

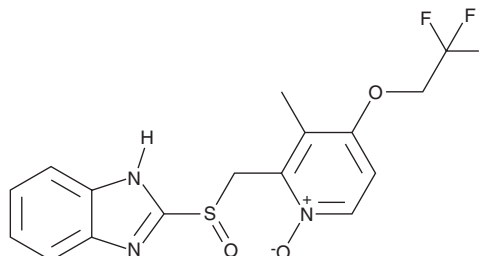
PRODUCT INFORMATION



Lansoprazole N-oxide

Item No. 35427

CAS Registry No.: 213476-12-1
Formal Name: 2-[[[3-methyl-1-oxido-4-(2,2,2-trifluoroethoxy)-2-pyridinyl]methyl]sulfinyl]-1H-benzimidazole
MF: C₁₆H₁₄F₃N₃O₃S
FW: 385.4
Purity: ≥98%
UV/Vis.: λ_{max}: 279 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

Lansoprazole N-oxide is supplied as a solid. A stock solution may be made by dissolving the lansoprazole N-oxide in the solvent of choice, which should be purged with an inert gas. Lansoprazole N-oxide is soluble in the organic solvent DMSO.

Description

Lansoprazole N-oxide is a potential impurity found in bulk preparations of the proton pump inhibitor lansoprazole.¹ It is a degradation product formed under acid or base stress.

Reference

1. Rao, P.V., Kumar, M.N., and Kumar, M.R. A novel, validated stability-indicating UPLC method for the estimation of lansoprazole and its impurities in bulk drug and pharmaceutical dosage forms. *Sci. Pharm.* **81**(1), 183-193 (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, 11/18/2024

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