

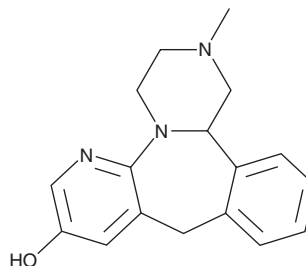
PRODUCT INFORMATION



8-hydroxy Mirtazapine

Item No. 35443

CAS Registry No.: 102335-57-9
Formal Name: 1,2,3,4,10,14b-hexahydro-2-methyl-pyrazino[2,1-a]pyrido[2,3-c][2]benzazepin-8-ol
Synonym: 8-hydroxy-6-Azamianserin
MF: C₁₇H₁₉N₃O
FW: 281.4
Purity: ≥95%
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

8-hydroxy Mirtazapine is supplied as a solid. A stock solution may be made by dissolving the 8-hydroxy mirtazapine in the solvent of choice, which should be purged with an inert gas. 8-hydroxy Mirtazapine is slightly soluble in DMSO and methanol (sonicated).

Description

8-hydroxy Mirtazapine is a metabolite of the antidepressant mirtazapine.¹ It is formed from mirtazapine by the cytochrome P450 (CYP) isoform CYP2D6.

Reference

1. Dahl, M.-L., Voortman, G., Alm, C., *et al.* *In vitro* and *in vivo* studies on the disposition of mirtazapine in humans. *Clin. Drug Investig.* **13**(1), 37-46 (1997).

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, 11/18/2022

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