

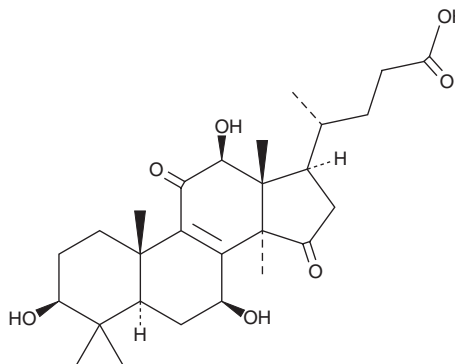
PRODUCT INFORMATION



Lucidenic Acid C

Item No. 43509

CAS Registry No.: 95311-96-9
Formal Name: (3 β ,5 α ,7 β ,12 β)-3,7,12-trihydroxy-4,4,14-trimethyl-11,15-dioxo-chol-8-en-24-oic acid
MF: C₂₇H₄₀O₇
FW: 476.6
Purity: $\geq 98\%$
Supplied as: A solid
Storage: -20°C
Stability: ≥ 4 years
Item Origin: Fungus/*Ganoderma* sp.



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

Laboratory Procedures

Lucidenic acid C is supplied as a solid. A stock solution may be made by dissolving the lucidenic acid C in the solvent of choice, which should be purged with an inert gas. Lucidenic acid C is slightly soluble (0.1-1 mg/ml) in DMSO and acetonitrile.

Description

Lucidenic acid C is a triterpene that has been found in *G. lucidum* and has anticancer activity.¹ It inhibits invasion of HepG2 cells induced by phorbol 12-myristate 13-acetate (PMA; Item No. 10008014) in a migration assay when used at a concentration of 50 μ M.

Reference

1. Weng, C.-J., Chau, C.-F., Chen, K.-D., *et al.* The anti-invasive effect of lucidenic acids isolated from a new *Ganoderma lucidum* strain. *Mol. Nutr. Food Res.* **51**(12), 1472-1477 (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, 06/25/2025

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