

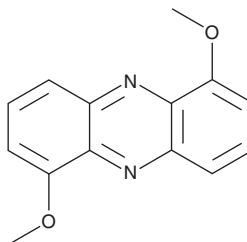
PRODUCT INFORMATION



1,6-Dimethoxyphenazine

Item No. 28061

CAS Registry No.: 13398-79-3
Formal Name: 1,6-dimethoxy-phenazine
Synonym: Crystalloiodinine B
MF: C₁₄H₁₂N₂O₂
FW: 240.3
Purity: ≥95%
Supplied as: A solid
Storage: -20°C
Stability: ≥4 years
Item Origin: Bacterium/*Streptomyces* sp.



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Laboratory Procedures

1,6-Dimethoxyphenazine is supplied as a solid. A stock solution may be made by dissolving the 1,6-dimethoxyphenazine in the solvent of choice, which should be purged with an inert gas. 1,6-Dimethoxyphenazine is soluble in ethanol, methanol, DMSO, and dimethyl formamide.

1,6-Dimethoxyphenazine is sparingly soluble in aqueous solutions. To enhance aqueous solubility, dilute the organic solvent solution into aqueous buffers or isotonic saline. If performing biological experiments, ensure the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. We do not recommend storing the aqueous solution for more than one day.

Description

1,6-Dimethoxyphenazine is a bacterial metabolite that has been found in *S. thioluteus*.¹ It is active against the fungus *T. mentagrophytes* (MIC = 25 µg/ml) and the bacteria *M. smegmatis*, *M. rhodochrous*, *N. asteroides*, and *N. coeliaca* (MICs = 25, 5, 5, and 20 µg/ml, respectively), as well as *Actinoplanes* species (MIC = 25 µg/ml).

Reference

- Gerber, N.N. Phenazines, phenoxazinones, and dioxopiperazines from *Streptomyces thioluteus*. *J. Org. Chem.* **32**(12), 4055-4057 (1967).

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, 10/20/2022

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