

PRODUCT INFORMATION



Roridin L2

Item No. 27374

CAS Registry No.: 85124-22-7
Formal Name: 12,13-epoxy-trichothec-9-ene-4,15-diol, 4β-[(2Z,4E)-6-[2-(2,5-dihydro-5-oxo-3-furanyl)ethoxy]-7-hydroxy-2,4-octadienoate]

MF: C₂₉H₃₈O₉

FW: 530.6

Purity: ≥70%

UV/Vis.: λ_{max}: 260 nm

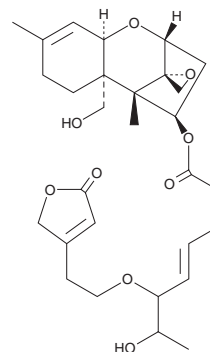
Supplied as: An oil

Storage: -20°C

Stability: ≥4 years

Item Origin: Fungus/Unidentified marine fungus

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



Laboratory Procedures

Roridin L2 is supplied as an oil. A stock solution may be made by dissolving the roridin L2 in the solvent of choice, which should be purged with an inert gas. Roridin L2 is soluble in ethanol, methanol, DMSO, and dichloromethane.

Description

Roridin L2 is a fungal metabolite that has been found in *M. roridum*.¹

Reference

1. Grove, J.F. The trichothecenes and their biosynthesis. *Fortschr. Chem. Org. Naturst.* **88**, 63-130 (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, 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