

PRODUCT INFORMATION



Tobramycin (sulfate)

Item No. 41453

CAS Registry No.: 49842-07-1

Formal Name: O-3-amino-3-deoxy- α -D-glucopyranosyl-(1 $\rightarrow$ 6)-O-[2,6-diamino-2,3,6-trideoxy- α -D-ribohexopyranosyl-(1 $\rightarrow$ 4)]-2-deoxy-D-streptamine, sulfuric acid

Synonyms: Deoxykanamycin B, Distobram, Gernebcin

MF: $C_{18}H_{37}N_5O_9 \cdot 5/2H_2SO_4$

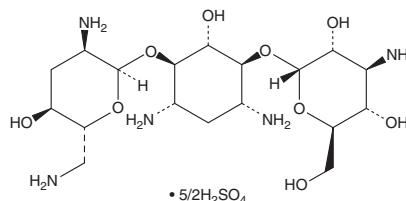
FW: 712.7

Purity: $\geq 98\%$

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

Tobramycin (sulfate) is supplied as a solid. Aqueous solutions of tobramycin (sulfate) can be prepared by directly dissolving the solid in aqueous buffers. Tobramycin (sulfate) is soluble (≥ 10 mg/ml) in PBS (pH 7.2). We do not recommend storing the aqueous solution for more than one day.

Description

Tobramycin is an aminoglycoside antibiotic.¹ It binds to the 30S ribosomal subunit, inhibiting protein synthesis. Tobramycin is active against a variety of Gram-positive and Gram-negative bacteria, including several strains of *S. aureus*, *P. aeruginosa*, *E. coli*, *K. pneumoniae*, and *Enterobacter* (mean MICs = 0.2, 0.6, 1.2, 0.8, and 0.8 μ g/ml, respectively). It reduces the number of colony-forming units (CFUs) in several mouse models of thigh infection.² Formulations containing tobramycin have been used in the treatment of Gram-positive and Gram-negative bacterial infections.

References

1. Neu, H.C. Tobramycin: An overview. *J. Infect. Dis.* **134**(Suppl. 1), S3-S19 (1976).
2. Fantin, B., Leggett, J., Ebert, S., *et al.* Correlation between in vitro and in vivo activity of antimicrobial agents against gram-negative bacilli in a murine infection model. *Antimicrob. Agents Chemother.* **35**(7), 1413-1422 (1991).

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, 04/18/2025

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