

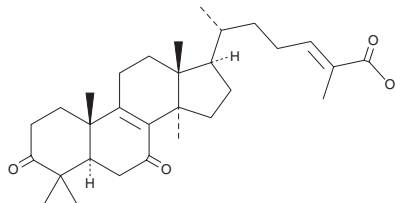
PRODUCT INFORMATION



Ganoderic Acid DM

Item No. 34536

CAS Registry No.: 173075-45-1
Formal Name: 3,7-dioxo-lanosta-8,24E-dien-26-oic acid
MF: $C_{30}H_{44}O_4$
FW: 468.7
Purity: $\geq 98\%$
UV/Vis.: λ_{max} : 217, 251 nm
Supplied as: A solid
Storage: $-20^{\circ}C$
Stability: ≥ 4 years
Item Origin: Plant/Unspecified sp.



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

Laboratory Procedures

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

Description

Ganoderic acid DM is a triterpenoid that has been found in *G. lucidum* and has diverse biological activities.¹⁻³ It inhibits 5 α -reductase and HMG-CoA reductase (IC_{50} s = 10.6 and 9.5 μ M, respectively) and binds to the ligand-binding domain of the androgen receptor in a fluorescence polarization assay.^{1,2} Ganoderic acid suppresses osteoclastogenesis of RAW 264.7 cells when used at a concentration of 12.5 μ M and is cytotoxic to K562 leukemia and PC3 prostate cancer cells (IC_{50} s = 18.8 and 81.6 μ M, respectively). It reduces ear edema induced by phorbol 12-myristate 13-acetate (TPA; Item No. 10008014) in mice (ID_{50} = 0.08 mg/ear).³

References

1. Liu, J., Shiono, J., Shimizu, K., *et al.* Ganoderic acid DM: Anti-androgenic osteoclastogenesis inhibitor. *Bioorg. Med. Chem. Lett.* **19(8)**, 2154-2157 (2009).
2. Wang, K., Bao, L., Xiong, W., *et al.* Lanostane triterpenes from the Tibetan medicinal mushroom *Ganoderma leucocontextum* and their inhibitory effects on HMG-CoA reductase and α glucosidase. *J. Nat. Prod.* **78(8)**, 1977-1989 (2015).
3. Akihisa, T., Nakamura, Y., Tagata, M., *et al.* Anti-inflammatory and anti-tumor-promoting effects of triterpene acids and sterols from the fungus *Ganoderma lucidum*. *Chem. Biodivers.* **4(2)**, 224-231 (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, 12/06/2022

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