

PRODUCT INFORMATION



Omeprazole sulfone-d₃

Item No. 26784

CAS Registry No.: 2749628-17-7
Formal Name: 6-methoxy-2-[[[4-methoxy-3,5-dimethyl-2-pyridinyl)methyl]sulfonyl]-1H-benzimidazole-4,5,7-d₃

Synonyms: OMEP sulfone-d₃,
OMP sulfone-d₃,
OMZ sulfone-d₃

MF: C₁₇H₁₆D₃N₃O₄S
FW: 364.4

Chemical Purity: ≥95% (Omeprazole sulfone)

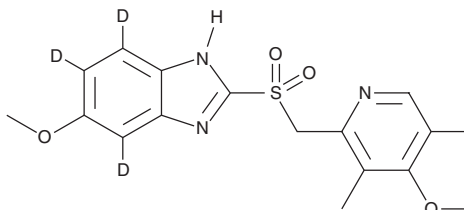
Deuterium

Incorporation: ≥99% deuterated forms (d₁-d₃); ≤1% d₀

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 sulfone-d₃ is intended for use as an internal standard for the quantification of omeprazole sulfone (Item No. 18882) by GC- or LC-MS. The accuracy of the sample weight in this vial is between 5% over and 2% under the amount shown on the vial. If better precision is required, the deuterated standard should be quantitated against a more precisely weighed unlabeled standard by constructing a standard curve of peak intensity ratios (deuterated versus unlabeled).

Omeprazole sulfone-d₃ is supplied as a solid. A stock solution may be made by dissolving the omeprazole sulfone-d₃ in the solvent of choice, which should be purged with an inert gas. Omeprazole sulfone-d₃ is slightly soluble in chloroform and methanol (warmed).

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.

References

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