

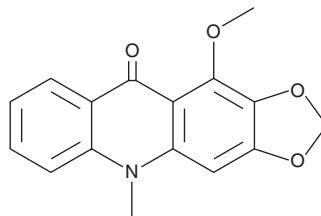
PRODUCT INFORMATION



Evoxanthine

Item No. 29437

CAS Registry No.: 477-82-7
Formal Name: 11-methoxy-5-methyl-1,3-dioxolo[4,5-b]acridin-10(5H)-one
Synonym: NSC 407812
MF: C₁₆H₁₃NO₄
FW: 283.3
Purity: ≥95%
Supplied as: A solid
Storage: -20°C
Stability: ≥4 years
Item Origin: Plant/*Evodia alata*



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

Laboratory Procedures

Evoxanthine is supplied as a solid. A stock solution may be made by dissolving the evoxanthine in the solvent of choice, which should be purged with an inert gas. Evoxanthine is soluble in the organic solvents ethanol, methanol, DMSO, and dimethyl formamide.

Description

Evoxanthine is an alkaloid that has been found in *O. renieri* and has antimalarial and anticancer activities.^{1,2} It is active against *P. falciparum* with an IC₅₀ value of 67.6 µg/ml.¹ Evoxanthine decreases proliferation of nine sensitive and drug-resistant cancer cell lines (IC₅₀s = 6.11-80.99 µM).²

References

1. Khalid, S.A., Farouk, A., Geary, T.G., *et al.* Potential antimalarial candidates from African plants: And *in vitro* approach using *Plasmodium falciparum*. *J. Ethnopharmacol.* **15(2)**, 201-209 (1986).
2. Kuete, V., Fouotsa, H., Mbaveng, A.T., *et al.* Cytotoxicity of a naturally occurring furoquinoline alkaloid and four acridone alkaloids towards multi-factorial drug-resistant cancer cells. *Phytomedicine* **22(10)**, 946-951 (2015).

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, 10/17/2022

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