

# PRODUCT INFORMATION



## Hypocrellin A

Item No. 29933

**CAS Registry No.:** 77029-83-5  
**Formal Name:** 1-acetyl-2,3-dihydro-2,6,11-trihydroxy-4,8,9,13-tetramethoxy-2-methyl-1H-cyclohepta[ghi]perylene-5,12-dione, stereoisomer

**MF:** C<sub>30</sub>H<sub>26</sub>O<sub>10</sub>

**FW:** 546.5

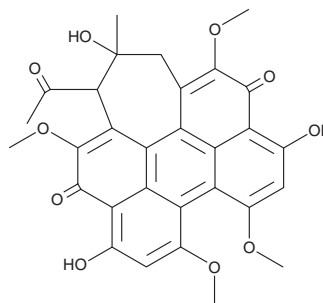
**Purity:** ≥98%

**Supplied as:** A solid

**Storage:** -20°C

**Stability:** ≥2 years

**Item Origin:** Fungi/*Hypocrella bambusae*



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

### Laboratory Procedures

Hypocrellin A is supplied as a solid. A stock solution may be made by dissolving the hypocrellin A in the solvent of choice, which should be purged with an inert gas. Hypocrellin A is soluble in ethanol, DMSO, and dimethyl formamide. The solubility of hypocrellin A in these solvents is 0.3, 3, and 2 mg/ml, respectively.

### Description

Hypocrellin A is a perylenequinoid that has been found in *Shiraia bambusicola* and has diverse biological activities.<sup>1,2</sup> It is active against *C. albicans*, *C. parapsilosis*, and *C. tropicalis* fungi (IC<sub>50</sub>s = 0.65, 1, and 85 µg/ml, respectively), as well as *S. aureus*, methicillin-resistant *S. aureus* (MRSA), *P. aeruginosa*, and *M. intracellulare* bacteria (IC<sub>50</sub>s = 3, 7, 10, and 2.5 µg/ml, respectively).<sup>1</sup> Hypocrellin A is also active against *L. donovani* promastigotes (IC<sub>50</sub> = 0.27 µg/ml). Upon photoactivation with UV radiation, hypocrellin A induces production of reactive oxygen species and apoptosis and reduces cell viability of A549 cancer cells.<sup>2</sup>

### References

1. Ma, G., Khan, S.I., Jacob, M.R., *et al.* Antimicrobial and antileishmanial activities of hypocrellins A and B. *Antimicrob. Agents Chemother.* **48(11)**, 4450-4452 (2004).
2. Qi, S., Guo, L., Yan, S., *et al.* Hypocrellin A-based photodynamic action induces apoptosis in A549 cells through ROS-mediated mitochondrial signaling pathway. *Acta Pharm. Sin. B.* **9(2)**, 279-293 (2019).

#### 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, 04/07/2020

#### 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