

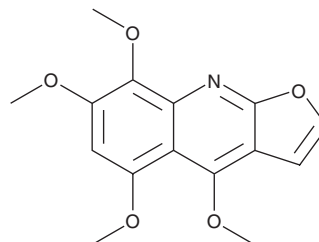
PRODUCT INFORMATION



5,7,8-Trimethoxydictamnine

Item No. 31418

CAS Registry No.: 521-43-7
Formal Name: 4,5,7,8-tetramethoxy-furo[2,3-b]quinoline
Synonyms: Acronycidine, NSC 30619
MF: C₁₅H₁₅NO₅
FW: 289.3
Purity: ≥95%
Supplied as: A powder
Storage: -20°C
Stability: ≥4 years
Item Origin: Plant/*Halfordia kendack*



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

Laboratory Procedures

5,7,8-Trimethoxydictamnine is supplied as a powder. A stock solution may be made by dissolving the 5,7,8-trimethoxydictamnine in the solvent of choice, which should be purged with an inert gas. 5,7,8-Trimethoxydictamnine is soluble in methanol and DMSO.

Description

5,7,8-Trimethoxydictamnine is a quinoline alkaloid that has been found in *Rutaceae* and has antimalarial activity.¹ It is active against chloroquine-sensitive and -resistant strains of *P. falciparum* (IC₅₀s =19.9 and 5.72 µg/ml, respectively).

Reference

1. Basco, L.K.M., S., Skaltsounis, A.-L., Ravvelomanantsoa, N., *et al.* In vitro activities of furoquinoline and acridone alkaloids against *Plasmodium falciparum*. *Antimicrob. Agents Chemother.* **38(5)**, 1169-1171 (1994).

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, 11/09/2022

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