

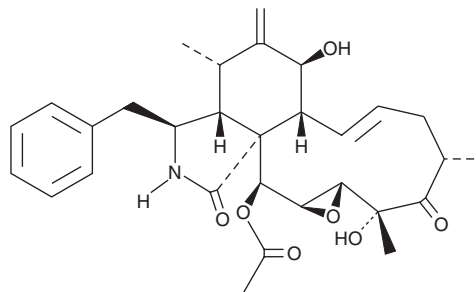
# PRODUCT INFORMATION



## 19,20-Epoxyctochalasin D

Item No. 28050

**CAS Registry No.:** 191349-10-7  
**Formal Name:** (1S,3S,3aR,4E,7S,9R,9aR,10aS,11S,11aR,14S,14aR)-11-(acetyloxy)-3,3a,6,7,9,9a,10a,11,14,14a-decahydro-3,9-dihydroxy-1,7,9-trimethyl-2-methylene-14-(phenylmethyl)-1H-oxireno[9,10]cycloundec[1,2-d]isoindole-8,12(2H,13H)-dione  
**MF:** C<sub>30</sub>H<sub>37</sub>NO<sub>7</sub>  
**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 D is supplied as a powder. A stock solution may be made by dissolving the 19,20-epoxyctochalasin D in the solvent of choice, which should be purged with an inert gas. 19,20-Epoxyctochalasin D is soluble in organic solvents such as ethanol, methanol, DMSO, and dimethyl formamide.

### Description

19,20-Epoxyctochalasin D is a fungal metabolite that has been found in the endophytic fungus *Nemania* sp. UM10M.<sup>1</sup> It is active against the chloroquine-sensitive and -resistant strains of *P. falciparum* (MIC = 0.4 ng/ml for both) without inducing cytotoxicity in Vero cells. It is cytotoxic to BT-549, LLC-PK11, and P388 cells (IC<sub>50</sub>s = 7.84, 8.4, and 0.16 μM, respectively) but not SK-MEL, KB, or SKOV3 cells up to a concentration of 10 μM.<sup>1,2</sup>

### References

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).
2. Shi, L.-M. and Zhan, Z.-J. Structural revision of 19, 20-epoxyctochalasin D and its cytotoxic activity. *J. Chem. Res.* **2007**(3), 144-145 (2007).

#### 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