

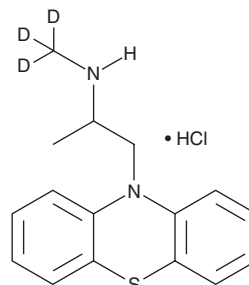
# PRODUCT INFORMATION



## N-demethyl Promethazine-d<sub>3</sub> (hydrochloride)

Item No. 39871

**CAS Registry No.:** 1330173-19-7  
**Formal Name:** N-(methyl-d<sub>3</sub>)-1-(10H-phenothiazin-10-yl)propan-2-amine, monohydrochloride  
**Synonyms:** Norpromethazine-d<sub>3</sub>, (±)-N-demethyl Promethazine-d<sub>3</sub>, rac-N-demethyl Promethazine-d<sub>3</sub>, Promethazine EP Impurity C-d<sub>3</sub>, N-desmethyl Promethazine-d<sub>3</sub>  
**MF:** C<sub>16</sub>H<sub>15</sub>D<sub>3</sub>N<sub>2</sub>S • HCl  
**FW:** 309.9  
**Chemical Purity:** ≥98% (N-demethyl promethazine)  
**Deuterium Incorporation:** ≥99% deuterated forms (d<sub>1</sub>-d<sub>3</sub>); ≤1% d<sub>0</sub>  
**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.

### Description

N-demethyl Promethazine-d<sub>3</sub> (hydrochloride) (Item No. 39871) is intended for use as an internal standard for the quantification of N-demethyl promethazine 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).

N-demethyl Promethazine is a phenothiazine metabolite.<sup>1</sup> It is a metabolite of promethazine (Item Nos. 21331 | 22450). N-demethyl Promethazine is also a degradation product of promethazine.<sup>2</sup> This product is intended for research and forensic applications.

### References

1. Blement, B. and Beckett, A.H. Metabolism of promethazine *in vitro*. Identificaton of N-oxidized products. *Xenobiotica* **11(9)**, 609-618 (1981).
2. Shearer, C.M., and Beckett, S.M. Promethazine hydrochloride. *Analyt. Profiles Drug. Subst.* **5**, 429-465 (1976).

#### 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, 03/08/2024

#### 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