

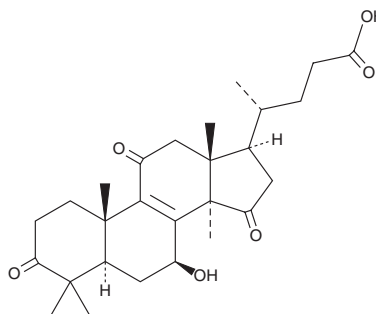
# PRODUCT INFORMATION



## Lucidenic Acid A

Item No. 30372

**CAS Registry No.:** 95311-94-7  
**Formal Name:** (5 $\alpha$ ,7 $\beta$ )-7-hydroxy-4,4,14-trimethyl-3,11,15-trioxo-chol-8-en-24-oic acid  
**Synonyms:** Lucidenic Acid A, Lucidenic Acid  $\alpha$   
**MF:** C<sub>27</sub>H<sub>38</sub>O<sub>6</sub>  
**FW:** 458.6  
**Purity:**  $\geq$ 95%  
**Supplied as:** A solid  
**Storage:** -20°C  
**Stability:**  $\geq$ 4 years  
**Item Origin:** Fungus/*Ganoderma* sp.



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

### Laboratory Procedures

Lucidenic acid A is supplied as a solid. A stock solution may be made by dissolving the lucidenic acid A in the solvent of choice, which should be purged with an inert gas. Lucidenic acid A is soluble in acetonitrile.

### Description

Lucidenic acid A is a triterpene originally isolated from *G. lucidum* and has anticancer activity.<sup>1,2</sup> It is cytotoxic to HepG2, HepG2 2.2.15', and P388 cancer cells (IC<sub>50</sub>s = 0.164, 0.205, and 17 nM, respectively). Lucidenic acid A inhibits increases in matrix metalloproteinase-2 (MMP-2) and MMP-9 activity and Matrigel™ invasion induced by phorbol 12-myristate 13-acetate (PMA; Item No. 10008014) in HepG2 cells.<sup>2</sup>

### References

1. Wu, T.-S., Shi, L.-S., and Kuo, S.-C. Cytotoxicity of *Ganoderma lucidum* triterpenes. *J. Nat. Prod.* **64**(8), 1121-1122 (2001).
2. Weng, C.-J., Chau, C.-F., Chen, K.-D., *et al.* The anti-invasive effect of lucidenic acids isolated from a new *Ganoderma lucidum* strain. *Mol. Nutr. Food Res.* **51**(12), 1472-1477 (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, 08/30/2023

#### 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