

PRODUCT INFORMATION



Teicoplanin Complex

Item No. 15511

CAS Registry No.: 61036-62-2

Synonyms: Antibiotic 8327A, Antibiotic MDL 507, MDL 507

MF: $C_{88}H_{95}Cl_2N_9O_{33}$ (for A2-1)

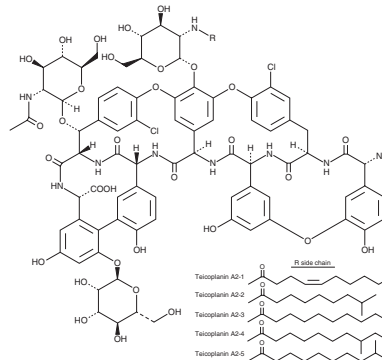
FW: 1,877.7 (for A2-1)

Purity: $\geq 95\%$

Supplied as: A crystalline solid

Storage: $-20^{\circ}C$

Stability: ≥ 4 years



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

Laboratory Procedures

Teicoplanin complex is supplied as a crystalline solid. A stock solution may be made by dissolving the teicoplanin complex in the solvent of choice. Teicoplanin complex is soluble in the organic solvent DMSO, which should be purged with an inert gas, at a concentration of approximately 0.15 mg/ml.

Further dilutions of the stock solution into aqueous buffers or isotonic saline should be made prior to performing biological experiments. Ensure that the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. Organic solvent-free aqueous solutions of teicoplanin complex can be prepared by directly dissolving the crystalline solid in aqueous buffers. The solubility of teicoplanin complex in PBS, pH 7.2, is approximately 0.25 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

Teicoplanin complex is a mixture of closely related metabolites produced by *A. teichomyceticus* that are effective against Gram-positive bacteria. Teicoplanin A2-2 (Item No. 16096), the major component of this complex, is a glycopeptide antibiotic structurally similar to vancomycin (Item No. 15327) that prevents bacterial cell-wall synthesis by binding and sequestering a cell-wall precursor with a D-alanine-containing peptide.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, 11/02/2022

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