

PRODUCT INFORMATION

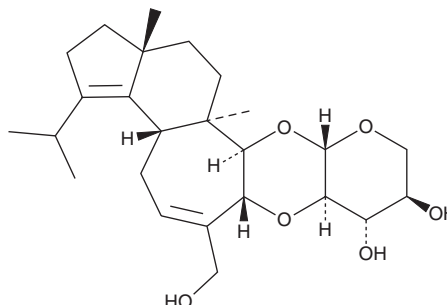


Erinacine C

Item No. 40079

CAS Registry No.: 156101-09-6
Formal Name: 1,2-O-[(3aR,5aR,6R,7R,10aR)-2,3,3a,4,5,5a,6,7,10,10a-decahydro-8-(hydroxymethyl)-3a,5a-dimethyl-1-(1-methylethyl)cyclohept[e]indene-6,7-diyl]-β-D-

Synonym: Erinacin C
MF: $C_{25}H_{38}O_6$
FW: 434.6
Purity: ≥95%
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

Erinacine C is supplied as a solid. A stock solution may be made by dissolving the erinacine C in the solvent of choice, which should be purged with an inert gas. Erinacine C is soluble (≥10 mg/ml) in ethanol and DMSO.

Description

Erinacine C is a cyathane diterpenoid that has been found in *H. erinaceum* and has diverse biological activities.¹⁻⁴ It inhibits transcription of the genes encoding the bacteriocins mutacin IV, mutacin V, and mutacin VI, which are *S. mutans* virulence factors, in reporter assays using *S. mutans*.² Erinacine C (1 mM) increases the production of nerve growth factor (NGF) in mouse astroglial cells.¹ It inhibits LPS-induced nitric oxide (NO) production in BV-2 microglia (IC₅₀ = 29.22 μM).³ Erinacine C (2 mg/kg) reduces mild traumatic brain injury-induced motor deficits and neuronal cell death in rats.⁴

References

1. Kawagishi, H., Shimada, A., Shirai, R., et al. Erinacines A, B and C, strong stimulators of nerve growth factor (NGF)-synthesis from the mycelia of *Hericium erinaceum*. *Tetrahedron Lett.* **35**(10), 1569-1572 (1994).
2. Premnath, P., Reck, M., Wittstein, K., et al. Screening for inhibitors of mutacin synthesis in *Streptococcus mutans* using fluorescent reporter strains. *BMC Microbiol.* **18**(1), 24 (2018).
3. Wei, J., Li, J.-Y., Feng, X.-L., et al. Unprecedented neoverrucosane and cyathane diterpenoids with anti-neuroinflammatory activity from cultures of the culinary-medicinal mushroom *Hericium erinaceus*. *Molecules* **28**(17), 6380 (2023).
4. K.F., L., Hsieh, Y.Y., Tung, S.Y., et al. The cerebral protective effect of novel erinacines from *Hericium erinaceus* Mycelium on in vivo mild traumatic brain injury animal model and primary mixed glial cells via Nrf2-dependent pathways. *Antioxidants (Basel)* **13**(3), 371 (2024).

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, 02/24/2025

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