

PRODUCT INFORMATION



Dihydrotetrodecamycin

Item No. 29725

CAS Registry No.: 166403-10-7
Formal Name: [3S-(3R*,5R*,6S*,6aR*,10aR*,11S*,13R*)]-6,6a,7,8,9,10,10a,11-octahydro-6,6a-dihydroxy-3,11,13-trimethyl-5,11-methanofuro[3,4-d][3]benzoxonin-1,12(3H,5H)-dione

Synonym: Streptonolid A

MF: C₁₈H₂₄O₆

FW: 336.4

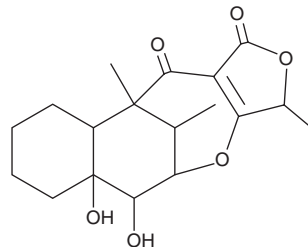
Purity: ≥70%

Supplied as: A solid

Storage: -20°C

Stability: ≥4 years

Item Origin: Bacteria/*Streptomyces zaomyceticus*



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

Laboratory Procedures

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

Description

Dihydrotetrodecamycin is a fungal metabolite originally isolated from *S. nashvillensis* that has antibacterial activity.^{1,2} It is active against the piscine pathogenic bacterium *P. piscicida*.

References

1. Tsuchida, T., Linuma, H., Nishida, C., *et al.* Tetrodecamycin and dihydrotetrodecamycin, new antimicrobial antibiotics against *Pasteurella piscicida* produced by *Streptomyces nashvillensis* MJ885-mF8. I. Taxonomy, fermentation, isolation, characterization and biological activities. *J. Antibiot. (Tokyo)* **48(10)**, 1104-1109 (1995).
2. Tsuchida, T., Linuma, H., Nakamura, K.T., *et al.* Derivatives of tetrodecamycin. *J. Antibiot. (Tokyo)* **48(11)**, 1330-1335 (1995).

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, 10/12/2022

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