

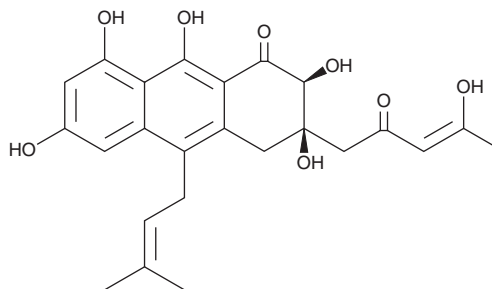
PRODUCT INFORMATION



Melicopine

Item No. 31636

CAS Registry No.: 568-01-4
Formal Name: 4,5-dimethoxy-11-methyl-1,3-dioxolo[4,5-c]acridin-6(11H)-one
MF: C₁₇H₁₅NO₅
FW: 313.3
Purity: ≥95%
Supplied as: A solid
Storage: -20°C
Stability: ≥4 years
Item Origin: Plant/*Melicope* sp.



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

Laboratory Procedures

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

Description

Melicopine is an acridone alkaloid that has been found in *Z. simulans* and has antimalarial and anticancer activities.¹ It is active against the chloroquine-sensitive 3D7 and chloroquine-resistant Dd2 strains of *P. falciparum* with IC₅₀ values of 29.7 and 33.7 µg/ml, respectively. Melicopine is cytotoxic to PC-3M and LNCaP prostate cancer, but not non-cancerous HEK293, cells (IC₅₀s = 47.9, 37.8, and >100 µg/ml, respectively).

Reference

1. Wang, C., Wan, J., Mei, Z., et al. Acridone alkaloids with cytotoxic and antimalarial activities from *Zanthoxylum simulans* Hance. *Pharmacogn. Mag.* **10(37)**, 73-76 (2014).

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, 11/10/2022

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