

PRODUCT INFORMATION



Shearamide A

Item No. 45506

CAS Registry No.: 212394-70-2
Formal Name: (6S,12S,15S,17aS,23S,26S,28aS)-12-((1H-indol-3-yl)methyl)-6-benzyl-15-((S)-sec-butyl)-23-((R)-1-hydroxyethyl)-26-isopropyloctadecahydro-1H,5H-dipyrrolo[1,2-a:1',2'-j][1,4,7,10,13,16,19,22]octaazacyclotetracosin-5,8,11,14,17,22,25,28(28aH)-octaone
Synonym: Cyclic(Pro-Val-Thr-Pro-Ile-Trp-Gly-Phe)

Peptide Sequence: c[PVTPIWGF]

MF: C₄₇H₆₃N₉O₉

FW: 898.1

Purity: ≥95%

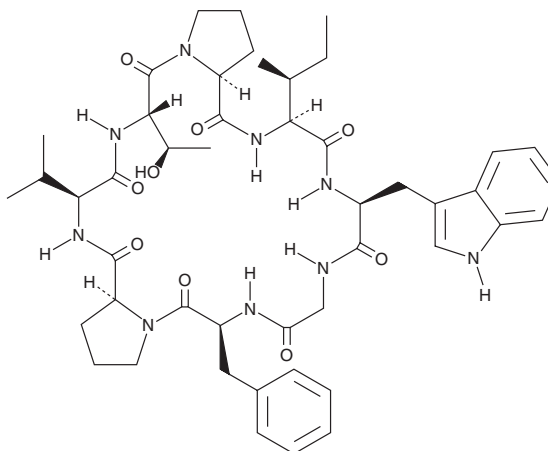
Supplied as: A solid

Storage: -20°C

Stability: ≥4 years

Item Origin: Fungus/*Talaromyces* sp.

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



Laboratory Procedures

Shearamide A is supplied as a solid. A stock solution may be made by dissolving the shearamide A in the solvent of choice, which should be purged with an inert gas. Shearamide A is slightly soluble (0.1-1 mg/ml) in methanol.

Description

Shearamide A is a cyclic octapeptide that has been found in *E. shearii* and has antiinsectan activity.¹ It reduces the growth rate of *H. zea* larvae by 35% when administered in the diet at 100 ppm.

Reference

1. Belofsky, G.N., Gloer, J.B., Wicklow, D.T., *et al.* Shearamide A: A new cyclic peptide from the ascostromata of *Eupenicillium shearii*. *Tetrahedron Lett.* **39(31)**, 5497-5500 (1998).

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, 04/13/2026

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