

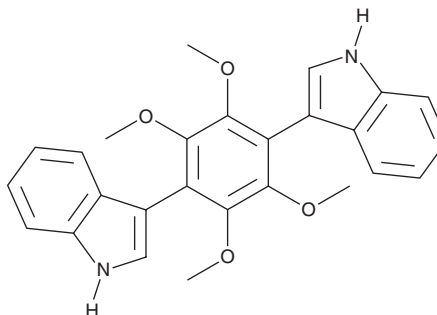
PRODUCT INFORMATION



Asterriquinol D dimethyl ether

Item No. 25080

CAS Registry No.: 287117-66-2
Formal Name: 3,3'-(2,3,5,6-tetramethoxy-1,4-phenylene)bis-1H-indole
MF: C₂₆H₂₄N₂O₄
FW: 428.5
Purity: ≥95%
Supplied as: A solid
Storage: -20°C
Stability: ≥4 years
Item Origin: Fungus/*Penicillium citreonigrum*



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

Laboratory Procedures

Asterriquinol D dimethyl ether is supplied as a solid. A stock solution may be made by dissolving the asterriquinol D dimethyl ether in the solvent of choice, which should be purged with an inert gas. Asterriquinol D dimethyl ether is soluble in DMSO.

Description

Asterriquinol D dimethyl ether is a fungal metabolite that has been found in *A. kumbius*.¹ It is cytotoxic to NS-1 mouse myeloma cells (IC₅₀ = 28 µg/ml) and has activity against *T. foetus in vitro* (IC₅₀ = 100 µg/ml).

Reference

1. Lacey, H.J., Vuong, D., Pitt, J.I., *et al.* Kumbicins A–D: Bis-Indolyl benzenoids and benzoquinones from an Australian soil fungus, *Aspergillus kumbius*. *Aust. J. Chem.* **69(2)**, 152-160 (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, 08/23/2023

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