

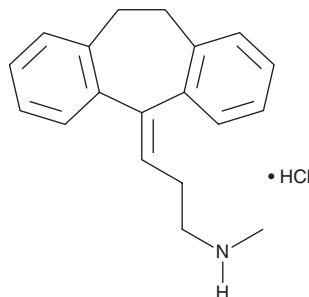
PRODUCT INFORMATION



Nortriptyline (hydrochloride)

Item No. 19041

CAS Registry No.: 894-71-3
Formal Name: 3-(10,11-dihydro-5H-dibenzo[a,d]cyclohepten-5-ylidene)-N-methyl-1-propanamine, monohydrochloride
MF: C₁₉H₂₁N • HCl
FW: 299.8
Purity: ≥98%
Supplied as: A neat solid
Storage: -20°C
Stability: ≥7 years



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

Description

Nortriptyline (hydrochloride) (Item No. 19041) is an analytical reference material categorized as an antidepressant.¹ Nortriptyline is an active metabolite of amitriptyline (Item Nos. 19031 | 19029).² It decreases immobility time in the forced swim test in rats and has antinociceptive activity in mice.^{1,3} Nortriptyline decreases nicotine self-administration in rats.⁴ Formulations containing nortriptyline have been used in the treatment of depression. This product is intended for analytical forensic applications. This product is also available as a general research tool (Item No. 15904).

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. Petersén, A., Wörtwein, G., Gruber, S.H., *et al.* Nortriptyline mediates behavioral effects without affecting hippocampal cytochrome in a genetic rat depression model. *Neurosci. Lett.* **451(2)**, 148-151 (2009).
2. Coutts, R.T., Bach, M.V., and Baker, G.B. Metabolism of amitriptyline with CYP2D6 expressed in a human cell line. *Xenobiotica* **27(1)**, 33-47 (1997).
3. Tavakoli, F., Hejazian, S.H., and Anoosheh, S.M. Antinociceptive activity of nortriptyline through the adrenergic system - an *in vivo* study in mice. *Int. J. Clin. Exp. Physiol.* **5(3)**, 120-122 (2018).
4. Wing, V.C., and Shoaib, M. Translating the smoking cessation properties of the antidepressant nortriptyline using reinforcing, discriminative and aversive stimulus effects of nicotine in rats. *Psychopharmacology (Berl.)* **219**, 847-857 (2012).

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, 11/01/2023

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