

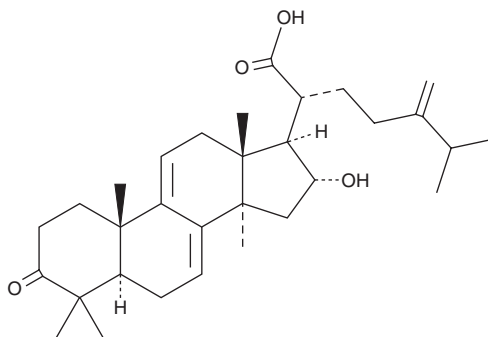
PRODUCT INFORMATION



Polyporenic Acid C

Item No. 44639

CAS Registry No.: 465-18-9
Formal Name: 16 α -hydroxy-24-methylene-3-oxo-lanosta-7,9(11)-dien-21-oic acid
MF: C₃₁H₄₆O₄
FW: 482.7
Purity: $\geq 98\%$
Supplied as: A solid
Storage: -20°C
Stability: ≥ 4 years
Item Origin: Fungus/*Wolfiporia cocos* (Schw.) Ryv.



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

Laboratory Procedures

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

Description

Polyporenic acid C is a triterpene that has been found in *P. cocos* and has diverse biological activities.¹⁻³ It is active against *M. phlei* and *M. smegmatis* when used at concentrations ranging from 0.78 to 2 $\mu\text{g/ml}$.¹ Polyporenic acid C (20, 60, and 80 μM) induces apoptosis in A549 lung cancer cells.² *In vivo*, polyporenic acid C (25 and 50 mg/kg) reduces hepatic lipid accumulation and inguinal white adipose tissue mitochondrial degeneration and increases the inguinal brown adipose tissue index in a mouse model of high-fat diet-induced steatosis.³

References

1. Marcus, S. Antibacterial activity of the triterpenoid acid (polyporenic acid C) and of ungulinic acid, metabolic products of *Polyporus benzoinus* (Wahl.) *Fr. Biochem J.* **50**(4), 516-517 (1952).
2. Ling, H., Zhou, L., Jia, X., *et al.* Polyporenic acid C induces caspase-8-mediated apoptosis in human lung cancer A549 cells. *Mol. Carcinog.* **48**(6), 498-507 (2009).
3. He, X., Tao, Y., Song, C., *et al.* Novel Parkin agonists from *Poria cocos* against dyslipidemia. *Fitoterapia* **182**, 106469 (2025).

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, 12/05/2025

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