

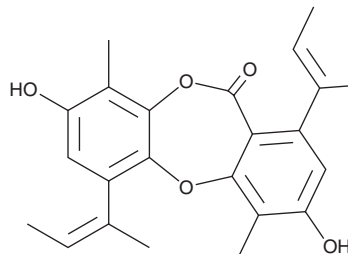
PRODUCT INFORMATION



Folipastatin

Item No. 24924

CAS Registry No.: 139959-71-0
Formal Name: 3,8-dihydroxy-4,9-dimethyl-1,6-bis[(1Z)-1-methyl-1-propen-1-yl]-11H-dibenzo[b,e][1,4]dioxepin-11-one
MF: C₂₃H₂₄O₅
FW: 380.4
Purity: ≥95%
Supplied as: A solid
Storage: -20°C
Stability: ≥4 years
Item Origin: Fungus/*Aspergillus unguis*



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

Laboratory Procedures

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

Description

Folipastatin is a fungal metabolite produced by *A. unguis* and an inhibitor of phospholipase A₂ (PLA₂; IC₅₀ = 39 μM).¹ It inhibits release of arachidonic acid (Item Nos. 90010 | 10006607) from rat polymorphonuclear leukocytes (IC₅₀ = 24 μM). Folipastatin also inhibits radioligand binding to L-type calcium channels in porcine heart membranes (IC₅₀ = 0.39 μM) but has no effect on contraction of taenia (IC₅₀ = >100 μM).

Reference

1. Hamano, K., Kinoshita-Okami, M., Hemmi, A., *et al.* Folipastatin, a new depsidone compound from *Aspergillus unguis* as an inhibitor of phospholipase A₂. Taxonomy, fermentation, isolation, structure determination and biological properties. *J. Antibiot. (Tokyo)* **45(8)**, 1195-1201 (1992).

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, 08/23/2023

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