

PRODUCT INFORMATION



Streptomycin (sulfate)

Item No. 21211

CAS Registry No.: 3810-74-0
Formal Name: O-2-deoxy-2-(methylamino)- α -L-glucopyranosyl-(1 \rightarrow 2)-O-5-deoxy-3-C-formyl- α -L-lyxofuranosyl-(1 \rightarrow 4)-N¹,N³-bis(aminoiminomethyl)-D-streptamine

MF: C₂₁H₃₉N₇O₁₂ • 1.5H₂SO₄

FW: 728.7

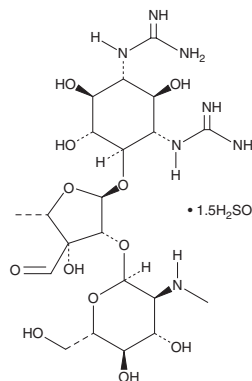
Purity: \geq 95%

UV/Vis.: λ_{max} : 231 nm

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

Streptomycin (sulfate) is supplied as a crystalline solid. A stock solution may be made by dissolving the streptomycin (sulfate) in the solvent of choice, which should be purged with an inert gas. Streptomycin (sulfate) is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of streptomycin (sulfate) in these solvents is approximately 2 and 13 mg/ml, respectively.

Further dilutions of the stock solution into aqueous buffers or isotonic saline should be made prior to performing biological experiments. Ensure that the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. Organic solvent-free aqueous solutions of streptomycin (sulfate) can be prepared by directly dissolving the crystalline solid in aqueous buffers. The solubility of streptomycin (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

Streptomycin is an aminoglycoside antibiotic that inhibits growth of both Gram-negative (MICs = 1 and 16 μ g/mL for *K. pneumoniae* and *S. marcescens*, respectively) and Gram-positive bacteria (MIC = 0.25 μ g/mL for *S. pneumoniae*).¹ It inhibits growth of *M. tuberculosis* H37Rv (MIC = 5 μ g/mL) and of susceptible strains from clinical isolates (MICs = \leq 0.125-1 μ g/mL).^{2,3} Streptomycin (150 mg/kg), in combination with isoniazid (Item No. 20378), rifapentine (Item No. 20307), and moxifloxacin (Item No. 14830), administered once per week over 6 months to *M. tuberculosis*-inoculated mice results in clearance of the pathogen from the lungs of the majority of treated mice.⁴ However, after 3 months without treatment, 58% of treated mice produce a positive lung culture. Streptomycin acts by inhibiting protein synthesis in prokaryotes by binding to the 30S ribosomal subunit.⁵⁻⁸ It has been used, in combination with penicillin G (Item No. 21615), in antibiotic cocktails to prevent bacterial growth in cell culture.⁹ Formulations containing streptomycin in combination with other antibiotics have been used to treat tuberculosis.

References

1. Syed, H.C., and Ravaoarino, M. LiF reduces MICs of antibiotics against clinical isolates of gram-positive and gram-negative bacteria. *Int. J. Microbiol.* **454065** (2012).
2. Jhamb, S.S., Goyal, A., and Singh, P.P. Determination of the activity of standard anti-tuberculosis drugs against intramacrophage *Mycobacterium tuberculosis*, *in vitro*: MGIT 960 as a viable alternative for BACTEC 460. *Braz. J. Infect. Dis.* **18(3)**, 336-340 (2014).

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, 12/01/2022

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

PRODUCT INFORMATION



3. Franzblau, S.G., Witzig, R.S., McLaughlin, J.C., *et al.* Rapid, low-technology MIC determination with clinical *Mycobacterium tuberculosis* isolates by using the microplate Alamar Blue assay. *J. Clin. Microbiol.* **36(2)**, 362-366 (1998).
4. Lounis, N., Bentoucha, A., Truffot-Pernot, C., *et al.* Effectiveness of once-weekly rifapentine and moxifloxacin regimens against *Mycobacterium tuberculosis* in mice. *Antimicrob. Agents Chemother.* **45(12)**, 3482-3486 (2001).
5. Luzzatto, L., Apirion, D., and Schlessinger, D. Mechanism of action of streptomycin in *E. coli*: Interruption of the ribosome cycle at the initiation of protein synthesis. *Proc. Natl. Acad. Sci. U.S.A.* **60(3)**, 873-880 (1968).
6. Moazed, D., and Noller, H.F. Interaction of antibiotics with functional sites in 16S ribosomal RNA. *Nature* **327(6121)**, 389-394 (1987).
7. Carter, A.P., Clemons, W.M., Brodersen, D.E., *et al.* Functional insights from the structure of the 30S ribosomal subunit and its interactions with antibiotics. *Nature* **407(6802)**, 340-348 (2000).
8. Sharma, D., Cukras, A.R., Rogers, E.J., *et al.* Mutational analysis of S12 protein and implications for the accuracy of decoding by the ribosome. *J. Mol. Biol.* **374(4)**, 1065-1076 (2007).
9. Varghese, D.S., Parween, S., Ardah, M.T., *et al.* Effects of aminoglycoside antibiotics on human embryonic stem cell viability during differentiation in vitro. *Stem Cells Int.* **2451927** (2017).

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