

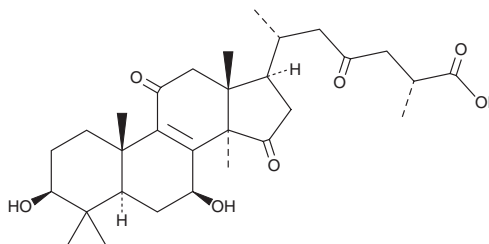
# PRODUCT INFORMATION



## Ganoderic Acid B

Item No. 30112

**CAS Registry No.:** 81907-61-1  
**Formal Name:** (3 $\beta$ ,7 $\beta$ ,25R)-3,7-dihydroxy-11,15,23-trioxo-lanost-8-en-26-oic acid  
**MF:** C<sub>30</sub>H<sub>44</sub>O<sub>7</sub>  
**FW:** 516.7  
**Purity:**  $\geq$ 98%  
**UV/Vis.:**  $\lambda_{\text{max}}$ : 255 nm  
**Supplied as:** A crystalline solid  
**Storage:** -20°C  
**Stability:**  $\geq$ 2 years  
**Item Origin:** Fungi/*Ganoderma lucidum*



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

### Laboratory Procedures

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

### Description

Ganoderic acid B is a triterpenoid that has been found in *G. lucidum* and has diverse biological activities.<sup>1-3</sup> It inhibits HIV-1 protease with an IC<sub>50</sub> value of 0.17 mM.<sup>1</sup> Ganoderic acid B is cytotoxic to HepG2 and Caco-2 cells (IC<sub>50</sub>s = 0.65 and 0.36 mg/ml, respectively).<sup>2</sup> It inhibits acetic acid-induced writhing in mice by 48.3% when administered at a dose of 5 mg/kg.<sup>3</sup>

### References

1. El-Mekawwy, S., Meselhy, M.R., Nakamura, N., *et al.* Anti-HIV-1 and anti-HIV-1-protease substances from *Ganoderma lucidum*. *Phytochemistry* **49(6)**, 1651-1657 (1998).
2. Ruan, W., Lim, A.H.H., Huang, L.G., *et al.* Extraction optimisation and isolation of triterpenoids from *Ganoderma lucidum* and their effect on human carcinoma cell growth. *Nat. Prod. Res.* **28(24)**, 2264-2272 (2014).
3. 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, 05/19/2020

#### 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