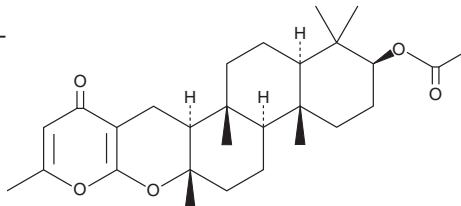


PRODUCT INFORMATION



Chevalone C Item No. 25084

CAS Registry No.: 1318025-77-2
Formal Name: (2S,4aR,4bR,6aS,12aS,12bR,14aR)-2-(acetyloxy)-2,3,4,4a,4b,5,6,6a,12,12a,12b,13,14,14a-tetradecahydro-1,1,4a,6a,9,12b-hexamethyl-1H,11H-phenanthro[2,1-b]pyrano[3,2-e]pyran-11-one
MF: C₂₈H₄₀O₅
FW: 456.6
Purity: ≥95%
Supplied as: A solid
Storage: -20°C
Stability: ≥4 years



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

Laboratory Procedures

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

Description

Chevalone C is a meroterpenoid fungal metabolite originally isolated from *E. chevalieri*.¹ It is active against *M. tuberculosis* H37Ra (MIC = 6.3 µg/ml) and is cytotoxic to BC1 human breast cancer cells (IC₅₀ = 8.7 µg/ml). Chevalone C inhibits the growth of multidrug-resistant isolates of *E. coli*, *S. aureus*, and *E. faecium* in a disc diffusion assay when used at a concentration of 15 µg/disc.² It also induces cell death in HCT116 colorectal carcinoma cells.³

References

1. Kanokmedhakul, K., Kanokmedhakul, S., Suwannatrai, R., *et al.* Bioactive meroterpenoids and alkaloids from the fungus *Eurotium chevalieri*. *Tetrahedron* **67(30)**, 5461-5468 (2011).
2. Gomes, N.M., Bessa, L.J., Buttachon, S., *et al.* Antibacterial and antibiofilm activities of tryptoquivalines and meroditerpenes isolated from the marine-derived fungi *Neosartorya paulistensis*, *N. laciniosa*, *N. tsunodae*, and the soil fungi *N. fischeri* and *N. siamensis*. *Mar. Drugs* **12(2)**, 822-839 (2014).
3. Prata-Sena, M., Ramos, A.A., Buttachon, S., *et al.* Cytotoxic activity of secondary metabolites from marine-derived fungus *Neosartorya siamensis* in human cancer cells. *Phytother. Res.* **30(11)**, 1862-1871 (2016).

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, 12/19/2022

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