

PRODUCT INFORMATION



Chrodrimanin H

Item No. 44759

CAS Registry No.: 1446801-59-7
Formal Name: 1,8S,9,9aR,10,12,13,13aS,13bS,14-decahydro-5,8-dihydroxy-2R,7aR,10,10,13a-pentamethyl-2H,4H-benzo[a]pyrano[3,4-j]xanthene-4,11(7aH)-dione

MF: C₂₅H₃₂O₆

FW: 428.5

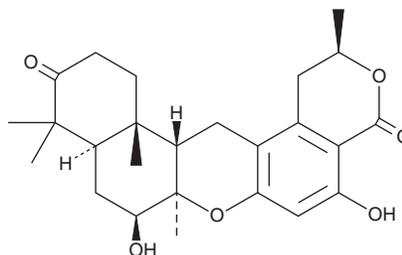
Purity: ≥95%

Supplied as: A solid

Storage: -20°C

Stability: ≥4 years

Item Origin: Fungi/*Penicillium* sp.



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

Laboratory Procedures

Chrodrimanin H is supplied as a solid. A stock solution may be made by dissolving the chrodrimanin H in the solvent of choice, which should be purged with an inert gas. Chrodrimanin H is soluble in methanol and DMSO.

Description

Chrodrimanin H is a meroterpenoid that has been isolated from *T. primulinus* and has anti-inflammatory activity.¹ It inhibits LPS-induced increases in nitric oxide (NO) production in RAW 264.7 macrophages when used at a concentration of 50 μM.

Reference

1. Chen, J.Q., Li, S., Fan, R.Z., *et al.* Talaesthanes A-C, three new meroterpenoids from the endophytic fungus *Talaromyces primulinus* H21. *Fitoterapia* **177**, 106085 (2024).

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