

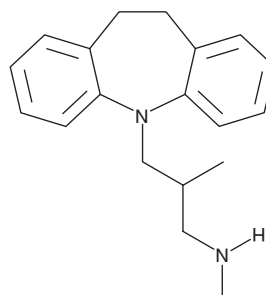
PRODUCT INFORMATION



N-desmethyl Trimipramine

Item No. 34982

CAS Registry No.: 2293-21-2
Formal Name: 10,11-dihydro-N, β -dimethyl-5H-dibenz[b,f]azepine-5-propanamine
Synonym: Monodesmethyltrimipramine, Nortrimipramine
MF: C₁₉H₂₄N₂
FW: 280.4
Purity: $\geq 95\%$
UV/Vis.: λ_{max} : 250 nm
Supplied as: A solution in methanol
Storage: -20°C
Stability: ≥ 2 years



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

Laboratory Procedures

N-desmethyl Trimipramine is supplied as a solution in methanol. To change the solvent, simply evaporate the N-desmethyl trimipramine under a gentle stream of nitrogen and immediately add the solvent of choice. Solvents such as chloroform can be used. N-desmethyl Trimipramine is slightly soluble in chloroform.

Description

N-desmethyl Trimipramine is an active metabolite of the tricyclic antidepressant trimipramine (Item No. 15921).¹ It inhibits the norepinephrine transporter (NET), serotonin (5-HT) transporter (SERT), and dopamine transporter (DAT; IC₅₀s = 2.92, 6.22, and 29.5 μM , respectively) as well as organic cation transporter 1 (OCT1) and OCT2 (IC₅₀s = 3.19 and 10.4 μM , respectively) in HEK293 cells expressing the human transporters.

Reference

1. Haenisch, B., Hiemke, C., and Bönisch, H. Inhibitory potencies of trimipramine and its main metabolites at human monoamine and organic cation transporters. *Psychopharmacology (Berl)* **217(2)**, 289-295 (2011).

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, 12/08/2022

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