

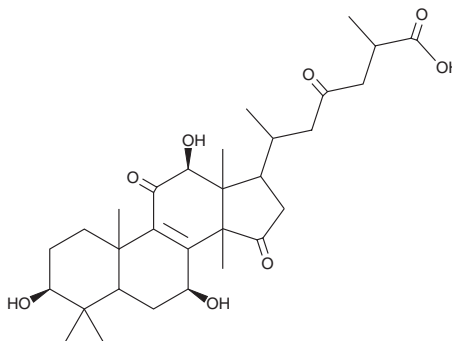
# PRODUCT INFORMATION



## Ganoderic Acid G

Item No. 34675

CAS Registry No.: 98665-22-6  
Formal Name: 3 $\beta$ ,7 $\beta$ ,12 $\beta$ -trihydroxy-11,15,23-trioxo-lanost-8-en-26-oic acid  
MF: C<sub>30</sub>H<sub>44</sub>O<sub>8</sub>  
FW: 532.7  
Purity:  $\geq$ 95% (mixture of isomers)  
UV/Vis.:  $\lambda_{\text{max}}$ : 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 G is supplied as a solid. A stock solution may be made by dissolving the ganoderic acid G in the solvent of choice, which should be purged with an inert gas. Ganoderic acid G is soluble in DMSO.

### Description

Ganoderic acid G is a triterpenoid that has been found in *G. lucidum* and has anticancer and anti-inflammatory activities.<sup>1,2</sup> It inhibits the proliferation of Meth A sarcoma cells (IC<sub>50</sub> = 6.8  $\mu$ g/ml).<sup>1</sup> Ganoderic acid G inhibits the production of nitric oxide (NO) in LPS-induced BV-2 microglial cells (IC<sub>50</sub> = 5.77  $\mu$ M).<sup>2</sup>

### References

1. Min, B.S., Gao, J.J., Nakamura, N., *et al.* Triterpenes from the spores of *Ganoderma lucidum* and their cytotoxicity against meth-A and LLC tumor cells. *Chem. Pharm. Bull. (Tokyo)* **48(7)**, 1026-1033 (2000).
2. Yang, J., Ting, X., Zou, L.-H., *et al.* Lanostane triterpenoids from *Ganoderma curtisii* and their NO production inhibitory activities of LPS-induced microglia. *Bioorg. Med. Chem. Lett.* **26(15)**, 3556-3561 (2018).

#### 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, 12/06/2022

#### 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