

PRODUCT INFORMATION

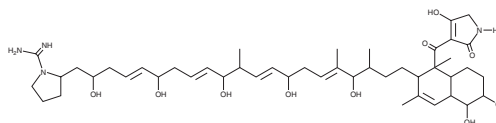


TPU-0037A

Item No. 25761

CAS Registry No.: 485815-59-6

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-11,17,19-trimethyl-4,8,12,16-heneicosatetraen-1-yl]-1-pyrrolidinecarboximidamide



MF: C₄₆H₇₂N₄O₁₀

FW: 841.1

Purity: ≥95%

Supplied as: A film

Storage: -20°C

Stability: ≥4 years

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

Laboratory Procedures

TPU-0037A is supplied as a film. A stock solution may be made by dissolving the TPU-0037A in the solvent of choice, which should be purged with an inert gas. TPU-0037A is soluble in organic solvents such as ethanol, methanol, DMSO, and dimethyl formamide.

TPU-0037A is sparingly soluble in aqueous solutions. To enhance aqueous solubility, dilute the organic solvent solution into aqueous buffers or isotonic saline. If performing biological experiments, ensure the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. We do not recommend storing the aqueous solution for more than one day.

Description

TPU-0037A is an antibiotic and a congener of lydicamycin (Item No. 20584).¹ It inhibits the growth of Gram-positive bacteria, including methicillin-resistant *S. aureus* (MRSA), *B. subtilis*, and *M. luteus* (MICs = 1.56-12.5 µg/ml), but not Gram-negative *E. coli*, *P. mirabilis*, *P. vulgaris*, or *P. aeruginosa* strains (MICs = >50 µg/ml).

Reference

1. Furumai, T., Eto, K., Sasaki, T., *et al.* TPU-0037-A, B, C and D, novel lydicamycin congeners with anti-MRSA activity from *Streptomyces platensis* TP-A0598. *J. Antibiot. (Tokyo)* **55**(10), 873-880 (2002).

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, 12/20/2022

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