

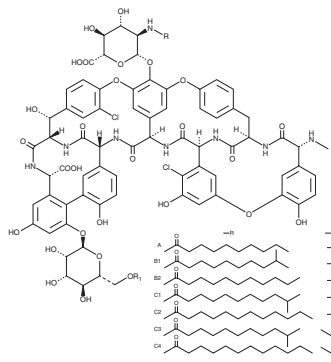
PRODUCT INFORMATION



Parvodicin Complex

Item No. 24036

CAS Registry No.: 187888-13-7
MF: $C_{83}H_{88}Cl_2N_8O_{29}$ (for C1)
FW: 1,732.5
Purity: $\geq 95\%$
Supplied as: A 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

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

Description

Parvodicin complex is a mixture of closely related lipophilic glycopeptides originally isolated from *A. parvosata*.¹ The parvodicin complex has activity against Gram-positive bacteria (MICs = 0.8-12.5 $\mu g/ml$) including methicillin-resistant strains of *S. aureus* (MRSA), *S. hemolyticus*, and *E. faecalis* (MICs = 0.1-25 $\mu g/ml$). Glycopeptide antibiotics act by preventing cell wall synthesis by binding and sequestering a cell wall precursor with a D-alanine-containing peptide.²

References

1. Christensen, S.B., Allaudeen, H.S., Burke, M.R., *et al.* Parvodicin, a novel glycopeptide from a new species, *Actinomadura parvosata*: Discovery, taxonomy, activity and structure elucidation. *J. Antibiot. (Tokyo)* **40(7)**, 970-990 (1987).
2. 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, 12/17/2022

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