

PRODUCT INFORMATION



Hericerin

Item No. 39057

CAS Registry No.: 140381-53-9

Formal Name: 5-[(2E)-3,7-dimethyl-2,6-octadien-1-yl]-2,3-dihydro-4-hydroxy-6-methoxy-2-(2-phenylethyl)-1H-isoindol-1-one

Synonym: Isohericerin

MF: $C_{27}H_{33}NO_3$

FW: 419.6

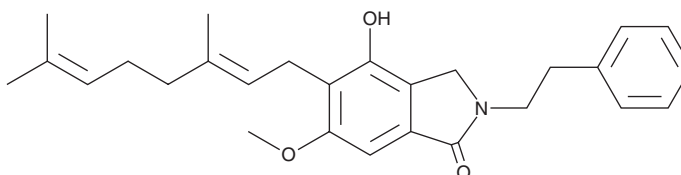
Purity: $\geq 98\%$

Supplied as: A solid

Storage: -20°C

Stability: ≥ 4 years

Item Origin: Synthetic



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

Laboratory Procedures

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

Description

Hericerin is a polyketide synthase-derived isoindolinone that has been found in *H. erinaceus* and has diverse biological activities.¹⁻³ It inhibits α -glucosidase ($IC_{50} = 5.3 \mu\text{M}$) and is cytotoxic to A549 lung and HeLa cervical cancer cells (IC_{50} s = 49 and 40.5 μM , respectively).² Hericerin induces neurite outgrowth in Neuro2a neuroblastoma cells and increases nerve growth factor (NGF) production in C6 glioma cells when used at concentrations of 25 and 50 μM , respectively.³

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 *Hericium erinaceus*. *Sci. Rep.* **7**(1), 10151 (2017).
2. Wang, K., Bao, L., Qi, Q., *et al.* Erinacerins C-L, isoindolin-1-ones with α -glucosidase inhibitory activity from cultures of the medicinal mushroom *Hericium erinaceus*. *J. Nat. Prod.* **78**(1), 146-154 (2015).
3. Ryu, S.H., Hong, S.M., Khan, Z., *et al.* Neurotrophic isoindolinones from the fruiting bodies of *Hericium erinaceus*. *Bioorg. Med. Chem. Lett.* **31**, 127714 (2021).

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, 05/07/2025

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