

PRODUCT INFORMATION



Hericene A

Item No. 45578

CAS Registry No.: 157207-54-0

Formal Name: hexadecanoic acid, [4-[(2E)-3,7-dimethyl-2,6-octadien-1-yl]-2-formyl-3-hydroxy-5-methoxyphenyl]methyl ester

Synonym: 5'-Deoxohericenone C

MF: C₃₅H₅₆O₅

FW: 556.8

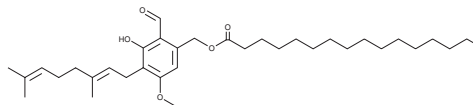
Purity: ≥98%

Supplied as: A solid

Storage: -20°C

Stability: ≥4 years

Item Origin: Fungus/*Heridium erinaceus*



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

Laboratory Procedures

Hericene A is supplied as a solid. A stock solution may be made by dissolving the hericene A in the solvent of choice, which should be purged with an inert gas. Hericene A is slightly soluble (0.1-1 mg/ml) in chloroform.

Description

Hericene A is a polyketide synthase-derived geranyl resorcyate that has been found in *H. erinaceus* and has diverse biological activities.¹⁻⁵ It is an inhibitor of α -glucosidase ($IC_{50} = 6.7 \mu M$).³ Hericene A (1-100 μM) decreases LPS-induced TNF- α , IL-6, and nitric oxide (NO) production in RAW 264.7 cells.⁴ It increases the axon length of primary rat hippocampal neurons and improves recognition memory in mice.⁵

References

1. Chen, J., Zeng, X., Yang, Y.L., *et al.* Genomic and transcriptomic analyses reveal differential regulation of diverse terpenoid and polyketides secondary metabolites in *Heridium erinaceus*. *Sci. Rep.* **7**(1), 10151 (2017).
2. Kobayashi, S., Hamada, Y., Yasumoto, T., *et al.* Total syntheses and endoplasmic reticulum stress suppressive activities of hericenones A-C and their derivatives. *Tetrahedron Lett.* **59**(18), 1733-1736 (2018).
3. Lee, S.K., Ryu, S.H., Turk, A., *et al.* Characterization of α -glucosidase inhibitory constituents of the fruiting body of lion's mane mushroom (*Heridium erinaceus*). *J. Ethnopharmacol.* **262**, 113197 (2020).
4. Xie, G., Tang, L., Xie, Y., *et al.* Secondary metabolites from *Heridium erinaceus* and their anti-inflammatory activities. *Molecules* **27**(7), 2157 (2022).
5. Martínez-Mármol, R., Chai, Y., Conroy, J.N., *et al.* Hericerin derivatives activates a pan-neurotrophic pathway in central hippocampal neurons converging to ERK1/2 signaling enhancing spatial memory. *J. Neurochem.* **165**(6), 791-808 (2023).

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, 06/03/2026

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