

PRODUCT INFORMATION

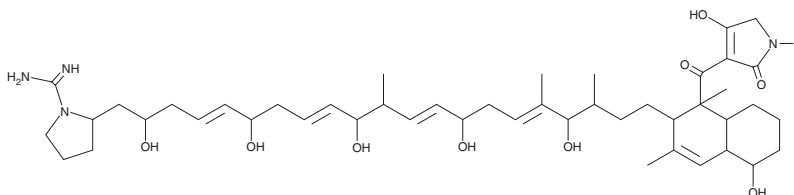


TPU-0037C

Item No. 24179

CAS Registry No.: 485815-61-0

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-hydroxy-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: $C_{46}H_{72}N_4O_9$

FW: 825.1

Purity: $\geq 95\%$

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

Description

TPU-0037C is a metabolite of the marine actinomycete *S. platensis* that is structurally similar to lydicamycin (Item No. 20584).¹ It is active against Gram-positive bacteria (MICs = 0.39-3.13 $\mu\text{g}/\text{ml}$), including methicillin-resistant *S. aureus* strains (MIC = 3.13 $\mu\text{g}/\text{ml}$), but is ineffective against Gram-negative bacteria (MICs = >50 $\mu\text{g}/\text{ml}$).

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, 11/16/2022

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