

# PRODUCT INFORMATION



## Teicoplanin A2-5

Item No. 20192

**CAS Registry No.:** 91032-38-1  
**Formal Name:** 34-O-[2-(acetylamino)-2-deoxy-β-D-glucopyranosyl]-22,31-dichloro-7-demethyl-64-O-demethyl-19-deoxy-56-O-[2-deoxy-2-[(9-methyl-1-oxodecyl)amino]-β-D-glucopyranosyl]-42-O-α-D-mannopyranosyl-ristomycin A aglycone

**MF:** C<sub>89</sub>H<sub>99</sub>Cl<sub>2</sub>N<sub>9</sub>O<sub>33</sub>

**FW:** 1,893.7

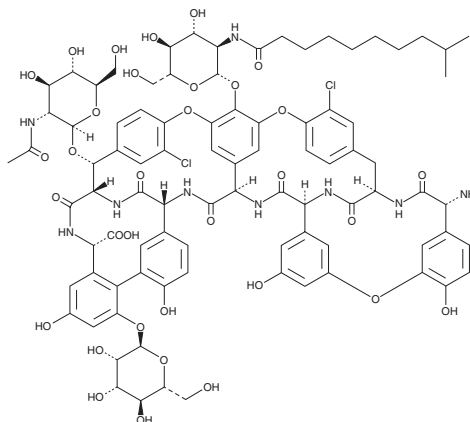
**Purity:** ≥98%

**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 A2-5 is supplied as a solid. A stock solution may be made by dissolving the teicoplanin A2-5 in the solvent of choice, which should be purged with an inert gas. Teicoplanin A2-5 is soluble in organic solvents such as ethanol, methanol, DMSO, and dimethyl formamide.

### Description

Teicoplanin A2-5 is a glycopeptide antibiotic produced by *A. teichomyceticus* that is broadly effective against Gram-positive bacteria *in vitro*.<sup>1</sup> It is the most non-polar metabolite in the family of teicoplanins.

### 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, 12/16/2022

#### 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](http://WWW.CAYMANCHEM.COM)