

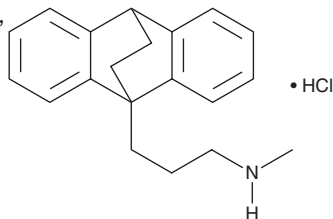
PRODUCT INFORMATION



Maprotiline (hydrochloride)

Item No. 32702

CAS Registry No.: 10347-81-6
Formal Name: N-methyl-9,10-ethanoanthracene-9(10H)-propanamine, monohydrochloride
Synonym: Ba 34276
MF: C₂₀H₂₃N • HCl
FW: 313.9
Purity: ≥98%
Supplied as: A neat solid
Storage: -20°C
Stability: ≥4 years



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

Description

Maprotiline (Item No. 32702) is an analytical reference material categorized as a tricyclic antidepressant.¹ Formulations containing maprotiline have been associated with overdoses.² This product is intended for use in analytical forensic applications. This product is also available as a general research tool (Item No. 15892).

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

References

1. Spencer, P.S. Review of the pharmacology of existing antidepressants. *Br. J. Clin. Pharmacol.* **4(Suppl 2)**, 57S-68S (1977).
2. Keller, T. and Zollinger, U. Gas chromatographic examination of postmortem specimens after maprotiline intoxication. *Forensic Sci. Int.* **88(2)**, 117-123 (1997).

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, 01/06/2021

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