

PRODUCT INFORMATION



19,20-Epoxyctochalasin C

Item No. 28025

CAS Registry No.: 189351-79-9
Formal Name: (3S,3aR,4E,7S,9R,9aR,10aS,11S,11aR,14S,14aR)-11-(acetyloxy)-6,7,9,9a,10a,11,14,14a-octahydro-3,9-dihydroxy-1,2,7,9-tetramethyl-14-(phenylmethyl)-3H-oxireno[9,10]cycloundec[1,2-d]isoindole-8,12(3aH,13H)-dione

MF: C₃₀H₃₇NO₇

FW: 523.6

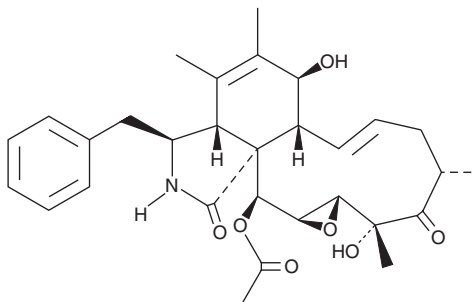
Purity: ≥95%

Supplied as: A powder

Storage: -20°C

Stability: ≥4 years

Item Origin: Fungus/*Geniculosporium* sp.



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

Laboratory Procedures

19,20-Epoxyctochalasin C is supplied as a powder. A stock solution may be made by dissolving the 19,20-epoxyctochalasin C in the solvent of choice, which should be purged with an inert gas. 19,20-Epoxyctochalasin C is soluble in organic solvents such as ethanol, methanol, DMSO, and dimethyl formamide.

Description

19,20-Epoxyctochalasin C is a fungal metabolite originally isolated from *Nemania* sp. UM10M.¹ It inhibits the growth of chloroquine-sensitive and -resistant strains of *P. falciparum* (IC₅₀s = 0.07 and 0.05 ng/ml, respectively) without inducing cytotoxicity in Vero cells (IC₅₀ = >4,760 ng/ml). 19,20-Epoxyctochalasin C is also phytotoxic, reducing growth of lettuce and bentgrass plants when used at a concentration of 1 mg/ml.

Reference

1. Kumarihamy, M., Ferreira, D., Croom, E.M., Jr., *et al.* Antiplasmodial and cytotoxic cytochalasins from an endophytic fungus, *Nemania* sp. UM10M, isolated from a diseased *Torreya taxifolia* leaf. *Molecules* **24**(4), E777 (2019).

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, 08/10/2023

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