

PRODUCT INFORMATION

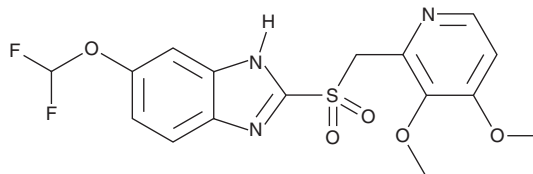


Pantoprazole sulfone

Item No. 22252

CAS Registry No.: 127780-16-9
Formal Name: 6-(difluoromethoxy)-2-[[[(3,4-dimethoxy-2-pyridinyl)methyl]sulfonyl]-1H-benzimidazole

MF: C₁₆H₁₅F₂N₃O₅S
FW: 399.4
Purity: ≥98%
UV/Vis.: λ_{max}: 287 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

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

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

Pantoprazole sulfone is a metabolite of the gastric H⁺/K⁺ ATPase pump inhibitor pantoprazole (Item No. 21345).¹ Pantoprazole is metabolized by the cytochrome P450 (CYP) isoforms CYP2C19 and CYP3A4 to form pantoprazole sulfone.

Reference

1. Shi, S., and Klotz, U. Proton pump inhibitors: An update of their clinical use and pharmacokinetics. *Eur. J. Clin. Pharmacol.* **64**(10), 935-951 (2008).

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