

PRODUCT INFORMATION



Chrodrimanin B

Item No. 27585

CAS Registry No.: 132196-54-4
Formal Name: (1R,2R,7aR,8S,9aR,13aS,13bS)-1-(acetyloxy)-1,8,9,9a,10,13a,13b,14-octahydro-5,8-dihydroxy-2,7a,10,10,13a-pentamethyl-2H,4H-benzo[a]pyrano[3,4-j]xanthene-4,11(7aH)-dione

Synonym: Thailandolide B

MF: C₂₇H₃₂O₈

FW: 484.5

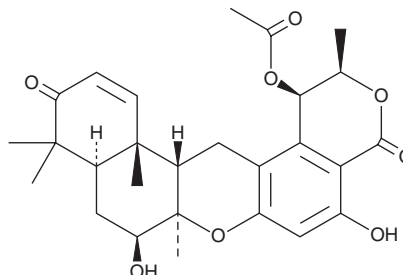
Purity: ≥95%

Supplied as: A solid

Storage: -20°C

Stability: ≥4 years

Item Origin: Fungus/*Penicillium* sp.



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

Laboratory Procedures

Chrodrimanin B is supplied as a solid. A stock solution may be made by dissolving the chrodrimanin B in the solvent of choice, which should be purged with an inert gas. Chrodrimanin B is soluble in organic solvents such as ethanol, methanol, and DMSO. The solubility of chrodrimanin B in ethanol is approximately 1 mg/ml.

Description

Chrodrimanin B is a meroterpenoid fungal metabolite that has been found in *Talaromyces* species.¹ It has insecticidal activity against third instar silkworm larvae (LD₅₀ = 10 µg/g of diet). Chrodrimanin B selectively inhibits the silkworm GABA receptor RDL (IC₅₀ = 1.13 nM) over human α1β2γ2 subunit-containing GABA receptors (IC₅₀ = 1.48 µM).² The structure of thailandolide B, originally described as a C-9 epimer of chrodrimanin B, was revised in 2012 and is identical to that of chrodrimanin B.¹

References

- Hayashi, H., Oka, Y., Kai, K., *et al.* A new meroterpenoid, chrodrimanin C, from YO-2 of *Talaromyces* sp. *Biosci. Biotech. Biochem.* **76**(4), 745-748 (2012).
- Xu, Y., Furutani, S., Ihara, M., *et al.* Meroterpenoid chrodrimanins are selective and potent blockers of insect GABA-gated chloride channels. *PLoS One* **10**(4), e0122629 (2015).

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, 08/09/2023

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