

PRODUCT INFORMATION

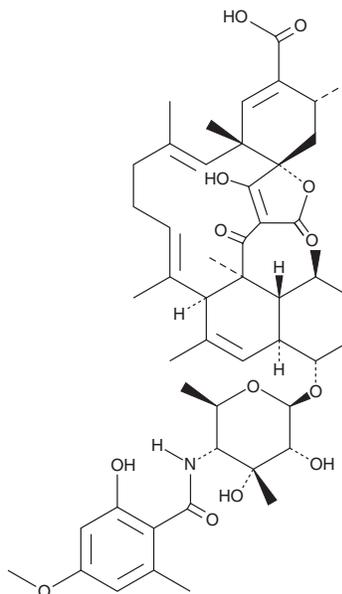


Tetromycin C5

Item No. 44801

CAS Registry No.: 3040055-61-3
Formal Name: (1S,4S,4aS,6aR,7E,11E,12aR,15R,16aS,19E,20aS,20bR)-4-(((2R,3R,4R,5R,6R)-3,4-dihydroxy-5-(2-hydroxy-4-methoxy-6-methylbenzamido)-4,6-dimethyltetrahydro-2H-pyran-2-yl)oxy)-21-hydroxy-1,6,7,11,12a,15,20a-heptamethyl-18,20-dioxo-1,2,3,4,4a,6a,9,10,12a,15,16,20,20a,20b-tetradecahydro-18H-16a,19-(metheno)benzo[b]naphtho[2,1-j][1]oxacyclotetradecine-14-carboxylic acid

Synonym: Glenthmycin I
MF: C₅₀H₆₅NO₁₃
FW: 888.1
Purity: ≥95%
Supplied as: A solid
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

Tetromycin C5 is supplied as a solid. A stock solution may be made by dissolving the tetromycin C5 in the solvent of choice, which should be purged with an inert gas. Tetromycin C5 is soluble in DMSO and methanol.

Description

Tetromycin C5 is a polyketide that has been found in *Streptomyces* and has antibiotic activity.¹ It is active against *S. aureus*, *E. faecalis*, methicillin-resistant *S. aureus* (MRSA), and vancomycin-resistant Enterococci (VRE; IC₅₀s = 10, 13, 10.2, and 7.3 μM, respectively).

Reference

- 1 Wu, T., Salim, A.A., Khalil, Z.G., et al. Glenthmycins A-M: Macrocyclic spiro-tetronate polyketide antibacterials from the Australian pasture plant-derived *Streptomyces* sp. CMB-PB041. *J. Nat. Prod.* **85**(6), 1641-1657 (2022).

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, 12/01/2025

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