

# PRODUCT INFORMATION



## Teicoplanin A3-1

Item No. 20193

**CAS Registry No.:** 93616-27-4  
**Formal Name:** 34-O-[2-(acetylamino)-2-deoxy-β-D-glucopyranosyl]-22,31-dichloro-7-demethyl-64-O-demethyl-19-deoxy-42-O-α-D-mannopyranosyl-ristomycin A aglycone

**Synonym:** Antibiotic L 17054

**MF:** C<sub>72</sub>H<sub>68</sub>Cl<sub>2</sub>N<sub>8</sub>O<sub>28</sub>

**FW:** 1,564.3

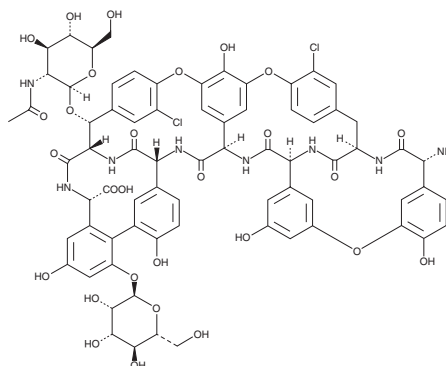
**Purity:** ≥98%

**Supplied as:** A solid

**Storage:** -20°C

**Stability:** ≥4 years

**Item Origin:** Bacterium/*Actinoplanes* sp.



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

### Laboratory Procedures

Teicoplanin A3-1 is supplied as a solid. A stock solution may be made by dissolving the teicoplanin A3-1 in the solvent of choice, which should be purged with an inert gas. Teicoplanin A3-1 is soluble in organic solvents such as ethanol, methanol, DMSO, and dimethyl formamide.

### Description

Teicoplanin A3-1 is a degradation product of teicoplanins A2-1 to 5 (Item Nos. 20187, 20188, 20189, 20192). Teicoplanins are glycopeptide antibiotics produced by *A. teichomyceticus* that are broadly effective against Gram-positive bacteria *in vitro*.<sup>1</sup>

### Reference

1. Somma, S., Gastaldo, L., and Corti, A. Teicoplanin, a new antibiotic from *Actinoplanes teichomyceticus* nov. sp. *Antimicrob. Agents Chemother.* **26(6)**, 917-923 (1984).

#### 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, 01/30/2023

#### 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