

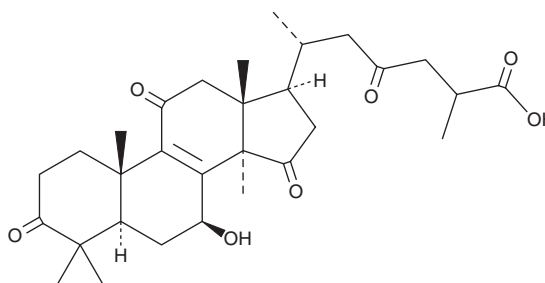
PRODUCT INFORMATION



Ganoderic Acid D

Item No. 26765

CAS Registry No.: 108340-60-9
Formal Name: 7 β -hydroxy-3,11,15,23-tetraoxo-lanost-8-en-26-oic acid
MF: C₃₀H₄₂O₇
FW: 514.7
Purity: $\geq 97\%$
UV/Vis.: λ_{max} : 254 nm
Supplied as: A crystalline 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

Ganoderic acid D is supplied as a crystalline solid. A stock solution may be made by dissolving the ganoderic acid D in the solvent of choice. Ganoderic acid D is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide, which should be purged with an inert gas. The solubility of ganoderic acid D in these solvents is approximately 30 mg/ml.

Ganoderic acid D is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, ganoderic acid D should first be dissolved in ethanol and then diluted with the aqueous buffer of choice. Ganoderic acid D has a solubility of approximately 0.25 mg/ml in a 1:3 solution of ethanol:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

Ganoderic acid D is an oxygenated triterpenoid that has been found in *G. lucidum*.¹ It inhibits glucose uptake and lactate production when used at concentrations ranging from 50 to 200 μM and decreases pyruvate and acetyl-coenzyme A production at concentrations of 100 and 200 μM in HT-29 and SW620 colon cancer cells. Ganoderic acid D increases protein levels of sirtuin 3 (SIRT3) in HT-29 and SW620 cells. It is cytotoxic to HepG2, HeLa, and Caco-2 cells with IC₅₀ values of 0.14, 0.18, and 0.02 $\mu\text{g/ml}$, respectively.²

References

1. Liu, Z., Li, L., and Xue, B. Effect of ganoderic acid D on colon cancer Warburg effect: Role of SIRT3/cyclophilin D. *Eur. J. Pharmacol.* **824**, 72-77 (2018).
2. Ruan, W., Lim, A.H.H., Huang, L.G., et al. Extraction optimisation and isolation of triterpenoids from *Ganoderma lucidum* and their effect on human carcinoma cell growth. *Nat. Prod. Res.* **28(24)**, 2264-2272 (2014).

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, 11/08/2022

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