

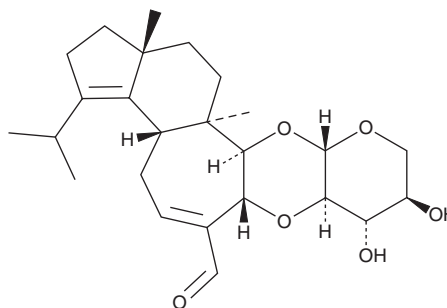
PRODUCT INFORMATION



Erinacine B

Item No. 40078

CAS Registry No.: 156101-10-9
Formal Name: 1,2-O-[(3aR,5aR,6R,7R,10aR)-8-formyl-2,3,3a,4,5,5a,6,7,10,10a-decahydro-3a,5a-dimethyl-1-(1-methylethyl)cyclohept[e]indene-6,7-diyl]-β-D-xylopyranose
Synonyms: Erinacin B, (-)-Erinacine B
MF: C₂₅H₃₆O₆
FW: 432.6
Purity: ≥98%
Supplied as: A solution in acetonitrile
Storage: -20°C
Stability: ≥1 year



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

Laboratory Procedures

Erinacine B is supplied as a solution in acetonitrile. To change the solvent, simply evaporate the acetonitrile under a gentle stream of nitrogen and immediately add the solvent of choice. Erinacine B is sparingly soluble (1-10 mg/ml) in methanol.

Description

Erinacine B is a cyathin diterpenoid fungal metabolite that has been found in *H. erinaceus*, also known as lion's mane mushroom.¹ It induces secretion of nerve growth factor (NGF) in isolated mouse astroglial cells when used at a concentration of 1 mM.

Reference

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).

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, 11/13/2024

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