

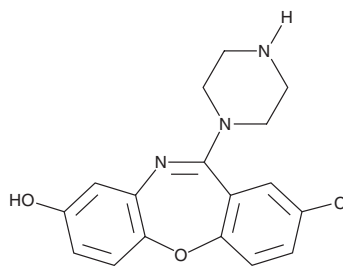
PRODUCT INFORMATION



8-hydroxy Amoxapine

Item No. 29192

CAS Registry No.: 61443-78-5
Formal Name: 2-chloro-11-(1-piperazinyl)-dibenz[b,f][1,4]oxazepin-8-ol
MF: C₁₇H₁₆ClN₃O₂
FW: 329.8
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 Amoxapine is supplied as a solid. A stock solution may be made by dissolving the 8-hydroxy amoxapine in the solvent of choice, which should be purged with an inert gas. 8-hydroxy Amoxapine is slightly soluble in DMSO.

Description

8-hydroxy Amoxapine is a metabolite of the tetracyclic antidepressant amoxapine (Item No. 17347).^{1,2}

References

1. Johnson, S.M., Nygard, G., and Khalil, S.K. Isocratic liquid chromatographic method for the determination of amoxapine and its metabolites. *J. Pharm. Sci.* **73(5)**, 696-699 (1984).
2. Tasset, J.J. and Pesce, A.J. Amoxapine in human overdose. *J. Anal. Toxicol.* **8(3)**, 124-128 (1984).

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, 10/03/2022

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