

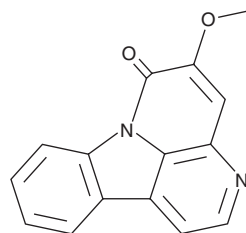
PRODUCT INFORMATION



5-Methoxycanthinone

Item No. 31342

CAS Registry No.: 15071-56-4
Formal Name: 5-methoxy-6H-indolo[3,2,1-de][1,5]naphthyridin-6-one
Synonyms: 5-Methoxycanthin-6-one, NSC 88929
MF: C₁₅H₁₀N₂O₂
FW: 250.3
Purity: ≥95%
Supplied as: A solid
Storage: -20°C
Stability: ≥4 years
Item Origin: Plant/*Quassia amara*



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

Laboratory Procedures

5-Methoxycanthinone is supplied as a solid. A stock solution may be made by dissolving the 5-methoxycanthinone in the solvent of choice, which should be purged with an inert gas. 5-Methoxycanthinone is soluble in organic solvents such as DMSO and methanol.

Description

5-Methoxycanthinone is an alkaloid that has been found in *L. floridana* and has anticancer properties.¹ It is cytotoxic to KB, A549, and HCT-8 cells (EC₅₀s = 2.9, 3.1, and 2.5 µg/ml, respectively). It is active against the *P. falciparum* mefloquine-resistant strain F32 and the multidrug and chloroquine-resistant strain K1 (IC₅₀s = 10.4 and 5.1 µg/ml, respectively).² However, it increases the parasite load by 68.4% in a mouse model of leishmaniasis when administered at a dose of 10 mg/kg per day.³

References

- Xu, Z., Chang, F.R., Wang, H.-K., et al. Anti-HIV agents 45 and antitumor agents 205. Two new sesquiterpenes, leitneridanins A and B, and the cytotoxic and anti-HIV principles from *Leitneria floridana*. *J. Nat. Prod.* **63**(12), 1712-1715 (2000).
- Cebrián-Torrejón, G., Spelman, K., Leblanc, K., et al. The antiplasmodium effects of a traditional South American remedy: *Zanthoxylum chiloperone* var. *Angustifolium* against chloroquine resistant and chloroquine sensitive strains of *Plasmodium falciparum*. *Revista Brasileira de Farmacognosia* **21**(4), 652-661 (2011).
- Ferreira, M.E., Rojas de Arias, A., Torres de Ortiz, S., et al. Leishmanicidal activity of two canthin-6-one alkaloids, two major constituents of *Zanthoxylum chiloperone* var. *angustifolium*. *J. Ethnopharmacol.* **80**(2-3), 199-202 (2002).

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, 08/18/2023

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