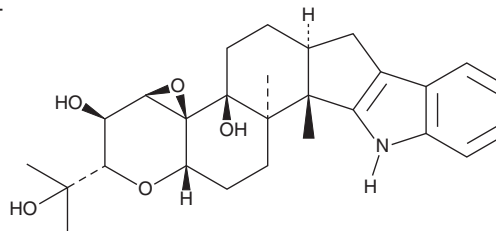


# PRODUCT INFORMATION



## Terpendole I Item No. 29443

**CAS Registry No.:** 167612-17-1  
**Formal Name:** (2S,3R,3aR,4aS,4bS,6aS,12bS,12cR,14aS)-3,3a,5,6,6a,7,12,12b,12c,13,14,14a-dodecahydro-2-(1-hydroxy-1-methylethyl)-12b,12c-dimethyl-2H,4bH-oxireno[4',4'a]-1-benzopyrano[5',6':6,7]indeno[1,2-b]indole-3,4b-diol  
**MF:** C<sub>27</sub>H<sub>35</sub>NO<sub>5</sub>  
**FW:** 453.6  
**Purity:** ≥95%  
**Supplied as:** A solid  
**Storage:** -20°C  
**Stability:** ≥4 years  
**Item Origin:** Fungus/Unidentified sp.



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

### Laboratory Procedures

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

### Description

Terpendole I is a fungal metabolite that has been found in *A. yamanashiensis*.<sup>1</sup> It is a weak inhibitor of acyl-coenzyme A:cholesterol acyltransferase (ACAT; IC<sub>50</sub> = 145 μM) and is active against the bacteria *B. cereus* and *B. subtilis* (MICs = 100 μg/ml for both) but not *S. aureus*, *P. aeruginosa*, or *K. pneumoniae* (MICs = >200 μg/ml for all) or the fungus *C. albicans* (MIC = 200 μg/ml).<sup>1,2</sup> It is cytotoxic to HeLa cells with an IC<sub>50</sub> value of 52.6 μM.<sup>3</sup>

### References

1. Tomoda, H., Tabata, N., Yang, D.-J., *et al.* Terpendoles, novel ACAT inhibitors produced by *Albophoma yamanashiensis*. III. Production, isolation and structure elucidation of new components. *J. Antibiot. (Tokyo)* **48**(8), 793-804 (1995).
2. Zhao, J.-C., Wang, Y.-L., Zhang, T.-Y., *et al.* Indole diterpenoids from the endophytic fungus *Drechmeria* sp. as natural antimicrobial agents. *Phytochemistry* **148**, 21-28 (2018).
3. Nagumo, Y., Motoyama, T., Hayashi, T., *et al.* Structure-activity relationships of terpendole E and its natural derivatives. *ChemistrySelect* **2**(4), 1533-1536 (2017).

#### 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](http://WWW.CAYMANCHEM.COM)