

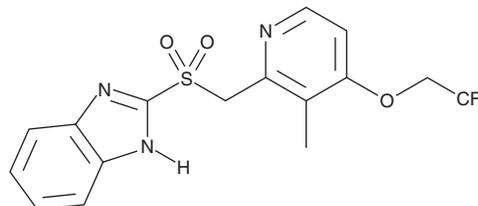
PRODUCT INFORMATION



Lansoprazole sulfone

Item No. 22569

CAS Registry No.: 131926-99-3
Formal Name: 2-[[[3-methyl-4-(2,2,2-trifluoroethoxy)-2-pyridinyl]methyl]sulfonyl]-1H-benzimidazole
Synonym: AG-1813
MF: C₁₆H₁₄F₃N₃O₃S
FW: 385.4
Purity: ≥98%
UV/Vis.: λ_{max}: 281 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

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

Lansoprazole sulfone is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, lansoprazole sulfone should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. Lansoprazole 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

Lansoprazole sulfone is a phase I metabolite of lansoprazole (Item Nos. 13627 | 21967 | 18235).¹ Lansoprazole is a proton pump inhibitor that inactivates the hydrogen/potassium-stimulated ATPase pumps in parietal cells, thus inhibiting gastric acid secretion and increasing intragastric pH.² Lansoprazole is metabolized by the cytochrome P450 (CYP) isoform CYP3A4 to form lansoprazole sulfone.¹

References

1. Meyer, U.A. Interaction of proton pump inhibitors with cytochromes P450: Consequences for drug interactions. *Yale J. Biol. Med.* **69(3)**, 203-209 (1996).
2. Schubert, M.L. Pharmacotherapy for acid/peptic disorders. *Yale J. Biol. Med.* **69(2)**, 197-201 (1996).

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, 10/24/2022

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