

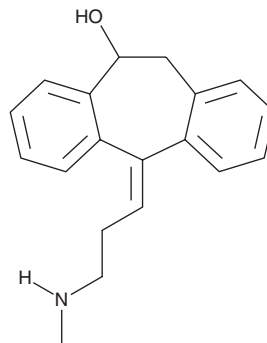
PRODUCT INFORMATION



cis-10-hydroxy Nortriptyline

Item No. 36659

CAS Registry No.: 47132-19-4
Formal Name: (5Z)-10,11-dihydro-5-[3-(methylamino)propylidene]-5H-dibenzo[a,d]cyclohepten-10-ol
Synonym: 10(Z)-hydroxy Nortriptyline
MF: C₁₉H₂₁NO
FW: 279.4
Purity: ≥90%
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.

Laboratory Procedures

cis-10-hydroxy Nortriptyline is supplied as a solid. A stock solution may be made by dissolving the cis-10-hydroxy nortriptyline in the solvent of choice, which should be purged with an inert gas. cis-10-hydroxy Nortriptyline is sparingly soluble (1-10 mg/ml) in chloroform and slightly soluble (0.1-1 mg/ml) in acetonitrile.

Description

cis-10-hydroxy Nortriptyline is an active metabolite of the tricyclic antidepressants amitriptyline (Item No. 15881) and nortriptyline (Item No. 15904).^{1,2} It is formed from amitriptyline and nortriptyline by the cytochrome P450 (CYP) isoform CYP2D6.¹ cis-10-hydroxy Nortriptyline inhibits norepinephrine reuptake in rat cortical slices (EC₅₀ = 160 nM).²

References

1. 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).
2. Bertilsson, L., Mellström, B., and Sjöqvist, F. Pronounced inhibition of noradrenaline uptake by 10-hydroxymetabolites of nortriptyline. *Life Sci.* **25(15)**, 1285-1292 (1979).

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, 01/24/2025

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