

# PRODUCT INFORMATION



## SNOB 1 Reagent

Item No. 17215

CAS Registry No.: 92924-45-3

Formal Name: (3aS,4S,6aR)-hexahydro-2-oxo-N-2-propen-1-yl-1H-thieno[3,4-d]imidazole-4-pentanamide

Synonym: S-Nitrosylated Binding 1 Reagent

MF:  $C_{13}H_{21}N_3O_2S$

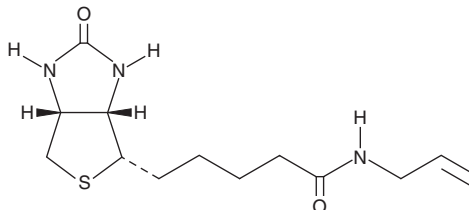
FW: 283.4

Purity:  $\geq 95\%$

Supplied as: A crystalline solid

Storage:  $-20^{\circ}\text{C}$

Stability:  $\geq 4$  years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

### Laboratory Procedures

SNOB 1 Reagent is supplied as a crystalline solid. A stock solution may be made by dissolving the SNOB 1 reagent in the solvent of choice, which should be purged with an inert gas. SNOB 1 Reagent is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of SNOB 1 reagent in these solvents is approximately 0.10, 10, and 5 mg/ml.

SNOB 1 Reagent is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, SNOB 1 reagent should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. SNOB 1 Reagent has a solubility of approximately 0.25 mg/ml in a 1:3 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

### Description

SNOB 1 Reagent is a biotinylated probe for detecting S-nitrosylated proteins in a single step. S-Nitrosylated binding (SNOB) proceeds rapidly on surface proteins or cell lysates under normal physiological conditions. Proteins tagged with SNOB 1 Reagent can be analyzed using avidin-linked probes (e.g., by immunoblot) or by mass spectrometry.

#### WARNING

THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

#### SAFETY DATA

This material should be considered hazardous until further information becomes available. Do not ingest, inhale, get in eyes, on skin, or on clothing. Wash thoroughly after handling. Before use, the user must review the [complete](#) Safety Data Sheet, which has been sent via email to your institution.

#### WARRANTY AND LIMITATION OF REMEDY

Buyer agrees to purchase the material subject to Cayman's Terms and Conditions. Complete Terms and Conditions including Warranty and Limitation of Liability information can be found on our website.

Copyright Cayman Chemical Company, 11/28/2022

#### CAYMAN CHEMICAL

1180 EAST ELLSWORTH RD  
ANN ARBOR, MI 48108 · USA

PHONE: [800] 364-9897  
[734] 971-3335

FAX: [734] 971-3640

CUSTSERV@CAYMANCHEM.COM  
WWW.CAYMANCHEM.COM