

PRODUCT INFORMATION



IT-143A

Item No. 30044

CAS Registry No.: 183485-32-7
Formal Name: 2-(10-hydroxy-3,5,7,9,11,13-hexamethyl-2,5,7,11,13-pentadecapentaenyl)-5,6-dimethoxy-3-methyl-4-pyridinol

MF: C₂₉H₄₃NO₄

FW: 469.7

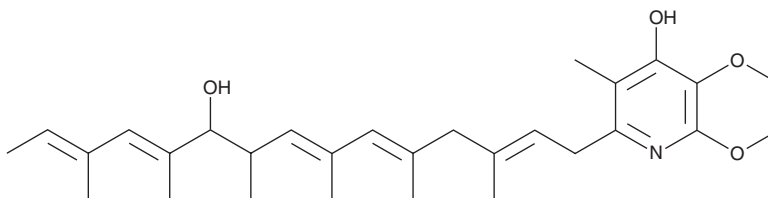
Purity: ≥95%

Supplied as: An oil

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

IT-143A is supplied as an oil. A stock solution may be made by dissolving the IT-143A in the solvent of choice, which should be purged with an inert gas. IT-143A is soluble in ethanol, methanol, DMSO, and dimethyl formamide.

Description

IT-143A is a bacterial metabolite originally isolated from *Streptomyces* sp. IT-143.¹ It is active against the bacterium *M. luteus* (MIC = 6.25 µg/ml) and the fungi *A. fumigatus* and *T. rubrum* *in vitro* (MICs = 12.5-25 µg/ml).

Reference

1. Urakawa, A., Sasaki, T., Yoshida, K., *et al.* IT-143-A and B, novel piericidin-group antibiotics produced by *Streptomyces* Sp. *J. Antibiot. (Tokyo)* **49(10)**, 1052-1055 (1996).

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, 10/25/2022

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