

PRODUCT INFORMATION



Dihydrostreptomycin (sulfate)

Item No. 27879

CAS Registry No.: 5490-27-7
Formal Name: O-2-deoxy-2-(methylamino)- α -L-glucopyranosyl-(1 \rightarrow 2)-O-5-deoxy-3-C-(hydroxymethyl)- α -L-lyxofuranosyl-(1 \rightarrow 4)-N¹,N³-bis(aminoiminomethyl)-D-streptamine, sulfate (2:3)

MF: C₂₁H₄₁N₇O₁₂ • 1.5H₂SO₄
FW: 730.7

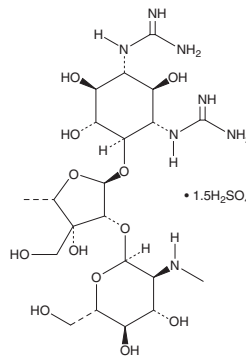
Purity: \geq 95%

Supplied as: A crystalline 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

Dihydrostreptomycin (sulfate) is supplied as a crystalline solid. Aqueous solutions of dihydrostreptomycin (sulfate) can be prepared by directly dissolving the crystalline solid in aqueous buffers. The solubility of dihydrostreptomycin (sulfate) in PBS, pH 7.2, is approximately 10 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

Dihydrostreptomycin is an aminoglycoside antibiotic and a derivative of streptomycin (Item No. 21211).¹ It is active against *B. subtilis*, *K. pneumoniae*, *S. marcescens*, *S. aureus*, *P. aeruginosa*, and *M. tuberculosis* (MICs = 0.07, 0.08, 0.09, 0.06, 0.8, and 5.7 μ g/ml, respectively). It interferes with bacterial protein synthesis by binding the 30S ribosomal subunit.² Dihydrostreptomycin enhances phagocytosis and intracellular killing of *E. coli* by isolated mouse peritoneal macrophages.³ It improves survival in a mouse model of *S. aureus* infection.⁴ Formulations containing dihydrostreptomycin have been used in the prevention and treatment of bacterial infections in livestock.

References

1. Donovan, R. and Rake, G. Studies on some biological aspects of dihydrostreptomycin. *J. Bacteriol.* **53**(2), 205-211 (1947).
2. Kaji, H. and Tanaka, Y. Binding of dihydrostreptomycin to ribosomal subunits. *J. Mol. Biol.* **32**(2), 221-230 (1968).
3. Adam, D., Staber, F., Belohradsky, B.H., et al. Effect of dihydrostreptomycin on phagocytosis of mouse-peritoneal macrophages in vitro. *Infect. Immun.* **5**(4), 537-541 (1972).
4. Waksman, S.A., Frankel, J., and Graessle, O. The in vivo activity of neomycin. *J. Bacteriol.* **58**(2), 229-237 (1949).

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, 11/01/2022

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