

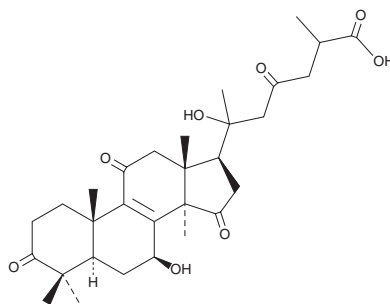
PRODUCT INFORMATION



Ganoderic Acid N

Item No. 34323

CAS Registry No.: 110241-19-5
Formal Name: 7 β ,20 ξ -dihydroxy-3,11,15,23-tetraoxo-lanost-8-en-26-oic acid
MF: C₃₀H₄₂O₈
FW: 530.7
Purity: $\geq$ 85%
UV/Vis.: λ_{max} : 252 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 N is supplied as a solid. A stock solution may be made by dissolving the ganoderic acid N in the solvent of choice, which should be purged with an inert gas. Ganoderic acid N is soluble in DMSO.

Description

Ganoderic acid N is a triterpene that has been found in *G. lingzhi* and has angiotensin-converting enzyme (ACE) inhibitory activity.¹ It inhibits ACE in a cell-free assay (IC₅₀ = 0.941 mM).

Reference

1. Hai-Bang, T. and Shimizu, K. Structure-activity relationship and inhibition pattern of reishi-derived (*Ganoderma lingzhi*) triterpenoids against angiotensin-converting enzyme. *Phytochem. Lett.* **12**, 243-247 (2015).

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, 08/25/2023

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