

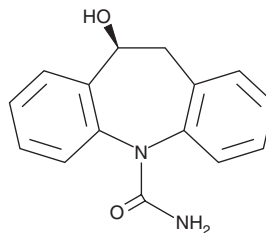
PRODUCT INFORMATION



Eslicarbazepine

Item No. 35008

CAS Registry No.: 104746-04-5
Formal Name: (10S)-10,11-dihydro-10-hydroxy-5H-dibenz[b,f]azepine-5-carboxamide
Synonyms: BIA 2-194, CGP 13751, (S)-Licarbazepine
MF: C₁₅H₁₄N₂O₂
FW: 254.3
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

Eslicarbazepine is supplied as a solid. A stock solution may be made by dissolving the eslicarbazepine in the solvent of choice, which should be purged with an inert gas. Eslicarbazepine is soluble in DMSO.

Description

Eslicarbazepine is a voltage-gated sodium channel (Na_v) inhibitor and an active metabolite of the prodrug eslicarbazepine acetate (Item No. 23656).^{1,2} It enhances slow activation of voltage-gated sodium channels, reducing channel availability.¹ Eslicarbazepine is protective against seizures induced by maximal electroshock (MES) in mice (ED₅₀ = 17.1 mg/kg).

References

- Soares-Da-Silva, P., Pires, N., Bonifácio, M.J., *et al.* Eslicarbazepine acetate for the treatment of focal epilepsy: An update on its proposed mechanisms of action. *Pharmacol. Res. Perspect.* **3**(2), e00124 (2015).
- Benes, J., Parada, A., Figueiredo, A.A., *et al.* Anticonvulsant and sodium channel-blocking properties of novel 10,11-dihydro-5H-dibenz[b,f]azepine-5-carboxamide derivatives. *J. Med. Chem.* **42**(14), 2582-2587 (1999).

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, 08/04/2023

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