

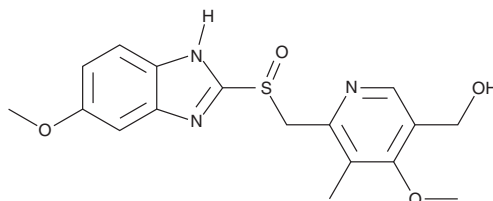
# PRODUCT INFORMATION



## 5-hydroxy Omeprazole

Item No. 10009028

**CAS Registry No.:** 92340-57-3  
**Formal Name:** 4-methoxy-6-[[[(6-methoxy-1H-benzimidazol-2-yl)sulfinyl]methyl]-5-methyl-3-pyridinemethanol  
**Synonyms:** 5-hydroxy OMEP, 5-hydroxy OMP, 5-hydroxy OMZ  
**MF:** C<sub>17</sub>H<sub>19</sub>N<sub>3</sub>O<sub>4</sub>S  
**FW:** 361.4  
**Purity:** ≥98%  
**UV/Vis.:** λ<sub>max</sub>: 205, 302 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

5-hydroxy Omeprazole is supplied as a solid. A stock solution may be made by dissolving the 5-hydroxy omeprazole in the solvent of choice, which should be purged with an inert gas. 5-hydroxy Omeprazole is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of 5-hydroxy omeprazole in these solvents is approximately 2, 10, and 20 mg/ml, respectively.

5-hydroxy Omeprazole is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, 5-hydroxy omeprazole should first be dissolved in DMF and then diluted with the aqueous buffer of choice. 5-hydroxy Omeprazole has a solubility of approximately 0.5 mg/ml in a 1:1 solution of DMF:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

### Description

5-hydroxy Omeprazole is a metabolite of the H<sup>+</sup>/K<sup>+</sup>-ATPase inhibitor omeprazole (Item No. 14880).<sup>1</sup> It is formed from omeprazole by the cytochrome P450 (CYP) isoform CYP2C19.

### Reference

1. Abelö, 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, 02/28/2023

#### 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