

PRODUCT INFORMATION



Hericenone D

Item No. 39785

CAS Registry No.: 137592-04-2
Formal Name: octadecanoic acid, [4-[(2E)-3,7-dimethyl-5-oxo-2,6-octadien-1-yl]-2-formyl-3-hydroxy-5-methoxyphenyl]methyl ester

MF: C₃₇H₅₈O₆

FW: 598.9

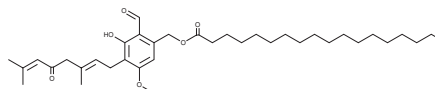
Purity: ≥98%

Supplied as: A solid

Storage: -20°C

Stability: ≥4 years

Item Origin: Fungus/*Hericium erinaceus*



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

Laboratory Procedures

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

Description

Hericenone D is a monoterpenoid originally isolated from *H. erinaceum*.¹

Reference

1. Kawagishi, H., Ando, M., Sakamoto, H., *et al.* Hericenones C, D and E, stimulators of nerve growth factor (NGF)-synthesis, from the mushroom *Hericium erinaceum*. *Tetrahedron Lett.* **35(32)**, 4561-4564 (1991).

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, 05/07/2026

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