

# PRODUCT INFORMATION



## Telomycin

Item No. 28137

CAS Registry No.: 19246-24-3

Formal Name: (3R)-L-β-aspartyl-L-seryl-L-threonyl-L-allothreonyl-L-alanylglycyl-(3S)-3-hydroxy-L-prolyl-α,β-didehydrotryptophyl-β-methyl-L-tryptophyl-(3S)-3-hydroxy-L-leucyl-3-hydroxy-L-proline, (11→3)-lactone

Synonym: NSC 235069

MF: C<sub>59</sub>H<sub>77</sub>N<sub>13</sub>O<sub>19</sub>

FW: 1,272.3

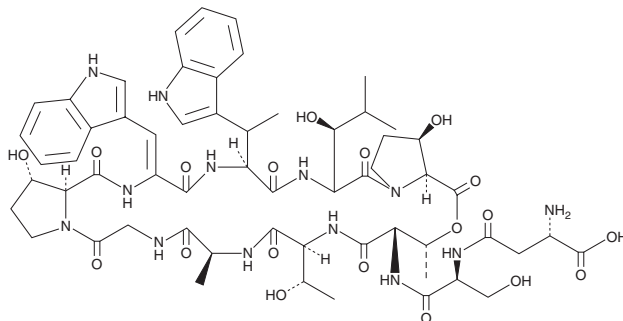
Purity: ≥95%

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

Telomycin is supplied as a solid. A stock solution may be made by dissolving the telomycin in the solvent of choice. Telomycin is soluble in organic solvents such as ethanol, methanol, DMSO, and dimethyl formamide, which should be purged with an inert gas. It is slightly soluble in water. We do not recommend storing the aqueous solution for more than one day.

### Description

Telomycin is a macrocyclic peptide lactone antibiotic originally isolated from *Streptomyces*.<sup>1</sup> It is active against Gram-positive bacteria, including *S. aureus*, methicillin-resistant *S. aureus* (MRSA), and *E. faecium* (MICs = 4, 2, and 16 µg/ml, respectively).<sup>2</sup> Telomycin eradicates *B. anserina* spirochetosis infection in chicks.<sup>3</sup>

### References

1. Sheehan, J.C., Mania, D., Nakamura, S., *et al.* The structure of telomycin. *J. Am. Chem. Soc.* **90**(2), 462-470 (1968).
2. Fu, C., Keller, L., Bauer, A., *et al.* Biosynthetic studies of telomycin reveal new lipopeptides with enhanced activity. *J. Am. Chem. Soc.* **137**(24), 7692-7705 (2015).
3. Packchanian, A. Chemotherapy of *Borrelia anserina* infection (spirochetosis) in chicks with amphotycin, kanamycin, telomycin, and tetracycline. *Antibiot. Chemother. (Northfield)* **10**, 731-739 (1960).

#### 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, 01/03/2023

#### 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