

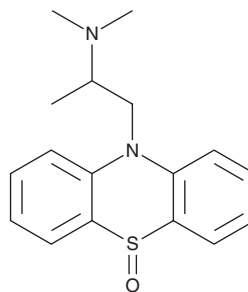
PRODUCT INFORMATION



Promethazine Sulfoxide

Item No. 39547

CAS Registry No.: 7640-51-9
Formal Name: N,N,α-trimethyl-10H-phenothiazine-10-ethanamine, 5-oxide
MF: C₁₇H₂₀N₂OS
FW: 300.4
Purity: ≥95%
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

Promethazine sulfoxide is supplied as a solid. A stock solution may be made by dissolving the promethazine sulfoxide in the solvent of choice, which should be purged with an inert gas. Promethazine sulfoxide is soluble in organic solvents such as methanol and chloroform.

Description

Promethazine sulfoxide is a metabolite of the histamine H₁ receptor antagonist promethazine (Item No. 16478).¹ It is formed from promethazine by the cytochrome P450 (CYP) isoform CYP2D6.²

References

1. Taylor, G. and Houston, J.B. Simultaneous determination of promethazine and two of its circulating metabolites by high-performance liquid chromatography. *J. Chromatogr.* **230(1)**, 194-198 (1982).
2. Nakamura, K., Yokoi, T., Inoue, K., et al. CYP2D6 is the principal cytochrome P450 responsible for metabolism of the histamine H1 antagonist promethazine in human liver microsomes. *Pharmacogenetics* **6(5)**, 449-457 (1996).

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, 10/25/2023

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