

PRODUCT INFORMATION



17-hydroxy Venturicidin A

Item No. 27508

CAS Registry No.: 113204-43-6
Formal Name: (1R,5R,6R,8S,9E,11R,15E,17R)-11-[[3-O-(aminocarbonyl)-2,6-dideoxy-β-D-arabino-hexopyranosyl]oxy]-1,7-dihydroxy-5-[(1R,3R,4S,5S)-4-hydroxy-1,3,5-trimethyl-6-oxooctyl]-6,8,16,18-tetramethyl-4,21-dioxabicyclo[15.3.1]heneicosa-9,15,18-trien-3-one

Synonym: YP 02259L-C

MF: C₄₁H₆₇NO₁₂

FW: 766.0

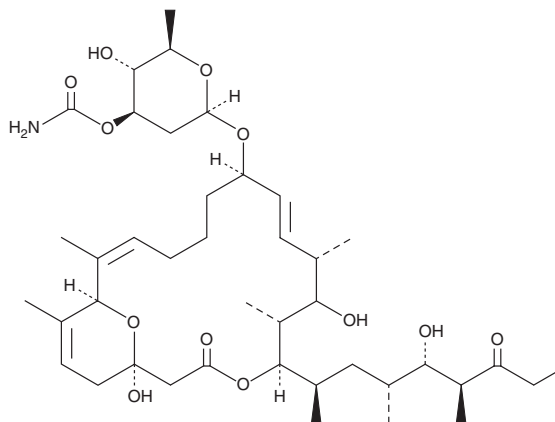
Purity: ≥95%

Supplied as: A solid

Storage: -20°C

Stability: ≥4 years

Item Origin: Bacterium/*Streptomyces* sp.



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

Laboratory Procedures

17-hydroxy Venturicidin A is supplied as a solid. A stock solution may be made by dissolving the 17-hydroxy venturicidin A in the solvent of choice, which should be purged with an inert gas. 17-hydroxy Venturicidin A is soluble in the organic solvent ethanol at a concentration of approximately 1 mg/ml. 17-hydroxy venturicidin A is also soluble in methanol and DMSO.

Description

17-hydroxy Venturicidin A is a macrolide fungal metabolite originally isolated from *Streptomyces*.¹ It has antibiotic activity against *M. luteus*, *B. subtilis*, and *S. aureus* and antifungal activity against *V. dahlia*, *Fusarium*, and *C. tropicalis* in a disc assay.

Reference

1. Fourati-Ben Fguira, L., Fotso, B., Ben Ameer-Mehdi, R., *et al.* Purification and structure elucidation of antifungal and antibacterial activities of newly isolated *Streptomyces* sp. strain US80. *Res. Microbiol.* **156**(3), 241-347 (2004).

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, 12/22/2022

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