

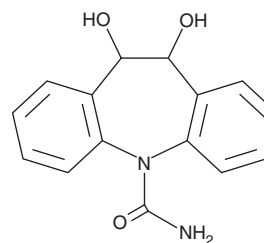
PRODUCT INFORMATION



10,11-dihydro-10,11-dihydroxy Carbamazepine

Item No. 35436

CAS Registry No.: 35079-97-1
Formal Name: 10,11-dihydro-10,11-dihydroxy-5H-dibenz[b,f]azepine-5-carboxamide
Synonyms: 10,11-dihydro-10,11-dihydroxy CBZ, CBZ-diol, Dihydroxycarbamazepine, Dihydroxycarbamazepine
MF: C₁₅H₁₄N₂O₃
FW: 270.3
Purity: ≥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

10,11-dihydro-10,11-dihydroxy Carbamazepine is supplied as a solid. A stock solution may be made by dissolving the 10,11-dihydro-10,11-dihydroxy carbamazepine in the solvent of choice, which should be purged with an inert gas. 10,11-dihydro-10,11-dihydroxy Carbamazepine is soluble in DMSO and methanol.

Description

10,11-dihydro-10,11-dihydroxy Carbamazepine is a metabolite of the anticonvulsant carbamazepine and the antiepileptic prodrug oxcarbazepine (Item No. 17340).¹ It is formed from carbamazepine via a carbamazepine 10,11-epoxide (Item No. 33908) intermediate by epoxide hydrolase and from oxcarbazepine via a 10,11-dihydro-10-hydroxy carbamazepine (Item No. 18467) intermediate.^{1,2} 10,11-dihydro-10,11-dihydroxy Carbamazepine has been found in wastewater effluent.³

References

1. Franceschi, L. and Furlanut, M. A simple method to monitor plasma concentrations of oxcarbazepine, carbamazepine, their main metabolites and lamotrigine in epileptic patients. *Pharmacol. Res.* **51**(4), 297-302 (2005).
2. Tybring, G., van Bahr, C., Bertilsson, L., et al. Metabolism of carbamazepine and its epoxide metabolite in human and rat liver *in vitro*. *Drug. Metab. Dispos.* **9**(6), 561-564 (1981).
3. Miao, X.-S., Yang, J.-J., and Metcalfe, C.D. Carbamazepine and its metabolites in wastewater and in biosolids in a municipal wastewater treatment plant. *Environ. Sci. Technol.* **39**(19), 7469-7475 (2005).

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