

PRODUCT INFORMATION



Lydicamycin

Item No. 20584

CAS Registry No.: 133352-27-9

Formal Name: 2-[21-[1-[(2,5-dihydro-4-hydroxy-2-oxo-1H-pyrrol-3-yl)carbonyl]-1,2,4a,5,6,7,8,8a-octahydro-5,6-dihydroxy-1,3-dimethyl-2-naphthalenyl]-2,6,10,14,18-pentahydroxy-5,11,17,19-tetramethyl-4,8,12,16-heneicosatetraen-1-yl]1-pyrrolidinecarboximidamide

MF: $C_{47}H_{74}N_4O_{10}$

FW: 855.10

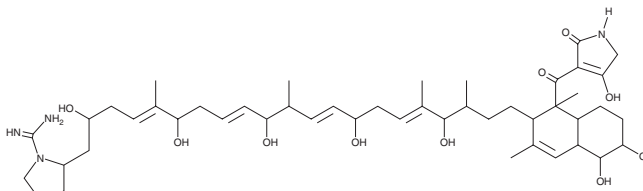
Purity: $\geq 95\%$

Supplied as: A film

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

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

Description

Lydicamycin is an antibiotic first isolated from *Streptomyces*. It is particularly cytotoxic against *S. aureus*, including the methicillin resistant 535 strain (MICs = 3.1-6.2 $\mu\text{g/ml}$).¹ Lydicamycin also inhibits the growth of *C. neoformans* (MIC = 25 $\mu\text{g/ml}$) but is ineffective against Gram-negative bacteria.¹

Reference

1. Hayakawa, Y., Kanamaru, N., Shimazu, A., *et al.* Lydicamycin, a new antibiotic of a novel skeletal type. I. Taxonomy, fermentation, isolation and biological activity. *J. Antibiot. (Tokyo)* **44**(3), 282-287 (1991).

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, 12/19/2022

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