

PRODUCT INFORMATION



BE-26263

Item No. 29070

CAS Registry No.: 147716-81-2
Formal Name: 7,8,11,12,15,16,23,24,27,28-decahydro-2,4,18,20-tetrahydroxy-7,11,15,23,27-pentamethyl-5H,9H,13H,21H,25H-dibenzo[k,u][1,5,9,15,19]pentaoxacyclotetracosin-5,9,13,21,25-pentone

MF: C₃₂H₃₈O₁₄

FW: 646.6

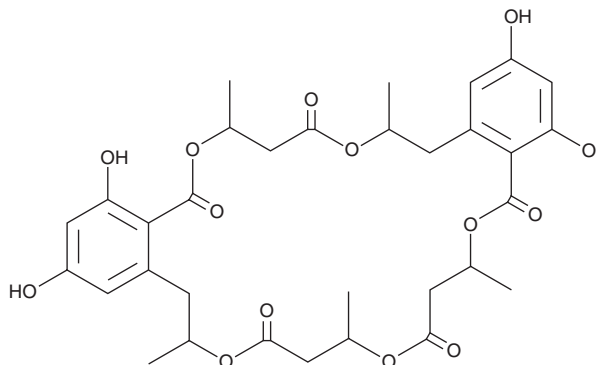
Purity: ≥70%

Supplied as: A solid

Storage: -20°C

Stability: ≥4 years

Item Origin: Fungus/*Pseudallescheria* sp.



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

Laboratory Procedures

BE-26263 is supplied as a solid. A stock solution may be made by dissolving the BE-26263 in the solvent of choice, which should be purged with an inert gas. BE-26263 is soluble in ethanol, methanol, dichloromethane, and DMSO.

Description

BE-26263 is a fungal metabolite that has been found in *S. apiospermum*.¹ It inhibits [¹²⁵I]-estradiol receptor binding with an IC₅₀ value of 0.96 μM.

Reference

1. Kondo, H., Kurama, M., Nakajima, S., *et al.* Estrogenic BE-26263 and its manufacture with *Scedosporium apiospermum*. *Banyu Pharma Co., Ltd. JP05032658* (1993).

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, 08/16/2023

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