## **INSTRUCTIONS**

# Sulfo-NHS Acetate



Number	Description
26777	Sulfo-NHS Acetate, 100mg
	Molecular Weight: 259.17

Storage: Upon receipt store at -20°C. Product shipped at ambient temperature.

## Introduction

Thermo Scientific<sup>TM</sup> Sulfo-NHS Acetate (Figure 1) is a protein modification reagent for blocking primary amines. The NHS-ester group of this reagent reacts with primary amines at pH 7.0-9.0, allowing modifications to occur in many standard non-aminecontaining buffers. Once reacted, the amine is irreversibly capped with an acyl group. For reversible amine blocking, Citraconic Anhydride (Product No. 20907) is used.

Sulfo-NHS Acetate is typically used to prevent polymerization when performing protein cross-linking reactions and when conjugating peptides to carrier proteins for immunogen production. Blocking amines on the peptide allows directed conjugation of carboxylic acids on the peptide to primary amines on the protein using EDC (Product No. 29980).

## Important Product Information

- Reconstitute Sulfo-NHS Acetate immediately before use. Do not prepare stock solutions for storage because the NHSester moiety readily hydrolyzes and becomes non-reactive. Discard any unused reconstituted reagent.
- Sulfo-NHS Acetate is moisture-sensitive. To avoid moisture condensation onto the product and subsequent hydrolysis of the reagent, the vial must be equilibrated to room temperature before opening.

## Procedure for Blocking Amines on Proteins or Peptides

#### A. Additional Materials Required

- Reaction Buffer: 0.1M sodium carbonate (NaHCO<sub>3</sub>) buffer, pH 8.5 or other amine-free buffer such as phosphate buffered saline (PBS; e.g., 100mM sodium phosphate, 0.15M NaCl; pH 7.2; Product No. 28372) or HEPES, pH 7.5-8
- Quenching Buffer (optional): 1M Tris•HCl, pH 7.5 (1M glycine or lysine may also be used) .
- Desalting Column (optional): Dextran Desalting Column, 5K MWCO (Product No. 43230) or Dialysis Unit (e.g., Thermo Scientific<sup>™</sup> Slide-A-Lyzer<sup>™</sup> Dialysis Cassette 0.1-0.5mL, 10K MWCO; Product No. 66385). For peptides larger than 2kDa, use Polyacrylamide Desalting Columns, 1.8K MWCO (Product No. 43426).

#### **B.** Procedure

Dissolve or buffer-exchange the protein or peptide sample into the Reaction Buffer. 1.

PO Box 117

Rockford, IL 61105 USA

Note: If the peptide is insoluble in the Reaction Buffer, dissolve peptide with DMSO or DMF and then dilute with an equal volume of Reaction Buffer.

2. Dissolve Sulfo-NHS Acetate in ultrapure water to 10mM (2.6mg/mL). Prepare a 25-fold molar excess of Sulfo-NHS Acetate to amino groups in the sample and add it to the reaction. Alternatively, if the number of amines in the sample is not known, adding an equal mass amount of the Sulfo-NHS Acetate to the sample will provide sufficient molar excess.

(815) 968-0747

(815) 968-7316 fax

Mix reaction and incubate for 1 hour at room temperature. 3.

Pierce Biotechnology

3747 N. Meridian Road



**Figure 1. Chemical** structure of Sulfo-NHS Acetate.





1503.2



4. If desired, add Quenching Buffer to quench any NHS esters that did not react or remove the non-reacted reagent by dialysis or desalting.

## Troubleshooting

Problem	Possible Cause	Solution
Amines not blocked	Reagent hydrolyzed	Completely equilibrate reagent vial to room temperature before opening to prevent condensation in the vial
		Dissolve reagent immediately before use – do not make a stock solution for storage
	Improper buffer was used	Use only non-amine containing buffers at pH 7-9 – at higher pH, hydrolysis is a competing reaction that reduces efficiency, and at lower pH, the reaction is extremely slow

## Additional Information Available on Our Website

- Tech Tip #18: Block amino groups to prevent polymer formation in peptide-carrier protein conjugations
- Tech Tip #3: Determine reactivity of NHS ester biotinylation and crosslinking reagents
- Tech Tip #43: Protein stability and storage
- Tech Tip #20: Dialysis: an overview

## **Related Thermo Scientific Products**

28372	BupH <sup>TM</sup> Phosphate Buffered Saline Packs, 40 packs, each pack results in 500 ml
28390	<b>BupH MES Buffer Saline Packs</b> , 10 packs, each yields 500mL of 0.1M 2-[morpholino]- ethanesulfonic acid, 0.9% NaCl, pH 4.7 when dissolved in 500mL deionized water (5L total)
77149	EDC (carbodiimide), 10mg
22980	EDC (carbodiimide), 5g
23031	Ethylenediamine Dihydrochloride, 10g
77171	Imject Bovine Serum Albumin (in MES buffer), 2mg
77110	<b>Imject Bovine Serum Albumin (in PBS)</b> , 5 × 20mg

#### **Product Reference**

Whaseon Lee-Kwon, *et al.* (2003). Lysophophatidic acid stimulates brush border Na+/H+ exchanger 3 (NHE3) activity by increasing its exocytosis by an NHE3 kinase A regulatory protein dominated mechanism. *J Biol Chem* **278**(**19**):16494-501.

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. © 2013 Thermo Fisher Scientific Inc. All rights reserved. All (other) trademarks are the property of Thermo Fisher Scientific Inc. and its subsidiaries. Printed in the USA.