

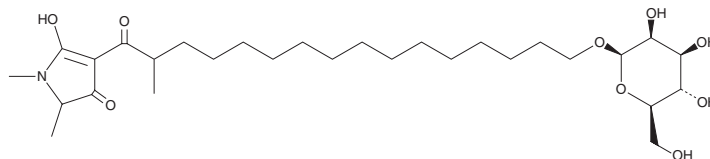
PRODUCT INFORMATION



Epicoccamide

Item No. 24176

CAS Registry No.: 606139-26-8
Formal Name: 1,2-dihydro-5-hydroxy-4-[16-(β-D-mannopyranosyloxy)-2-methyl-1-oxohexadecyl]-1,2-dimethyl-3H-pyrrol-3-one
MF: C₂₉H₅₁NO₉
FW: 557.7
Purity: ≥98%
Supplied as: A residue
Storage: -20°C
Stability: ≥4 years



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

Laboratory Procedures

Epicoccamide is supplied as a residue. A stock solution may be made by dissolving the epicoccamide in the solvent of choice, which should be purged with an inert gas. Epicoccamide is soluble in organic solvents such as ethanol, methanol, DMSO, and dimethyl formamide.

Epicoccamide is sparingly soluble in aqueous solutions. To enhance aqueous solubility, dilute the organic solvent solution into aqueous buffers or isotonic saline. If performing biological experiments, ensure the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. We do not recommend storing the aqueous solution for more than one day.

Description

Epicoccamide is a marine fungal metabolite originally isolated from *E. purpurascens*.¹ It is comprised of three subunits including a glycosidic subunit, a fatty acid, and a tetramic acid. Epicoccamide has no activity against various bacteria, fungi, and nematodes and is not cytotoxic.^{1,2}

References

1. Wright, A.D., Osterhage, C., and König, G.M. Epicoccamide, a novel secondary metabolite from a jellyfish-derived culture of *Epicoccum purpurascens*. *Org. Biomol. Chem.* **1(3)**, 507-510 (2003).
2. Wangun, H.V., Dahse, H.-M., and Hertweck, C. Epicoccamides B-D, glycosylated tetramic acid derivatives from an *Epicoccum* sp. associated with the tree fungus *Pholiota squarrosa*. *J. Nat. Prod.* **70(11)**, 1800-1803 (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, 12/17/2022

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