

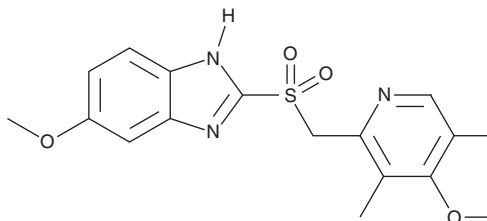
PRODUCT INFORMATION



Omeprazole sulfone

Item No. 18882

CAS Registry No.: 88546-55-8
Formal Name: 6-methoxy-2-[[[4-methoxy-3,5-dimethyl-2-pyridinyl)methyl]sulfonyl]-1H-benzimidazole
Synonyms: OMEP sulfone, OMP sulfone, OMZ sulfone
MF: C₁₇H₁₉N₃O₄S
FW: 361.4
Purity: ≥98%
UV/Vis.: λ_{max}: 273, 299 nm
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

Omeprazole sulfone is supplied as a crystalline solid. A stock solution may be made by dissolving the omeprazole sulfone in the solvent of choice, which should be purged with an inert gas. Omeprazole sulfone is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of omeprazole sulfone in ethanol is approximately 5 mg/ml and approximately 30 mg/ml in DMSO and DMF.

Omeprazole sulfone is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, omeprazole sulfone should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. Omeprazole sulfone has a solubility of approximately 0.5 mg/ml in a 1:1 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

Omeprazole sulfone is a metabolite of the inhibitor of gastric H⁺/K⁺ ATPase omeprazole (Item No. 14880).¹ It is formed from omeprazole by the cytochrome P450 (CYP) isoform CYP3A4.

Reference

1. Äbelö, A., Andersson, T.B., Antonsson, M., *et al.* Stereoselective metabolism of omeprazole by human cytochrome P450 enzymes. *Drug Metab. Dispos.* **28(8)**, 966-972 (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, 07/10/2023

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