

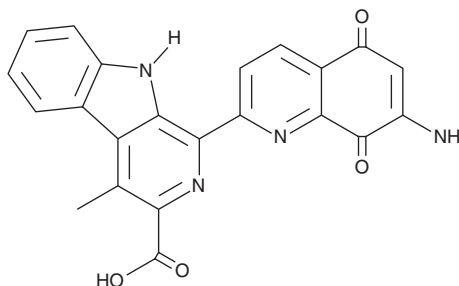
PRODUCT INFORMATION



Lavendamycin

Item No. 42360

CAS Registry No.: 81645-09-2
Formal Name: 1-(7-amino-5,8-dihydro-5,8-dioxo-2-quinoliny)-4-methyl-9H-pyrido[3,4-b]indole-3-carboxylic acid
Synonym: NSC 322370
MF: C₂₂H₁₄N₄O₄
FW: 398.4
Purity: ≥90%
Supplied as: A solid
Storage: -20°C
Stability: ≥4 years
Item Origin: Bacterium/*Actinoplanes* sp.



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

Description

Lavendamycin is a bacterial metabolite that has been found in *S. lavendulae* and has antimicrobial and anticancer activities.^{1,2} It is active against various bacteria, including *S. pneumoniae*, *S. aureus*, *E. coli*, *P. mirabilis*, and *P. vulgaris* (MICs = 0.13-1 µg/ml), and various fungi, including *T. rubrum*, *T. mentagrophytes*, and *M. canis* (MICs = 0.5 µg/ml for all).¹ Lavendamycin is cytotoxic to P388 murine leukemia, MKN45 gastric carcinoma, and WiDr colon adenocarcinoma cells (IC₅₀s = 0.06, 0.1, 0.09 µg/ml, respectively).² *In vivo*, lavendamycin (0.4-6.4 mg/kg) increases median survival time in a P388 murine leukemia model.¹

References

1. Balitz, D.M., Bush, J.A., Bradner, W.T., et al. Isolation of lavendamycin, a new antibiotic from *Streptomyces lavendulae*. *J. Antibiot. (Tokyo)* **35**(3), 259-265 (1982).
2. Abe, N., Nakakita, Y., Nakamura, T., et al. Novel cytotoxic compounds, oxopropalines from *Streptomyces* sp. G324 producing lavendamycin. I. Taxonomy of the producing organism, fermentation, isolation and biological activities. *J. Antibiot. (Tokyo)* **46**(11), 1672-1677 (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, 07/02/2025

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