

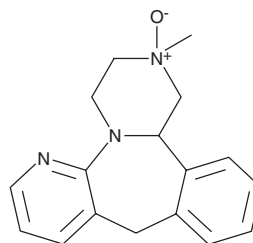
PRODUCT INFORMATION



Mirtazapine N-oxide

Item No. 34214

CAS Registry No.: 155172-12-6
Formal Name: 1,2,3,4,10,14b-hexahydro-2-methyl-pyrazino[2,1-a]pyrido[2,3-c][2]benzazepine, 2-oxide
MF: $C_{17}H_{19}N_3O$
FW: 281.4
Purity: $\geq 95\%$ (mixture of diastereomers)
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

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

Description

Mirtazapine N-oxide is a metabolite of mirtazapine.¹ It is formed from mirtazapine by the cytochrome P450 (CYP) isoforms CYP1A2 and CYP3A4 in human liver microsomes.

Reference

1. Störmer, E., von Moltke, L.L., Shader, R.I., *et al.* Metabolism of the antidepressant mirtazapine in vitro: Contribution of cytochromes P-450 1A2, 2D6, and 3A4. *Drug Metab. Dispos.* **28(10)**, 1168-1175 (2000).

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, 11/18/2024

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