

PRODUCT INFORMATION



Teicoplanin A2-3

Item No. 20188

CAS Registry No.: 91032-36-9
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-[(1-oxodecyl)amino]-β-D-glucopyranosyl]-42-O-α-D-mannopyranosyl-ristomycin A aglycone

MF: C₈₈H₉₇Cl₂N₉O₃₃

FW: 1,879.7

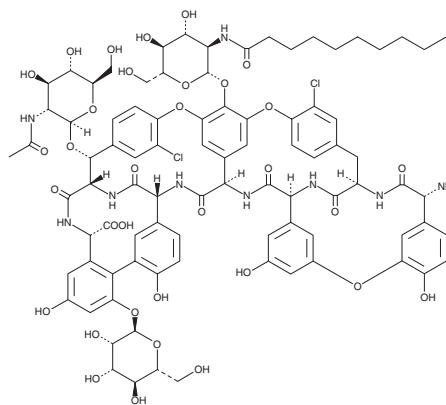
Purity: ≥85%

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 A2-3 is supplied as a solid. A stock solution may be made by dissolving the teicoplanin A2-3 in the solvent of choice, which should be purged with an inert gas. Teicoplanin A2-3 is soluble in ethanol, methanol, DMSO, and dimethyl formamide.

Description

Teicoplanin A2-3 is a major component of the teicoplanin complex (Item No. 15511), a glycopeptide antibiotic produced by *A. teichomyceticus* that is broadly effective against Gram-positive bacteria *in vitro*.¹

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, 12/12/2022

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