

PRODUCT INFORMATION



A-83016F

Item No. 23240

CAS Registry No.: 142435-72-1
Formal Name: [2R-[2a,2[S*[1S*[2R*(2E,4E,6E),3R*,5R*),2S*,3E,5E]],3β,4β,6β (1E,3Z)]-7-[tetrahydro-3-hydroxy-5-[2-methoxy-1,3-dimethyl-7-[[1-oxo-2-[tetrahydro-2,3,4-trihydroxy-5,5-dimethyl-6-(1,3-pentadienyl)-2H-pyran-2-yl]butyl]amino]-3,5-heptadienyl]-2-furanyl]-2,4,6-heptatrienoic acid

Synonym: Antibiotic A-83016F

MF: C₃₇H₅₅NO₁₀

FW: 673.8

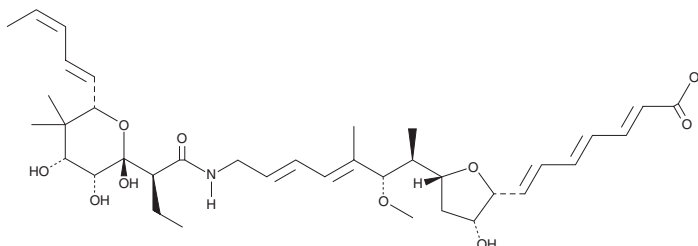
Purity: ≥70%

Supplied as: A solid

Storage: -20°C

Stability: ≥4 years

Item Origin: *Streptomyces* sp



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

Laboratory Procedures

A-83016F is supplied as a solid. A stock solution may be made by dissolving the A-83016F in the solvent of choice. A-83016F is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide.

Description

A-83016F is an aurodox antibiotic isolated from the actinomycete species A83016.¹ It has weak activity against *S. pneumoniae* PARK, *S. pyogenes*, and *H. influenzae* with MIC values of 16, 32, and 64 µg/ml, respectively, in a broth dilution assay.

Reference

1. Smitka, T.A., Bonjouklian, R., Perun, T.J., Jr., et al. A83016F, a new member of the aurodox family. *J. Antibiot. (Tokyo)* 45(4), 433-443 (1992).

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, 11/30/2022

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