

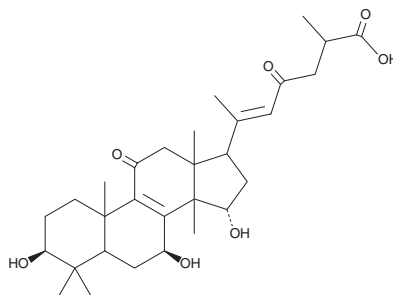
PRODUCT INFORMATION



Ganoderenic Acid C

Item No. 30366

CAS Registry No.: 100665-42-7
Formal Name: (3 β ,7 β ,15 α ,20E)-3,7,15-trihydroxy-11,23-dioxo-lanosta-8,20(22)-dien-26-oic acid
MF: C₃₀H₄₄O₇
FW: 516.7
Purity: $\geq 95\%$
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

Ganoderenic acid C is supplied as a solid. A stock solution may be made by dissolving the ganoderenic acid C in the solvent of choice, which should be purged with an inert gas. Ganoderenic acid C is soluble in methanol.

Description

Ganoderenic acid C is a triterpene that has been found in *G. lucidum*.¹ It is cytotoxic to H460 cancer cells (IC₅₀ = 93 μ M).

Reference

1. Komoda, Y., Nakamura, H., Ishihara, S., *et al.* Structures of new terpenoid constituents of *Ganoderma lucidum* (Fr.) KARST (Polyporaceae). *Chem. Pharm. Bull.* **33(11)**, 4829-4835 (1985).

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, 08/30/2023

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