

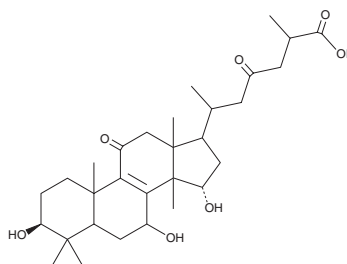
PRODUCT INFORMATION



Ganoderic Acid C2

Item No. 30371

CAS Registry No.: 98296-48-1
Formal Name: 3 β ,7,15 α -trihydroxy-11,23-dioxo-lanost-8-en-26-oic acid
MF: C₃₀H₄₆O₇
FW: 518.7
Purity: \geq 95%
Supplied as: A solid
Storage: -20°C
Stability: \geq 4 years
Item Origin: Fungus/*Ganoderma* sp.



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

Laboratory Procedures

Ganoderic acid C2 is supplied as a solid. A stock solution may be made by dissolving the ganoderic acid C2 in the solvent of choice, which should be purged with an inert gas. Ganoderic acid C2 is soluble in the organic solvent methanol.

Description

Ganoderic acid C2 is a triterpene that has been found in *G. lucidum*.¹ It inhibits histamine release from isolated rat mast cells induced by concanavalin A (Item No. 14951) or compound 48/80 (Item No. 22173) when used at a concentration of 2 mg/ml.

Reference

1. Kohda, H., Tokumoto, W., Sakamoto, K., *et al.* The biologically active constituents of *Ganoderma lucidum* (Fr.) Karst. histamine release-inhibitory triterpenes. *Chem. Pharm. Bull.* **33**(4), 1367-1374 (1985).

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