

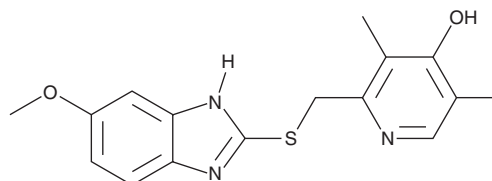
PRODUCT INFORMATION



4-hydroxy Omeprazole sulfide

Item No. 35448

CAS Registry No.: 103876-98-8
Formal Name: 2-[[[(6-methoxy-1H-benzimidazol-2-yl)thio]methyl]-3,5-dimethyl-4-pyridinol
Synonyms: 4-hydroxy OMEP sulfide, 4-hydroxy OMP sulfide, 4-hydroxy OMZ sulfide
MF: C₁₆H₁₇N₃O₂S
FW: 315.4
Purity: ≥98%
Supplied as: A 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

4-hydroxy Omeprazole sulfide is supplied as a solid. A stock solution may be made by dissolving the 4-hydroxy omeprazole sulfide in the solvent of choice, which should be purged with an inert gas. 4-hydroxy Omeprazole sulfide is soluble in DMSO.

Description

4-hydroxy Omeprazole sulfide is a metabolite of the proton pump inhibitor omeprazole (Item No. 14880).¹ It has been found in wastewater effluent.²

References

1. Hoffmann, K.J. Identification of the main urinary metabolites of omeprazole after an oral dose to rats and dogs. *Drug Metab. Dispos.* **14**(3), 341-348 (1986).
2. Hernández, F., Ibáñez, M., Gracia-Lor, E., *et al.* Retrospective LC-QTOF-MS analysis searching for pharmaceutical metabolites in urban wastewater. *J. Sep. Sci.* **34**(24), 3517-3526 (2011).

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, 09/30/2022

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