

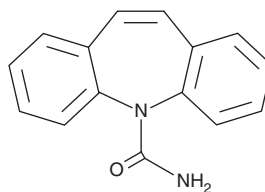
PRODUCT INFORMATION



Carbamazepine

Item No. 15422

CAS Registry No.: 298-46-4
Formal Name: 5H-dibenz[b,f]azepine-5-carboxamide
Synonyms: CBZ, NSC 169864
MF: C₁₅H₁₂N₂O
FW: 236.3
Purity: ≥98%
UV/Vis.: λ_{max}: 214, 238, 285 nm
Supplied as: A crystalline 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

Carbamazepine is supplied as a crystalline solid. A stock solution may be made by dissolving the carbamazepine in the solvent of choice, which should be purged with an inert gas. Carbamazepine is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of carbamazepine in ethanol is approximately 3 mg/ml and approximately DMSO in DMF and 25 mg/ml.

Carbamazepine is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, carbamazepine should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. Carbamazepine has a solubility of approximately 0.5 mg/ml in a 1:1 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

Carbamazepine is an inhibitor of sodium channels.¹ It preferentially inhibits inactivated voltage-gated sodium channel Na_v1.7 and Na_v1.3 (IC₅₀s = 406 and 900 μM, respectively) over resting (closed) Na_v1.7 and Na_v1.3 channels in whole-cell patch-clamp assays using HEK293 cells (IC₅₀s = 1,584 and 2,464 μM, respectively). Carbamazepine (100 μM) inhibits sodium currents in a rat sciatic nerve patch-clamp assay *ex vivo*.² It reduces levels of the endogenous catecholamine dopamine in the striatum and hippocampus, as well as the levels of the dopamine precursor L-DOPA (Item No. 13248) in the striatum, in rats when administered at a dose of 100 mg/kg.³ Dietary administration of carbamazepine (0.5% w/w) reduces amphetamine-induced hyperactivity in mice.⁴ It decreases immobility time in the forced swim test in mice when administered in the diet at 0.75% w/w. Formulations containing carbamazepine have been used in the treatment of seizures, trigeminal neuralgia, and bipolar disorder.

References

1. Sheets, P.L., Heers, C., Stoehr, T., *et al.* Differential block of sensory neuronal voltage-gated sodium channels by lacosamide [(2R)-2-(acetylamino)-N-benzyl-3-methoxypropanamide], lidocaine, and carbamazepine. *J. Pharmacol. Exp. Ther.* **326**(1), 89-99 (2008).
2. Schwarz, J.R. and Grigat, G. Phenytoin and carbamazepine: Potential- and frequency-dependent block of Na currents in mammalian myelinated nerve fibers. *Epilepsia* **30**(3), 286-294 (1989).
3. Okada, M., Hirano, T., Mizuno, K., *et al.* Biphasic effects of carbamazepine on the dopaminergic system in rat striatum and hippocampus. *Epilepsy Res.* **28**(2), 143-153 (1997).
4. Kara, N.Z., Karpel, O., Toker, L., *et al.* Chronic oral carbamazepine treatment elicits mood-stabilising effects in mice. *Acta Neuropsychiatr.* **26**(1), 29-34 (2014).

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, 08/28/2023

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