

# 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:** C<sub>17</sub>H<sub>19</sub>N<sub>3</sub>O<sub>4</sub>S

**FW:** 361.4

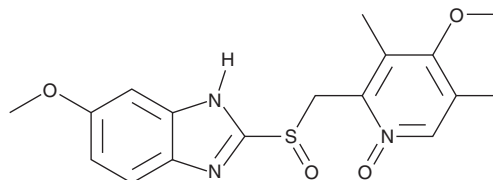
**Purity:** ≥98%

**UV/Vis.:** λ<sub>max</sub>: 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).<sup>1</sup> It is also a potential impurity found in commercial preparations of omeprazole.<sup>2</sup>

### 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, 11/09/2022

#### 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