

PRODUCT INFORMATION



Esomeprazole

Item No. 40851

CAS Registry No.: 119141-88-7

Formal Name: 6-methoxy-2-[(S)-[(4-methoxy-3,5-dimethyl-2-pyridinyl)methyl]sulfinyl]-1H-benzimidazole

Synonyms: (-)-Omeprazole, (S)-Omeprazole

MF: C₁₇H₁₉N₃O₃S

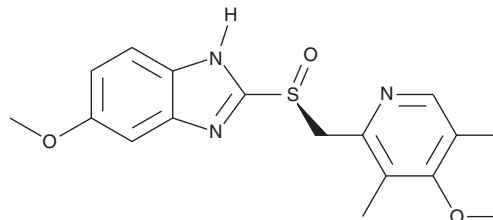
FW: 345.4

Purity: ≥98%

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

Esomeprazole is supplied as a crystalline solid. A stock solution may be made by dissolving the esomeprazole in the solvent of choice, which should be purged with an inert gas. Esomeprazole is sparingly soluble (1-10 mg/ml) in ethanol and DMSO.

Description

Esomeprazole is an inhibitor of the gastric H⁺/K⁺-ATPase and the (S) enantiomer of omeprazole (Item No. 14880).¹ It reversibly inhibits the activity of the H⁺/K⁺-ATPase in an enzyme assay. Esomeprazole (30 mg/kg, s.c.) inhibits histamine-induced gastric acid secretion in rats. Formulations containing esomeprazole in complex with magnesium have been used as proton pump inhibitors (PPIs) in the treatment of gastroesophageal reflux disease (GERD), *H. pylori* eradication, and hypersecretory conditions, as well as in the prevention of non-steroidal anti-inflammatory drug-related gastric ulcer.

Reference

1. Shin, J.M., and Sachs, G. Restoration of acid secretion following treatment with proton pump inhibitors. *Gastroenterology* **123**(5), 1588-1597 (2002).

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, 03/27/2024

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