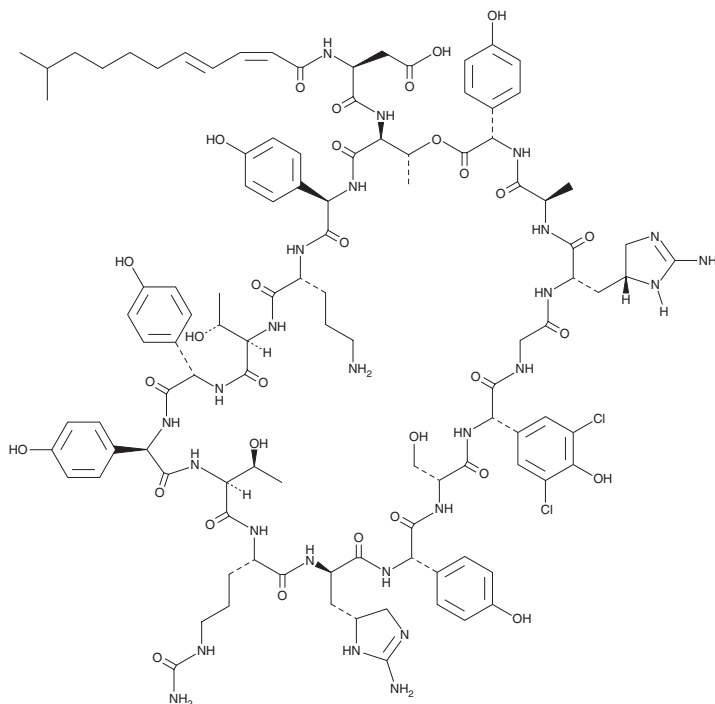


PRODUCT INFORMATION



Enramycin A Item No. 40847

CAS Registry No.: 34438-27-2
Formal Name: (2S)-N-[(2Z,4E)-10-methyl-1-oxo-2,4-undecadien-1-yl]-L- α -aspartyl-L-threonyl-(2R)-2-(4-hydroxyphenyl)glycyl-D-ornithyl-D-allothreonyl-(2S)-2-(4-hydroxyphenyl)glycyl-(2R)-2-(4-hydroxyphenyl)glycyl-L-allothreonyl-N⁵-(aminocarbonyl)-L-ornithyl-3-[(5R)-2-amino-4,5-dihydro-1H-imidazol-5-yl]-D-alanyl-(2S)-2-(4-hydroxyphenyl)glycyl-D-seryl-(2S)-2-(3,5-dichloro-4-hydroxyphenyl)glycylglycyl-3-[(5R)-2-amino-4,5-dihydro-1H-imidazol-5-yl]-L-alanyl-D-alanyl-2-(4-hydroxyphenyl)glycine (17 $\rightarrow$ 2)-lactone
Synonym: Enduramycin A
MF: C₁₀₇H₁₃₈Cl₂N₂₆O₃₁
FW: 2,355.3
Purity: $\geq$ 95%
Supplied as: A solid
Storage: -20°C
Stability: $\geq$ 4 years



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

Laboratory Procedures

Enramycin A is supplied as a solid. A stock solution may be made by dissolving the enramycin A in the solvent of choice, which should be purged with an inert gas. Enramycin A is slightly soluble (0.1-1 mg/ml) in methanol. It is also slightly soluble (0.1-1 mg/ml) in water. We do not recommend storing the aqueous solution for more than one day.

Description

Enramycin A is a lipopeptide antibiotic that has been found in *S. fungicidicus*.¹ It also inhibits avian myeloblastosis virus reverse transcriptase by 50% when used at a concentration of 10 μ g/ml in a cell-free assay.² Formulations containing enramycin A have been used as feed additives to prevent bacterial necrotic enteritis in livestock.

References

1. Iwasaki, H., Horii, S., Asai, M., *et al.* Enduracidin, a new antibiotic. VIII. Structures of enduracidins A and B. *Chem. Pharm. Bull.* **21**(6), 1184-1191 (1973).
2. Inouye, Y., Take, Y., Nakamura, S., *et al.* Screening for inhibitors of avian myeloblastosis virus reverse transcriptase and effect on the replication of AIDS-virus. *J. Antibiot. (Tokyo)* **40**(1), 100-104 (1987).

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, 04/29/2024

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