

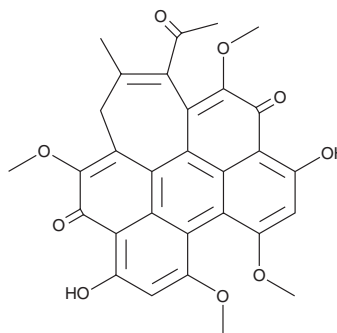
PRODUCT INFORMATION



Hypocrellin B

Item No. 28544

CAS Registry No.: 123940-54-5
Formal Name: 3-acetyl-6,11-dihydroxy-4,8,9,13-tetramethoxy-2-methyl-1H-cyclohepta[ghi]perylene-5,12-dione
MF: $C_{30}H_{24}O_9$
FW: 528.5
Purity: $\geq 98\%$
UV/Vis.: λ_{\max} : 336, 461, 549 nm
Supplied as: A crystalline solid
Storage: -20°C
Stability: ≥ 4 years
Item Origin: Fungus/*Hypocrella bambusae*



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

Laboratory Procedures

Hypocrellin B is supplied as a crystalline solid. A stock solution may be made by dissolving the hypocrellin B in the solvent of choice, which should be purged with an inert gas. Hypocrellin B is soluble in DMSO and dimethyl formamide.

Description

Hypocrellin B is a fungal metabolite that has been found in *H. bambusae* and has anticancer and antibacterial activities.^{1,2} It acts as a sonosensitizer, decreasing the viability of MDA-MB-231 breast cancer cells and inducing apoptosis in the presence of ultrasound exposure, effects that can be inhibited by the caspase inhibitor z-VAD-FMK.¹ It also increases the production of reactive oxygen species (ROS) in MDA-MB-231 cells when used at a concentration of 2.5 μM with concomitant ultrasound exposure. Hypocrellin B (20 μM), in combination with ultrasound sonication, is active against methicillin-resistant *S. aureus* (MRSA).²

References

1. Jia, Y., Wang, X., Liu, Q., *et al.* Sonodynamic action of hypocrellin B triggers cell apoptosis of breast cancer cells involving caspase pathway. *Ultrasonics* **73**, 154-161 (2017).
2. Leung, A.W.N., Ip, M., Xu, C.S., *et al.* Sonodynamic bactericidal efficacy of hypocrellin A and B against methicillin-resistant *Staphylococcus aureus*. *Hong Kong Med. J.* **23 Suppl 5(4)**, 36-37 (2017).

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, 08/16/2023

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