

PRODUCT INFORMATION



IT-143B

Item No. 29960

CAS Registry No.: 183485-34-9
Formal Name: 2-[(2E,5E,7E,9R,10R,11E,13E)-10-hydroxy-3,7,9,11,13-pentamethyl-2,5,7,11,13-pentadecapentaen-1-yl]-5,6-dimethoxy-3-methyl-4-pyridinol

MF: $C_{28}H_{41}NO_4$

FW: 455.6

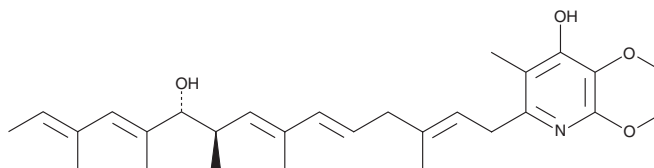
Purity: $\geq 95\%$

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

Description

IT-143B is a bacterial metabolite that has been found in *S. iakyrus*.¹ It inhibits the proliferation of OS-RC-2 and ACHN cancer cells ($\text{IC}_{50}\text{s} = 22$ and $98 \mu\text{M}$, respectively).

Reference

1. Li, K., Liang, Z., Chen, W., *et al.* Iakyrindins A-D, antiproliferative piericidin analogues bearing a carbonyl group or cyclic skeleton from *Streptomyces iakyrus* SCSIO NS104. *J. Org. Chem.* **84**(19), 12626-12631 (2019).

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