

PRODUCT INFORMATION



Enniatin A

Item No. 17456

CAS Registry No.: 2503-13-1
Formal Name: cyclo[(2R)-2-hydroxy-3-methylbutanoyl-N-methyl-L-isoleucyl-(2R)-2-hydroxy-3-methylbutanoyl-N-methyl-L-isoleucyl-(2R)-2-hydroxy-3-methylbutanoyl-N-methyl-L-isoleucyl]

MF: C₃₆H₆₃N₃O₉

FW: 681.9

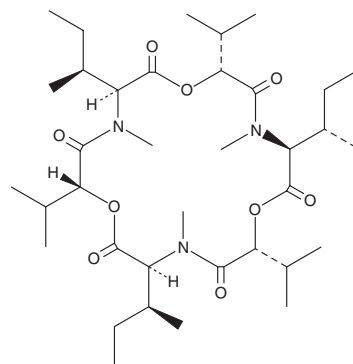
Purity: ≥99%

Supplied as: A solid

Storage: -20°C

Stability: ≥4 years

Item Origin: Fungus/*Fusarium* sp.



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

Laboratory Procedures

Enniatin A is supplied as a solid. A stock solution may be made by dissolving the enniatin A in the solvent of choice. Enniatin A is soluble in organic solvents such as ethanol, methanol, DMSO, and dimethyl formamide, which should be purged with an inert gas.

Description

Enniatins are cyclohexadepsipeptides commonly isolated from fungi that are known to have antibiotic properties and to induce apoptosis in several cancer lines.¹ Many function as ionophores, forming pores in cellular membranes to allow selective ion transport.^{1,2} Enniatin A is one of four major analogs in the enniatin complex (Item No. 9002040). It has been shown to moderately inhibit acyl-CoA:cholesterol acyltransferase activity in rat liver microsomes with an IC₅₀ value of 22 μM.³ Enniatin A also demonstrates anthelmintic properties against *N. brasiliensis*, *T. spiralis*, and *H. spumosa*.¹

References

1. Sy-Cordero, A.A., Pearce, C.J., and Oberlies, N.H. Revisiting the enniatins: A review of their isolation, biosynthesis, structure determination and biological activities. *J. Antibiot. (Tokyo)* **65(11)**, 541-549 (2012).
2. Kamyar, M.R., Rawnduzi, P., Studenik, C.R., et al. Investigation of the electrophysiological properties of enniatins. *Arch. Biochem. Biophys.* **429(2)**, 215-223 (2004).
3. Tomoda, H., Huang, X.H., Cao, J., et al. Inhibition of acyl-CoA: Cholesterol acyltransferase activity by cyclodepsipeptide antibiotics. *J. Antibiot. (Tokyo)* **45(10)**, 1626-1632 (1992).

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, 08/08/2023

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