2014 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.

Pierce Biotechnology	PO Box 117	(815) 968-0747	www.thermoscientific.com/pierce
3747 N. Meridian Road	Rockford, IL 61105 USA	(815) 968-7316 fax	