

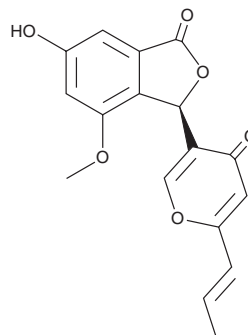
PRODUCT INFORMATION



6-O-Demethylvermistatin

Item No. 44763

CAS Registry No.: 1048672-98-5
Formal Name: (3R)-6-hydroxy-4-methoxy-3-[4-oxo-6-(1E)-1-propen-1-yl-4H-pyran-3-yl]-1(3H)-isobenzofuranone
Synonym: 6-Desmethylvermistatin
MF: C₁₇H₁₄O₆
FW: 314.3
Purity: ≥95%
Supplied as: A solid
Storage: -20°C
Stability: ≥4 years
Item Origin: Fungi/*Talaromyces* sp.



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

Laboratory Procedures

6-O-Demethylvermistatin is supplied as a solid. A stock solution may be made by dissolving the 6-O-demethylvermistatin in the solvent of choice, which should be purged with an inert gas. 6-O-Demethylvermistatin is soluble in methanol and DMSO.

Description

6-O-Demethylvermistatin is a fungal metabolite that has been found in *Penicillium*.¹

Reference

1. Liu, Y., Xia, G., Li, H., et al. Vermistatin derivatives with α -glucosidase inhibitory activity from the mangrove endophytic fungus *Penicillium* sp. HN29-3B1. *Planta Med.* **80**(11), 912-917 (2014).

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, 11/20/2025

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