

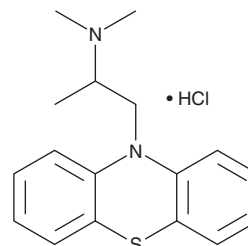
PRODUCT INFORMATION



Promethazine (hydrochloride)

Item No. 22450

CAS Registry No.: 58-33-3
Formal Name: N,N,α-trimethyl-10H-phenothiazine-10-ethanamine, monohydrochloride
Synonym: NSC 231688
MF: C₁₇H₂₀N₂S • HCl
FW: 320.9
Purity: ≥98%
Supplied as: A neat solid
Storage: -20°C
Stability: ≥1 year



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

Description

Promethazine (Item No. 22450) is an analytical reference material categorized as a phenothiazine.¹ Promethazine has been associated with misuse in conjunction with opioids or sedatives.^{1,2} Formulations containing promethazine have been used in the treatment of various allergy-related conditions, anesthesia- or surgery-related nausea and vomiting, and to induce sedation. This product is intended for analytical forensic applications. This product is also available as a general research tool (Item No. 16478).

This product is qualified as a Reference Material that has been manufactured and tested to ISO/IEC 17025 and ISO 17034 international standards.

References

1. Tsay, M.E., Procopio, G., Anderson, B.D., *et al.* Abuse and intentional misuse of promethazine reported to US Poison Centers: 2002 to 2012. *J. Addict. Med.* **9(3)**, 233-237 (2015).
2. Chiappini, S., Schifano, F., Corkery, J.M., *et al.* Beyond the 'purple drank': Study of promethazine abuse according to the European Medicines Agency adverse drug reaction reports. *J. Psychopharmacol.* **35(6)**, 681-692 (2021).

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, 02/27/2024

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