

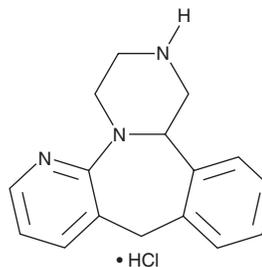
PRODUCT INFORMATION



Desmethyl Mirtazapine (hydrochloride)

Item No. 34261

CAS Registry No.: 1188265-41-9
Formal Name: 1,2,3,4,10,14b-hexahydro-pyrazino[2,1-a]pyrido[2,3-c][2]benzazepine, monohydrochloride
Synonyms: N-Desmethyl Mirtazapine, Normirtazapine
MF: C₁₆H₁₇N₃ • HCl
FW: 287.8
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.

Laboratory Procedures

Desmethyl mirtazapine (hydrochloride) is supplied as a solid. A stock solution may be made by dissolving the desmethyl mirtazapine (hydrochloride) in the solvent of choice, which should be purged with an inert gas. Desmethyl mirtazapine (hydrochloride) is soluble in the organic solvent DMSO.

Description

Desmethyl mirtazapine is a metabolite of the antidepressant mirtazapine.¹ It is formed from mirtazapine by the cytochrome P450 (CYP) isoform CYP3A4.

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, 11/07/2022

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