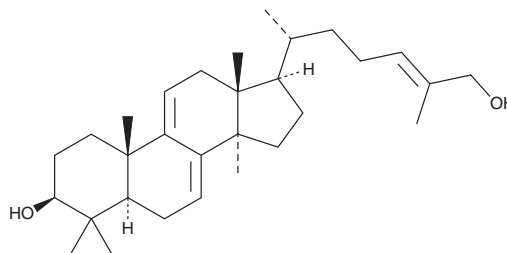


PRODUCT INFORMATION

Ganoderol B

Item No. 34324

CAS Registry No.: 104700-96-1
Formal Name: lanosta-7,9(11),24E-triene-3 β ,26-diol
Synonyms: Ganoderadiol, (+)-Ganoderol B
MF: C₃₀H₄₈O₂
FW: 440.7
Purity: $\geq 95\%$
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

Ganoderol B is supplied as a solid. A stock solution may be made by dissolving the ganoderol B in the solvent of choice, which should be purged with an inert gas. Ganoderol B is soluble in DMSO.

Description

Ganoderol B is a triterpenoid that has been found in *G. lucidum* and has diverse biological activities.¹⁻⁵ It inhibits angiotensin-converting enzyme (ACE) and α -glucosidase (IC₅₀s = 220 and 119.8 μ M, respectively).^{1,2} Ganoderol B is cytotoxic to HGC-27 gastric, HeLa cervical, and A549 lung cancer cells (IC₅₀s = 16.15, 12.64, and 9.77 μ M, respectively).³ It prevents herpes simplex virus 1 (HSV-1) infection in Vero cells (EC₅₀ = 68 μ M).⁴ Ganoderol B (0.001 and 0.1 mg/kg) reduces testosterone-induced increases in ventral prostate weight in castrated rats.⁵

References

1. Morigiwa, A., Kitabatake, K., Fujimoto, Y., *et al.* Angiotensin converting enzyme-inhibitory triterpenes from *Ganoderma lucidum*. *Chem. Pharm. Bull. (Tokyo)* **34**(7), 3025-3028 (1986).
2. Fatmawati, S., Shimizu, K., and Kondo, R. Ganoderol B: A potent α -glucosidase inhibitor isolated from the fruiting body of *Ganoderma lucidum*. *Phytomedicine* **18**(12), 1053-1055 (2011).
3. Su, H.-G., Zhou, Q.-M., Guo, L., *et al.* Lanostane triterpenoids from *Ganoderma luteomarginatum* and their cytotoxicity against four human cancer cell lines. *Phytochemistry* **156**, 89-95 (2018).
4. Mothana, R.A.A., Ali, N.A.A., Jansen, R., *et al.* Antiviral lanostanoid triterpenes from the fungus *Ganoderma pfeifferi*. *Fitoterapia* **74**(1-2), 177-180 (2003).
5. Liu, J., Shimizu, K., Konishi, F., *et al.* The anti-androgen effect of ganoderol B isolated from the fruiting body of *Ganoderma lucidum*. *Bioorg. Med. Chem.* **15**(14), 4966-4972 (2007).

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