INSTRUCTIONS



2052.4

Halt Protease Inhibitor Cocktail, EDTA-Free

87785 78437 78439

Number	Description
87785	Halt Protease Inhibitor Cocktail, EDTA-Free (100X), 1mL
78437	Halt Protease Inhibitor Cocktail, EDTA-Free (100X), 5mL, sufficient to treat 500mL of sample
78439	Halt Protease Inhibitor Cocktail, EDTA-Free (100X), 10mL, sufficient to treat 1L of sample

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

Introduction

The Thermo ScientificTM HaltTM Protease Inhibitor Cocktail, EDTA-Free protects protein from degradation by endogenous proteases released during protein extraction and purification. This ready-to-use cocktail contains several potent inhibitors (Table 1) in DMSO against the major classes of proteases including aspartic acid, cysteine and serine.

Table 1. Protease inhibitors included in the cocktail.		
Inhibitor	1X Concentration	
AEBSF	1mM	
Aprotinin	800nM	
Bestatin	50µM	
E64	15μM	
Leupeptin	20μΜ	
Pepstatin A	10µM	

Important Product Information

- Halt Protease Inhibitor Cocktail, EDTA-Free is a broad spectrum protease inhibitor and does not contain phosphatase inhibitors. Each protease inhibitor in the cocktail is listed in Table 1 at the 1X working concentration.
- This protease inhibitor cocktail is supplied at a 100X concentration in DMSO and is generally effective when used at a 1X final concentration; however, if a sample contains particularly high levels of proteases, the effective cocktail concentration might require optimization.
- This cocktail might freeze during storage, which does not affect its function or performance. If frozen, equilibrate the bottle to room temperature until the cocktail becomes a liquid.

Procedure

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

- 1. Equilibrate the bottle to room temperature.
- 2. Vortex the bottle to ensure a homogeneous suspension.
- 3. Immediately before use, add 10µL of Halt Protease Inhibitor Cocktail per milliliter of sample directly to the lysis buffer or extract to produce a 1X final concentration.



Troubleshooting

Problem	Possible Cause	Solution
Poor inhibition of protease activity	Sample contains high levels of proteases	Add sufficient cocktail to produce a
		2X or 3X final concentration
Cocktail does not provide	Inhibitor for a specific family is not a	Add individual inhibitors to the
sufficient protection against a	component of the cocktail	samples for the specific protease
particular protease family	A specific inhibitor is not at the needed	Use the cocktail at a higher
	concentration	concentration

Related Thermo Scientific Products

78425	Halt Protease Inhibitor Single-Use Cocktail, EDTA-free (100X), $100\mu L \times 24$ microtubes
88661	Pierce TM Protease Inhibitor Tablets, EDTA-free, 30 tablets
88660	Pierce Protease Inhibitor Tablets, 30 tablets
78440	Halt Protease and Phosphatase Inhibitor Cocktail (100X), 1mL
78441	Halt Protease and Phosphatase Inhibitor Cocktail, EDTA-free (100X), 1mL
78444	Halt Protease and Phosphatase Inhibitor Cocktail (100X), $5 \times 1mL$
78445	Halt Protease and Phosphatase Inhibitor Cocktail (100X), EDTA-free, $5 \times 1 mL$
78428	Halt Phosphatase Inhibitor Single-Use Cocktail (100X), $100\mu L \times 24$ microtubes
88664	Pierce Protease and Phosphatase Inhibitor Tablets, EDTA-free, 20 tablets
66380	Slide-A-Lyzer TM Dialysis Cassettes 10K MWCO, 0.5-3 mL
89882	Zeba TM Spin Desalting Columns, 0.5mL, 25 columns
23225	Pierce BCA Protein Assay Kit
23236	Coomassie Plus (Bradford) Assay Kit
78248	B-PER™ Bacterial Protein Extraction Reagent, 500mL
78990	Y-PER TM Yeast Protein Extraction Reagent, 500mL
78833	NE-PER [™] Nuclear and Cytoplasmic Extraction Kit
78501	M-PER TM Mammalian Protein Extraction Reagent, 250mL

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.