

PRODUCT INFORMATION



N-demethyl Promethazine (hydrochloride)

Item No. 39870

CAS Registry No.: 60113-77-1
Formal Name: N, α -dimethyl-10H-phenothiazine-10-ethanamine, monohydrochloride
Synonyms: DMPZ, Norpromethazine, (\pm)-N-demethyl Promethazine, *rac*-N-demethyl Promethazine, Promethazine EP Impurity C, N-desmethyl Promethazine

MF: C₁₆H₁₈N₂S • HCl

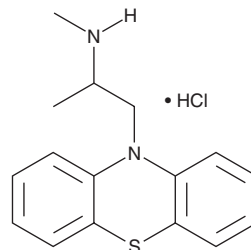
FW: 306.9

Purity: \geq 98%

Supplied as: A solid

Storage: -20°C

Stability: \geq 4 years



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

Description

N-demethyl Promethazine (hydrochloride) (Item No. 39870) is an analytical reference standard categorized as a phenothiazine metabolite.¹ N-demethyl Promethazine is a metabolite of promethazine (Item Nos. 21331 | 22450). This product is intended for research and forensic applications.

Reference

1. Campbell, C., Cornthwaite, H., and Watterson, J. Oxidation of selected phenothiazine drugs during sample preparation: Effects of varying extraction conditions on the extent of oxidation. *J. Anal. Toxicol.* **42**(2), 99-114 (2018).

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