

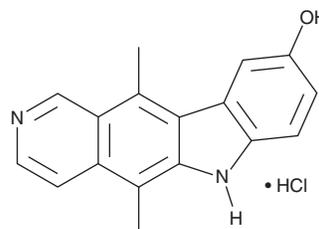
PRODUCT INFORMATION



9-Hydroxyellipticine (hydrochloride)

Item No. 17921

CAS Registry No.: 52238-35-4
Formal Name: 5,11-dimethyl-6H-pyrido[4,3-b]carbazol-9-ol, monohydrochloride
Synonyms: 9-OH-Ellipticine, NSC 210717
MF: C₁₇H₁₄N₂O • HCl
FW: 298.8
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

9-Hydroxyellipticine (hydrochloride) is supplied as a solid. A stock solution may be made by dissolving the 9-hydroxyellipticine (hydrochloride) in the solvent of choice, which should be purged with an inert gas. 9-Hydroxyellipticine (hydrochloride) is slightly soluble in DMSO (warmed) and methanol.

Description

9-Hydroxyellipticine is a derivative of ellipticine (Item No. 18742) with diverse biological activities.¹⁻³ It inhibits aroclor-induced activation of aniline hydroxylase, aminopyrine N-demethylase, and 7-ethoxycoumarin O-deethylase in rat liver microsomes (K_i s = 3.5, 0.6, and 0.74 μ M, respectively).¹ 9-Hydroxyellipticine inhibits the growth of L1210 murine leukemia cells (IC_{50} = 3 nM) *in vitro* and increases survival in an L1210 mouse leukemia model.^{2,3} It inhibits carrageenan-induced edema and UV-induced erythema in guinea pigs.⁴

References

1. Lesca, P., Rafidinarivo, E., Lecointe, P., *et al.* A class of strong inhibitors of microsomal monooxygenases: The ellipticines. *Chem. Biol. Interact.* **24(2)**, 189-197 (1979).
2. Monnot, M., Mauffret, O., Simon, V., *et al.* DNA-drug recognition and effects on topoisomerase II-mediated cytotoxicity. A three-mode binding model for ellipticine derivatives. *J. Biol. Chem.* **266(3)**, 1820-1829 (1991).
3. Le Pecq, J.B., Grosse, C., Dat-Xuong, N., *et al.* Antitumor activity of 9-hydroxyellipticine (NSC 210717) ON L1210 mouse leukemia and the effect of route of injection. *Cancer Res.* **36(9 pt. 1)**, 3067-3076 (1976).
4. Cros, J., Thibault, A., and Dat-Xuong, N. Anti-inflammatory effect of hydroxy-9-ellipticine. *C. R. Acad. Sci. Hebd. Seances Acad. Sci. D.* **281(15)**, 1139-1142 (1975).

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, 10/24/2022

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