

PRODUCT INFORMATION



Omeprazole N-oxide

Item No. 35430

CAS Registry No.: 176219-04-8
Formal Name: 6-methoxy-2-[[[4-methoxy-3,5-dimethyl-1-oxido-2-pyridinyl)methyl]sulfinyl]-1H-benzimidazole

Synonyms: OMEP N-oxide, OMP N-oxide, OMZ N-oxide

MF: C₁₇H₁₉N₃O₄S

FW: 361.4

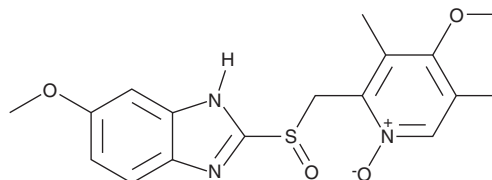
Purity: ≥98%

UV/Vis.: λ_{max}: 272, 303 nm

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

Omeprazole N-oxide is supplied as a solid. A stock solution may be made by dissolving the omeprazole N-oxide in the solvent of choice, which should be purged with an inert gas. Omeprazole N-oxide is soluble in the organic DMSO at a concentration of approximately 1 mg/ml.

Description

Omeprazole N-oxide is a metabolite of the gastric proton pump inhibitor omeprazole (Item No. 14880).¹ It is also a potential impurity found in commercial preparations of omeprazole.²

References

1. Seshadri, R.K., Raghavaraju, T.V., and Chakravarthy, I.E. A single gradient stability-indicating reversed-phase LC method for the estimation of impurities in omeprazole and domperidone capsules. *Sci. Pharm.* **81**(2), 437-438 (2013).
2. Shirasaka, Y., Sager, J.E., Lutz, J.D., et al. Inhibition of CYP2C19 and CYP3A4 by omeprazole metabolites and their contribution to drug-drug interactions. *Drug Metab. Dispos.* **41**(7), 1414-1424 (2013).

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, 11/22/2024

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