

PRODUCT INFORMATION



Enopeptin A

Item No. 17092

CAS Registry No.: 139601-96-0
Formal Name: (4R)-N-[(2E,4E,6E,8E,10E)-12-[(2-hydroxy-5-oxo-1-cyclopenten-1-yl)amino]-1,12-dioxo-2,4,6,8,10-dodecapentaen-1-yl]-L-phenylalanyl-L-seryl-L-prolyl-L-proline, (6→2)-lactone

Synonyms: BRN 537924, NSC 657143

MF: C₄₇H₅₇N₇O₁₁

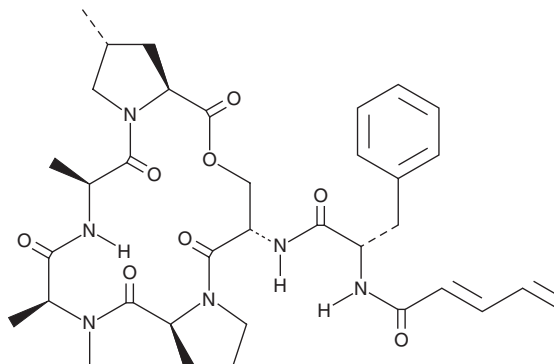
FW: 896.0

Purity: ≥95%

Supplied as: A solid

Storage: -20°C

Stability: ≥4 years



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

Laboratory Procedures

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

Description

Enopeptin A, originally isolated from a culture broth of *Streptomyces* sp. RK-1051, is a depsipeptide antibiotic that contains two unusual amino acids (N-methylalanine and 4-methylproline) and features a pentaenone side chain.^{1,2} It is effective against Gram-positive bacteria, including methicillin-resistant *S. aureus* (MIC = 25 µg/ml), and Gram-negative bacteria, including mutant forms of *E. coli* and *P. aeruginosa* (MICs = 200 µg/ml); however, it is not inhibitory to fungi.¹

References

1. Osada, H., Yano, T., Koshino, H., *et al.* Enopeptin A, a novel depsipeptide antibiotic with anti-bacteriophage activity. *J. Antibiot.* **44(12)**, 1463-1466 (1991).
2. Young, J.-J., Jung, L.-J., Liu, W.-T., *et al.* Solution conformation of enopeptin A, a depsipeptide antibiotic, using 2D NMR and restrained molecular dynamics studies. *J. Antibiot.* **47(8)**, 922-931 (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, 10/11/2022

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