

PRODUCT INFORMATION



Kanamycin B

Item No. 34561

CAS Registry No.: 4696-76-8
Formal Name: O-3-amino-3-deoxy- α -D-glucopyranosyl-(1 \rightarrow 6)-O-[2,6-diamino-2,6-dideoxy- α -D-glucopyranosyl-(1 \rightarrow 4)]-2-deoxy-D-streptamine

Synonyms: Bekanamycin, Kanendomycin, Nebramycin V

MF: C₁₈H₃₇N₅O₁₀

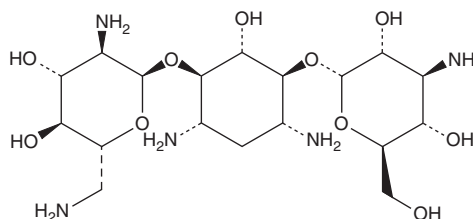
FW: 483.5

Purity: \geq 95%

Supplied as: A solid

Storage: -20°C

Stability: \geq 4 years



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

Laboratory Procedures

Kanamycin B is supplied as a solid. Aqueous solutions of kanamycin B can be prepared by directly dissolving the solid in aqueous buffers. The solubility of kanamycin B in PBS (pH 7.2) is approximately 1 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

Kanamycin B is an aminoglycoside antibiotic.¹ It is active against *S. aureus*, *E. coli*, *P. aeruginosa*, and *P. rettgeri* (MICs = 0.78-25 μ g/ml). Kanamycin B binds to the ribosomal decoding A site to interfere with bacterial protein synthesis.²

References

1. Umezawa, H., Umezawa, S., Tsuchiya, T., *et al.* 3',4'-Dideoxy-kanamycin B active against kanamycin-resistant *Escherichia coli* and *Pseudomonas aeruginosa*. *J. Antibiot. (Tokyo)* **24(7)**, 485-487 (1971).
2. Salian, S., Matt, T., Akbergenov, R., *et al.* Structure-activity relationships among the kanamycin aminoglycosides: Role of ring I hydroxyl and amino groups. *Antimicrob. Agents Chemother.* **56(12)**, 6104-6108 (2012).

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, 05/17/2023

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