

# PRODUCT INFORMATION



## Previridicatumtoxin

Item No. 29069

**CAS Registry No.:** 1379585-81-5  
**Formal Name:** (4aS,5S,12aS)-6-[(2E)-3,7-dimethyl-2,6-octadien-1-yl]-1,4,4a,5,12,12a-hexahydro-3,4a,5,10,11,12a-hexahydroxy-8-methoxy-1,12-dioxo-2-naphthacene-carboxamide

**MF:** C<sub>30</sub>H<sub>33</sub>NO<sub>10</sub>

**FW:** 567.6

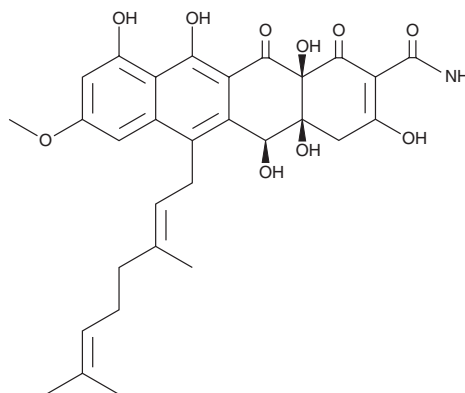
**Purity:** ≥70%

**Supplied as:** A solid

**Storage:** -20°C

**Stability:** ≥4 years

**Item Origin:** Fungus/*Emmonsia* sp.



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

### Laboratory Procedures

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

### Description

Previridicatumtoxin is a fungal metabolite that has been found in *P. aethiopicum* and has diverse biological activities.<sup>1,2</sup> It is an intermediate in the biosynthesis of the mycotoxin viridicatumtoxin (Item No. 25480). Previridicatumtoxin is active against methicillin-resistant *S. aureus* (MRSA) and vancomycin-resistant *E. faecalis* (IC<sub>50</sub>s = 4.4 and 4.8 μM, respectively), as well as *C. albicans* and *S. cerevisiae* (MIC = 32 μg/ml for both).<sup>1,2</sup> It is cytotoxic to NCI H460, KB-3-1, and SW620 cancer cells (IC<sub>50</sub>s = 5.3, 4.1, and 6 μM, respectively).<sup>2</sup>

### References

1. Chooi, Y.H., Wang, P., Fang, J., *et al.* Discovery and characterization of a group of fungal polycyclic polyketide prenyltransferases. *J. Am. Chem. Soc.* **134**(22), 9428-9437 (2012).
2. Shang, Z., Salim, A.A., Khalil, Z., *et al.* Viridicatumtoxins: Expanding on a rare tetracycline antibiotic scaffold. *J. Org. Chem.* **80**(24), 12501-12508 (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, 08/16/2023

#### 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