

PRODUCT INFORMATION



Shearinine B

Item No. 45476

CAS Registry No.: 163136-25-2
Formal Name: (2R,4bS,6aS,9aS,15bS,15cR,17aS)-5,6,6a,7,9,9a,10,12,15,15b,15c,16,17,17a-tetradecahydro-4b-hydroxy-2-(1-hydroxy-1-methylethyl)-10,10,12,12,15b,15c-hexamethyl-2H-1-benzopyrano[5',6':6,7]indeno[1,2-b]pyrano[4',3':3,4]cyclopent [1,2-f]indol-3(4bH)-one

MF: C₃₇H₄₇NO₅

FW: 585.8

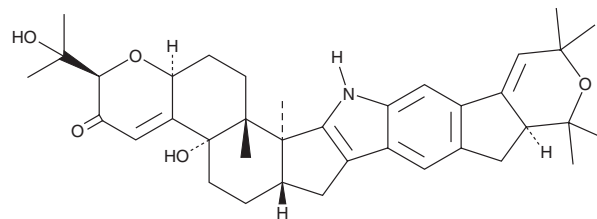
Purity: ≥95%

Supplied as: A solid

Storage: -20°C

Stability: ≥1 year

Item Origin: Bacterium/*Penicillium* sp.



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

Laboratory Procedures

Shearinine B is supplied as a solid. A stock solution may be made by dissolving the shearinine B in the solvent of choice, which should be purged with an inert gas. Shearinine B is soluble in methanol and DMSO.

Description

Shearinine B is an indole diterpene originally isolated from *E. shearii* that has antiinsectan activity.¹ Shearinine B (100 ppm in the diet) reduces the growth rate of *H. zea* larvae and the feeding rate of *C. hemipterus* larvae by 94 and 50%, respectively. It also induces 84% mortality in *S. frugiperda* larvae in a leaf disc assay.

Reference

1. Belofsky, G.N. and Gloer, J.B. Antiinsectan alkaloids: Shearinines A-C and a new paxilline derivative from the ascostromata of *Eupenicillium shearii*. *Tetrahedron* **51**(14), 3959-3968 (1995).

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, 05/22/2026

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