

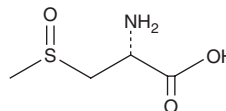
PRODUCT INFORMATION



S-Methyl-L-Cysteine-S-oxide

Item No. 17600

CAS Registry No.: 6853-87-8
Formal Name: (methylsulfinyl)-D-alanine
Synonyms: S-Methylcysteine sulfoxide, NSC 226572
MF: C₄H₉NO₃S
FW: 151.2
Purity: ≥98%
Supplied as: A crystalline solid
Storage: -20°C
Stability: ≥4 years



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

Laboratory Procedures

S-Methyl-L-cysteine-S-oxide is supplied as a crystalline solid. Aqueous solutions of S-methyl-L-cysteine-S-oxide can be prepared by directly dissolving the crystalline solid in aqueous buffers. The solubility of S-methyl-L-cysteine-S-oxide in PBS (pH 7.2) is approximately 10 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

S-Methyl-L-cysteine-S-oxide is an analog of alliin (Item No. 14001) found in some cruciferous vegetables including cabbage, turnip, cauliflower, and kale.¹ This odorless amino acid is converted through alliinase activity into a volatile thiosulfinate.²

References

1. Synge, R.L.M. and Wood, J.C. (+)-(S-Methyl-L-cysteine S-oxide) in cabbage. *Biochem. J.* **64(2)**, 252-259 (1956).
2. Krest, I., Glodek, J., and Keusgen, M. Cysteine sulfoxides and alliinase activity of some Allium species. *J. Agric. Food Chem.* **48(8)**, 3753-3760 (2000).

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, 12/02/2022

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