

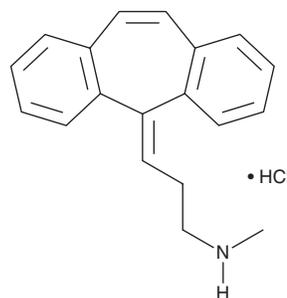
PRODUCT INFORMATION



N-Desmethylocyclobenzaprine (hydrochloride)

Item No. 15906

CAS Registry No.: 438-59-5
Formal Name: 3-(5H-dibenzo[a,d]cyclohepten-5-ylidene)-N-methyl-1-propanamine, monohydrochloride
Synonyms: 10,11-Dehydronortriptyline, N 7068, Ro 4-6011
MF: C₁₉H₁₉N • HCl
FW: 297.8
Purity: ≥98%
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-Desmethylocyclobenzaprine (hydrochloride) (Item No. 15906) is an analytical reference standard categorized as an antidepressant.¹ N-Desmethylocyclobenzaprine is an active metabolite of amitriptyline (Item Nos. 19031 | 19029) and cyclobenzaprine (Item Nos. ISO60170 | 15885).¹⁻³ This product is intended for research and forensic applications.

References

1. Petersen, P.V., Lassen, N., Hansen, V., *et al.* Pharmacological studies of a new series of bicyclic thymoleptics. *Acta Pharmacol. Toxicol. (Copenh)* **24(2)**, 121-133 (1966).
2. Gupta, R.N., Molnar, G., Hill, R.E., *et al.* Estimation of amitriptyline and its metabolites in serum and urine by GLC using nitrogen-specific detector. *Clin. Biochem.* **9(5)**, 247-251 (1976).
3. Belvedere, G., Rovei, V., Pantarotto, C., *et al.* Identification of cyclobenzaprine-10, 11-epoxide and other metabolites after incubation of cyclobenzaprine with rat liver microsomes. *Xenobiotica* **5(12)**, 765-771 (2008).

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, 05/23/2025

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