

PRODUCT INFORMATION



Beauvericin A

Item No. 31773

CAS Registry No.: 165467-50-5
Formal Name: (3S,6R,9S,12R,15S,18R)-4,10,16-trimethyl-6,12-bis(1-methylethyl)-18-[(1S)-1-methylpropyl]-3,9,15-tris(phenylmethyl)-1,7,13-trioxo-4,10,16-triazacyclooctadecane-2,5,8,11,14,17-hexone

MF: C₄₆H₅₉N₃O₉

FW: 798.0

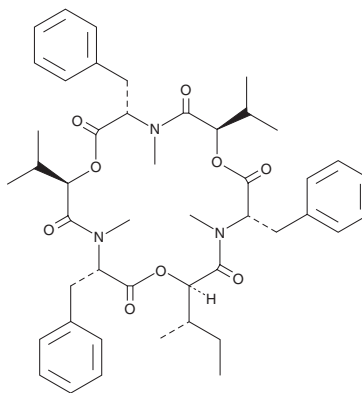
Purity: ≥95%

Supplied as: A solid

Storage: -20°C

Stability: ≥4 years

Item Origin: Fungus/*Beauveria bassiana*



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

Laboratory Procedures

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

Description

Beauvericin A is a cyclodepsipeptide and derivative of beauvericin (Item No. 11426) originally isolated from *B. bassiana* that has diverse biological activities.¹⁻³ It is active against *M. tuberculosis* (MIC = 25 µg/ml) and *P. falciparum* (IC₅₀ = 12 µg/ml).² Beauvericin A is toxic to brine shrimp (LD₁₀₀ = 32 µg/ml).³

References

1. Gupta, S., Montillor, C., and Hwang, Y.-S. Isolation of novel beauvericin analogues from the fungus *Beauveria bassiana*. *J. Nat. Prod.* **58(5)**, 733-738 (1995).
2. Nilanonta, C., Isaka, M., Kittakoop, P., et al. Antimycobacterial and antiplasmodial cyclodepsipeptides from the insect pathogenic fungus *Paecilomyces tenuipes* BCC 1614. *Planta Med.* **66(8)**, 756-758 (2000).
3. Shi, S., Li, Y., Ming, Y., et al. Biological activity and chemical composition of the endophytic fungus *Fusarium* sp. TP-G1 obtained from the root of *Dendrobium officinale* Kimura et Migo. *Rec. Nat. Prod.* **12(6)**, 549-556 (2018).

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, 08/11/2023

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