

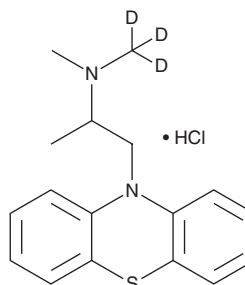
PRODUCT INFORMATION



Promethazine-d₃ (hydrochloride)

Item No. 38299

CAS Registry No.: 1435933-86-0
Formal Name: N,α-dimethyl-N-(methyl-d₃)-10H-phenothiazine-10-ethanamine, monohydrochloride
MF: C₁₇H₁₇D₃N₂S • HCl
FW: 323.9
Chemical Purity: ≥98% (Promethazine)
Deuterium Incorporation: ≥99% deuterated forms (d₁-d₃); ≤1% d₀
Supplied as: A crystalline solid
Storage: -20°C
Stability: ≥4 years



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

Description

Promethazine-d₃ (Item No. 38299) is intended for use as an internal standard for the quantification of promethazine (Item Nos. 21331 | 22450) by GC- or LC-MS. The accuracy of the sample weight in this vial is between 5% over and 2% under the amount shown on the vial. If better precision is required, the deuterated standard should be quantitated against a more precisely weighed unlabeled standard by constructing a standard curve of peak intensity ratios (deuterated *versus* unlabeled).

Promethazine is categorized as a phenothiazine.¹ It 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 research and forensic applications.

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