

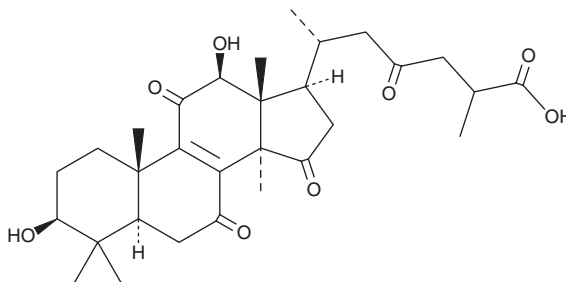
PRODUCT INFORMATION



Ganoderic Acid C6

Item No. 33241

CAS Registry No.: 105742-76-5
Formal Name: 3 β ,12 β -dihydroxy-7,11,15,23-tetraoxo-lanost-8-en-26-oic acid
Synonym: Compound C6
MF: C₃₀H₄₂O₈
FW: 530.7
Purity: \geq 98%
UV/Vis.: λ_{max} : 253 nm
Supplied as: A 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

Ganoderic acid C6 is supplied as a solid. A stock solution may be made by dissolving the ganoderic acid C6 in the solvent of choice, which should be purged with an inert gas. Ganoderic acid C6 is soluble in the organic solvent methanol.

Description

Ganoderic acid C6 is a triterpene that has been found in *G. lucidum*.¹

Reference

1. 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, 08/30/2023

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