

PRODUCT INFORMATION



Glenthmycin L

Item No. 45513

Formal Name: (1S,4S,4aS,6aR,7E,11E,12aR,15R,16aS,19E,20aS,20bR)-21-hydroxy-1,6,7,11,14,15,20a-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]oxacyclotetradecin-4-yl acetate

MF: C₃₅H₄₆O₆

FW: 562.7

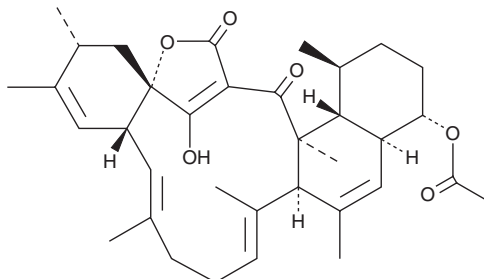
Purity: ≥95%

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

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

Description

Glenthmycin L is a macrocyclic spirotetronate originally isolated from *Streptomyces* that has antibacterial activity.¹ It is active against *S. aureus*, methicillin-resistant *S. aureus* (MRSA), and *E. faecalis* (IC₅₀S = 1.2, 0.9, and 3.7 μM, respectively).

Reference

1. Wu, T., Salim, A.A., Khalil, Z.G., *et al.* Glenthmycins A-M: Macrocyclic spirotetronate 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, 05/21/2026

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