

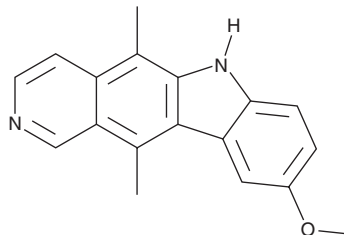
PRODUCT INFORMATION



9-Methoxyellipticine

Item No. 31631

CAS Registry No.: 10371-86-5
Formal Name: 9-methoxy-5,11-dimethyl-6H-pyrido[4,3-b]carbazole
Synonyms: Methoxy-9-ellipticine, NSC 69187
MF: C₁₈H₁₆N₂O
FW: 276.3
Purity: ≥95%
UV/Vis.: λ_{max}: 245, 276, 293 nm
Supplied as: A solid
Storage: -20°C
Stability: ≥4 years
Item Origin: Animal/*Bleekeria* sp.



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Laboratory Procedures

9-Methoxyellipticine is supplied as a solid. A stock solution may be made by dissolving the 9-methoxyellipticine in the solvent of choice, which should be purged with an inert gas. 9-Methoxyellipticine is soluble in DMSO.

Description

9-Methoxyellipticine is an alkaloid and a derivative of the DNA intercalating agent ellipticine (Item No. 18742) that has been found in *O. maculata* and has anticancer activity.^{1,2} It intercalates into DNA and is an inhibitor of wild-type c-Kit and c-Kit containing the D816V activating mutation (c-Kit^{D816V}; IC₅₀s = 0.8 and 0.6 μM, respectively).^{2,3} 9-Methoxyellipticine inhibits proliferation of Ba/F3 cells in the presence of the c-Kit ligand SCF and Ba/F3 cells expressing c-Kit^{D816V} (IC₅₀s = 0.5 and 0.3 μM, respectively).² It also increases survival in a variety of rodent tumor models.¹

References

1. Svoboda, G.H., Poore, G.A., and Montfort, M.L. Alkaloids of *Ochrosia maculata* Jacq. (*Ochrosia borbonica* Gmel.). Isolation of the alkaloids and study of the antitumor properties of 9-methoxyellipticine. *J. Pharm. Sci.* **57**(10), 1720-1725 (1968).
2. Vendôme, J., Letard, S., Martin, F., et al. Molecular modeling of wild-type and D816V c-Kit inhibition based on ATP-competitive binding of ellipticine derivatives to tyrosine kinases. *J. Med. Chem.* **48**(20), 6194-6201 (2005).
3. Festy, B., Poisson, J., and Paoletti, C. A new DNA intercalating drug: Methoxy-9-ellipticine. *FEBS Lett.* **17**(2), 321-323 (1971).

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, 11/14/2022

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