

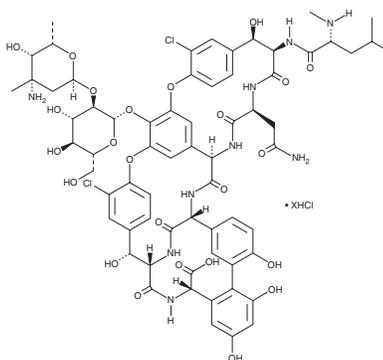
PRODUCT INFORMATION



Vancomycin (hydrochloride)

Item No. 15327

CAS Registry No.: 1404-93-9
Synonym: Lyphocin
MF: $C_{66}H_{75}Cl_2N_9O_{24} \cdot XHCl$
FW: 1,449.3
Purity: $\geq 98\%$
UV/Vis.: λ_{max} : 282 nm
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

Vancomycin (hydrochloride) is supplied as a crystalline solid. Aqueous solutions of vancomycin (hydrochloride) can be prepared by directly dissolving the crystalline solid in aqueous buffers. The solubility of vancomycin (hydrochloride) in PBS, pH 7.2, is approximately 5 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

Vancomycin is a glycopeptide antibiotic identified for its utility in the treatment of gram-positive bacteria, including penicillin-resistant staphylococci.¹ It acts by binding to bacterial cell wall protein precursors, interfering with continuing protein synthesis.² Vancomycin resistance in staphylococcal species has recently emerged as a clinical issue.³

References

1. Levine, D.P. Vancomycin: A history. *Clin. Infect. Dis.* **42(Suppl 1)**, S5-S12 (2006).
2. Williamson, M.P. and Williams, D.H. Hydrophobic interactions affect hydrogen bond strengths in complexes between peptides and vancomycin or ristocetin. *Eur. J. Biochem.* **138(2)**, 345-348 (1984).
3. Srinivasan, A., Dick, J.D., and Perl, T.M. Vancomycin resistance in staphylococci. *Clin. Microbiol. Rev.* **15(3)**, 430-438 (2002).

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, 10/03/2022

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