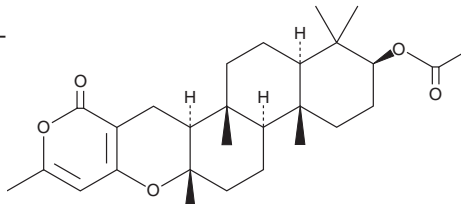


PRODUCT INFORMATION



Chevalone B Item No. 25083

CAS Registry No.: 1318025-75-0
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[3,4-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 B is supplied as a solid. A stock solution may be made by dissolving the chevalone B in the solvent of choice, which should be purged with an inert gas. Chevalone B is slightly soluble (0.1-1 mg/ml) in acetonitrile.

Description

Chevalone B is a meroterpenoid originally isolated from the fungus *E. chevalieri*.¹ It is cytotoxic to KB and NCI-H187 cells with IC₅₀ values of 2.9 and 9.8 µg/ml, respectively.

Reference

1. Kanokmedhakul, K., Kanokmedhakul, S., Suwannatrai, R., *et al.* Bioactive meroterpenoids and alkaloids from the fungus *Eurotium chevalieri*. *Tetrahedron* **67**(30), 5461-5468 (2011).

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, 01/23/2025

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