

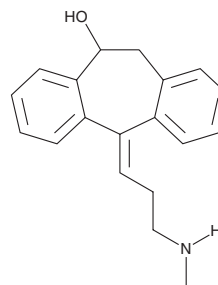
# PRODUCT INFORMATION



## *trans*-10-hydroxy Nortriptyline

Item No. 36660

**CAS Registry No.:** 47132-16-1  
**Formal Name:** (5E)-10,11-dihydro-5-[3-(methylamino)propylidene]-5H-dibenzo[a,d]cyclohepten-10-ol  
**Synonym:** 10(E)-hydroxy Nortriptyline  
**MF:** C<sub>19</sub>H<sub>21</sub>NO  
**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

*trans*-10-hydroxy Nortriptyline is supplied as a solid. A stock solution may be made by dissolving the *trans*-10-hydroxy nortriptyline in the solvent of choice, which should be purged with an inert gas. *trans*-10-Hydroxy nortriptyline is soluble in organic solvents such as ethanol and DMSO. The solubility of *trans*-10-hydroxy nortriptyline in these solvents is approximately 3 and 1 mg/ml, respectively. *trans*-10-hydroxy Nortriptyline is slightly soluble in dimethyl formamide.

### Description

*trans*-10-hydroxy Nortriptyline is an active metabolite of the tricyclic antidepressants amitriptyline (Item No. 15881) and nortriptyline (Item No. 15904).<sup>1,2</sup> It is formed from amitriptyline and nortriptyline by the cytochrome P450 (CYP) isoform CYP2D6.<sup>1</sup> *trans*-10-hydroxy Nortriptyline inhibits norepinephrine reuptake in rat cortical slices and isolated human plasma (EC<sub>50</sub>s = 160 and 176 nM, respectively).<sup>2</sup>

### 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, I., 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, 12/20/2022

#### 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