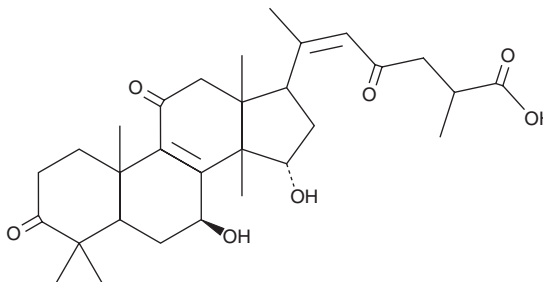


PRODUCT INFORMATION

Ganoderenic Acid A

Item No. 27685

CAS Registry No.: 100665-40-5
 Formal Name: (7 β ,15 α ,20E)-7,15-dihydroxy-3,11,23-trioxo-lanosta-8,20(22)-dien-26-oic acid
 MF: C₃₀H₄₂O₇
 FW: 514.7
 Purity: \geq 95%
 UV/Vis.: λ_{max} : 247 nm
 Supplied as: A crystalline solid
 Storage: -20°C
 Stability: \geq 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

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

Description

Ganoderenic acid A is a triterpenoid originally isolated from *G. lucidum* that has enzyme inhibitory activities.¹⁻³ It is an inhibitor of β -glucuronidase (IC₅₀ = 0.12 mg/ml in rat liver microsomes).² Ganoderenic acid A (100 mg/kg) reduces increases in serum aspartate aminotransferase (AST), alanine aminotransferase (ALT), and lactate dehydrogenase (LDH) in a rat model of liver injury induced by carbon tetrachloride (CCl₄). It also inhibits aldose reductase by 54.2% when used at a concentration of 100 μ g/ml.³

References

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).
2. Kim, D.-H., Shim, S.-B., Kim, N.-J., *et al.* β -Glucuronidase-inhibitory activity and hepatoprotective effect of *Ganoderma lucidum*. *Biol. Pharm. Bull.* **22**(2), 162-164 (1999).
3. Fatmawati, S., Shimizu, K., and Kondo, R. Inhibition of aldose reductase *in vitro* by constituents of *Ganoderma lucidum*. *Planta Med.* **76**(15), 1691-1693 (2010).

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, 04/25/2024

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