

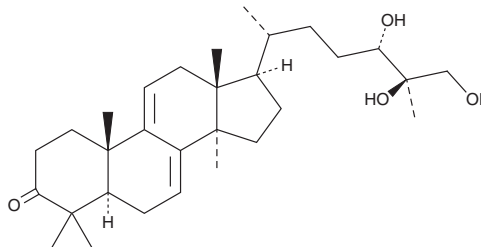
PRODUCT INFORMATION



Ganodermanontriol

Item No. 34325

CAS Registry No.: 106518-63-2
Formal Name: 24S,25R,26-trihydroxy-lanosta-7,9(11)-dien-3-one
MF: $C_{30}H_{48}O_4$
FW: 472.7
Purity: $\geq 90\%$
UV/Vis.: λ_{max} : 244, 253 nm
Supplied as: A solid
Storage: -20°C
Stability: ≥ 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

Ganodermanontriol is supplied as a solid. A stock solution may be made by dissolving the ganodermanontriol in the solvent of choice, which should be purged with an inert gas. Ganodermanontriol is soluble in the organic solvent chloroform.

Description

Ganodermanontriol is a triterpene that has been found in *G. lucidum* and has diverse biological activities.¹⁻⁴ It inhibits the activity of HIV-1 protease ($IC_{50} = 70 \mu\text{M}$), as well as reduces the activity of 5α -reductase by 32% when used at a concentration of $667 \mu\text{M}$.^{1,2} Ganodermanontriol (25 and $50 \mu\text{M}$) inhibits proliferation, colony formation, invasion, and migration of MDA-MB-231 human breast cancer cells.³ It reduces hepatocyte necrosis and increases in hepatic TNF- α and IL-6 levels in a mouse model of LPS/D-galactosamine-induced acute liver injury when administered at a dose of 5 mg/kg.⁴

References

1. Min, B.-S., Nakamura, N., Miyashiro, H., *et al.* Triterpenes from the spores of *Ganoderma lucidum* and their inhibitory activity against HIV-1 protease. *Chem. Pharm. Bull.* **46(10)**, 1607-1612 (1998).
2. Liu, J., Kurashiki, K., Shimizu, K., *et al.* Structure-activity relationship for inhibition of 5α -reductase by triterpenoids isolated from *Ganoderma lucidum*. *Bioorg. Med. Chem.* **14(24)**, 8654-8660 (2006).
3. Jiang, J., Jedinak, A., and Silva, D. Ganodermanontriol (GDNT) exerts its effect on growth and invasiveness of breast cancer cells through the down-regulation of CDC20 and uPA. *Biochem. Biophys. Res. Commun.* **415(2)**, 325-329 (2011).
4. Hu, Z., Du, R., Xiu, L., *et al.* Protective effect of triterpenes of *Ganoderma lucidum* on lipopolysaccharide-induced inflammatory responses and acute liver injury. *Cytokine* **127**, 154917 (2020).

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