

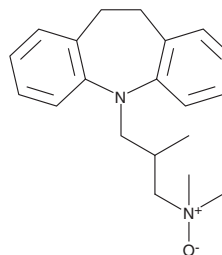
PRODUCT INFORMATION



Trimipramine N-oxide

Item No. 35414

CAS Registry No.: 14171-70-1
Formal Name: 10,11-dihydro-N,N,β-trimethyl-5H-dibenz[b,f]azepine-5-propanamine, N-oxide
MF: C₂₀H₂₆N₂O
FW: 310.4
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

Trimipramine N-oxide is supplied as a solid. A stock solution may be made by dissolving the trimipramine N-oxide in the solvent of choice, which should be purged with an inert gas. Trimipramine N-oxide is soluble in DMSO.

Description

Trimipramine N-oxide is an active metabolite of the tricyclic antidepressant trimipramine (Item No. 15921).¹ It is formed via N-oxidation of trimipramine by the cytochrome P450 (CYP) isoforms CYP2D6, CYP2C9, CYP2C19, CYP1A2, and CYP3A4. Trimipramine N-oxide selectively inhibits the serotonin transporter (SERT) over organic cation transporter 1 (OCT1) and OCT2, the dopamine transporter (DAT), and the norepinephrine transporter (NET) in HEK293 cells expressing the human transporters (IC₅₀s = 3.59, 9.35, 27.4, 9.4, and 11.7 μM, respectively).

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, 11/19/2024

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