

PRODUCT INFORMATION



FR901379

Item No. 29648

CAS Registry No.: 144371-88-0
Formal Name: 1-[(4R,5R)-4,5-dihydroxy-N²-(1-oxohexadecyl)-L-ornithine]-4-[(4S)-4-hydroxy-4-[4-hydroxy-3-(sulfoxy)phenyl]-L-threonine]-pneumocandin A₀

MF: C₅₁H₈₂N₈O₂₁S

FW: 1,175.3

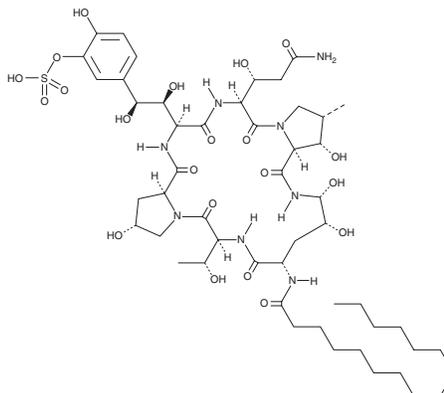
Purity: ≥90%

Supplied as: A solid

Storage: -20°C

Stability: ≥4 years

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



Laboratory Procedures

FR901379 is supplied as a solid. A stock solution may be made by dissolving the FR901379 in the solvent of choice, which should be purged with an inert gas. FR901379 is slightly soluble in methanol.

Description

FR901379 is an echinocandin-like antifungal lipopeptide.¹ It is active against *C. albicans*, *C. krusei*, *C. tropicalis*, *C. utilis*, *A. fumigatus*, and *A. niger* fungi (IC₅₀s = <0.003-1.9 μg/ml). FR901379 is protective against *C. albicans* infection in mice (ED₅₀ = 1.1 mg/kg). It also reduces the number of pulmonary cysts and trophozoites in a mouse model of *P. carinii* infection.

Reference

1. Fujie, A., Iwamoto, T., Sato, B., *et al.* FR131535, a novel water-soluble echinocandin-like lipopeptide: Synthesis and biological properties. *Bioorg. Med. Chem. Lett.* **11(3)**, 399-402 (2001).

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, 12/09/2022

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