

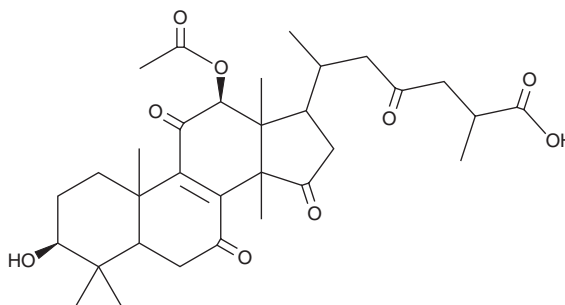
PRODUCT INFORMATION



Ganoderic Acid H

Item No. 29703

CAS Registry No.: 98665-19-1
Formal Name: 12 β -(acetyloxy)-3 β -hydroxy-7,11,15,23-tetraoxo-lanost-8-en-26-oic acid
MF: C₃₂H₄₄O₉
FW: 572.7
Purity: $\geq 95\%$
UV/Vis.: λ_{max} : 255 nm
Supplied as: A crystalline solid
Storage: -20°C
Stability: ≥ 4 years
Item Origin: Fungus/*Ganoderma lucidum*



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

Laboratory Procedures

Ganoderic acid H is supplied as a crystalline solid. A stock solution may be made by dissolving the ganoderic acid H in the solvent of choice, which should be purged with an inert gas. Ganoderic acid H is soluble in the organic solvent DMSO.

Description

Ganoderic acid H is a triterpene that has been found in *G. lucidum* and has diverse biological activities.¹⁻⁴ It inhibits angiotensin-converting enzyme (ACE; IC₅₀ = 26 μM).¹ Ganoderic acid H also inhibits HIV-1 protease with an IC₅₀ value of 200 μM .² It reduces colony formation and migration of MDA-MB-231 breast cancer cells in a concentration-dependent manner.³ *In vivo*, ganoderic acid H (1, 3, and 5 mg/kg) inhibits acetic acid-induced writhing in mice.⁴

References

1. Morigiwa, A., Kitabatake, K., Fujimoto, Y., *et al.* Angiotensin converting enzyme-inhibitory triterpenes from *Ganoderma lucidum*. *Chem. Pharm. Bull. (Tokyo)* **34**(7), 3025-3028 (1986).
2. El-Mekkawy, S., Meselhy, M.R., Nakamura, N., *et al.* Anti-HIV-1 and anti-HIV-1-protease substances from *Ganoderma lucidum*. *Phytochemistry* **49**(6), 1651-1657 (1998).
3. Jiang, J., Grieb, B., Thyagarajan, A., *et al.* Ganoderic acids suppress growth and invasive behavior of breast cancer cells by modulating AP-1 and NF- κ B signaling. *Int. J. Mol. Med.* **21**(5), 577-584 (2008).
4. Koyama, K., Imaizumi, T., Akiba, M., *et al.* Antinociceptive components of *Ganoderma lucidum*. *Planta Med.* **63**(3), 224-227 (1997).

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, 08/16/2023

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