

PRODUCT INFORMATION



A 26771B

Item No. 29593

CAS Registry No.: 56448-20-5
Formal Name: butanedioic acid, 1-[(3E,6S,16R)-16-methyl-2,5-dioxooxacyclohexadec-3-en-6-yl] ester

MF: C₂₀H₃₀O₇

FW: 382.5

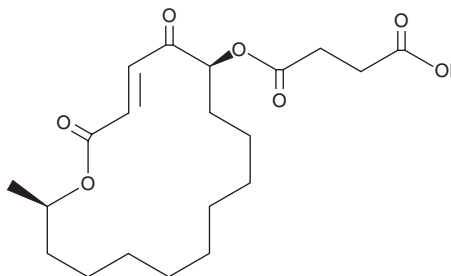
Purity: ≥95%

Supplied as: A solid

Storage: -20°C

Stability: ≥4 years

Item Origin: Fungus/*Penicillium* sp.



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

Laboratory Procedures

A 26771B is supplied as a solid. A stock solution may be made by dissolving the A 26771B in the solvent of choice, which should be purged with an inert gas. A 26771B is soluble in organic solvents such as ethanol, methanol, DMSO, and dimethyl formamide.

Description

A 26771B is a macrolide antibiotic originally isolated from *P. turbatum*.¹ It is active against *S. aureus*, *S. faecalis*, *V. coli*, *M. gallisepticum*, *M. granularum*, *M. synoviae*, *M. hyosynoviae*, *M. hyopneumoniae*, *E. amylovora*, *P. multocida*, and *X. phaseoli* bacteria (MICs = <0.78-50 µg/ml). A 26771B is active against various methicillin-resistant *S. aureus* (MRSA) strains (MICs = 6-16 µg/ml).² A 26771B is also active against *C. tropicalis*, *T. mentagrophytes*, *B. cinerea*, *C. ulmi*, and *V. albo-atrum fungi* (MICs = 6.25-100 µg/ml).¹

References

1. Michel, K.H., Demarco, P.V., and Nagarjan, R. The isolation and structure elucidation of macrocyclic lactone antibiotic, A26771B. *J. Antibiot. (Tokyo)* **30(7)**, 571-575 (1977).
2. Stierle, A.A., Stierele, D.B., Decato, D., et al. The berkeleylactones, antibiotic macrolides from fungal coculture. *J. Nat. Prod.* **80(4)**, 1150-1160 (2017).

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, 08/15/2023

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