

PRODUCT INFORMATION

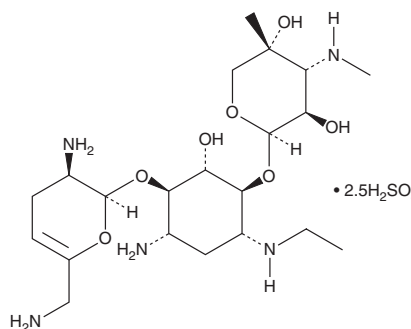


Netilmicin (sulfate)

Item No. 26666

CAS Registry No.: 56391-57-2
Formal Name: O-3-deoxy-4-C-methyl-3-(methylamino)-β-L-arabinopyranosyl-(1→6)-O-[2,6-diamino-2,3,4,6-tetradeoxy-α-D-glycero-hex-4-enopyranosyl-(1→4)]-2-deoxy-N¹-ethyl-D-streptomine, sulfate (2:5)

Synonyms: Netillin, Netilyn
MF: C₂₁H₄₁N₅O₇ • 2.5H₂SO₄
FW: 720.8
Supplied as: A crystalline solid
Storage: -20°C
Stability: ≥4 years



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

Laboratory Procedures

Netilmicin (sulfate) is supplied as a crystalline solid. Aqueous solutions of netilmicin (sulfate) can be prepared by directly dissolving the crystalline solid in aqueous buffers. The solubility of netilmicin (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

Netilmicin is an aminoglycoside antibiotic.¹ It is active against *S. aureus*, *N. gonorrhoeae*, *E. coli*, *P. mirabilis*, and *P. aeruginosa* (MICs = 0.4, 3.1, 0.8, 1.6, and 0.4 µg/ml, respectively) as well as *Enterococci*, *Enterobacter*, *Citrobacter*, and *Klebsiella* species (MICs = 3.1, 0.4, 0.8, and ≤0.2 µg/ml, respectively). Netilmicin induces nephrotoxicity in rats when administered at doses ranging from 50 to 150 mg/kg.² It also induces neuromuscular block, decreasing twitch tension in isolated diaphragm and soleus muscles in rabbits (ED₅₀s = 18.5 and 62.2 mg/kg, respectively).³ Formulations containing netilmicin were previously used in the intravenous treatment of severe bacterial infections.

References

1. Sanders, C.C., Sanders W.E., J., and Goering, R.V. In vitro studies with Sch 21420 and Sch 22591: Activity in comparison with six other aminoglycosides and synergy with penicillin against enterococci. *Antimicrob. Agents Chemother.* **14(2)**, 178-184 (1978).
2. Bowman, R.L., Silverblatt, F.J., and Kaloyanides, G.J. Comparison of the nephrotoxicity of netilmicin and gentamicin in rats. *Antimicrob. Agents Chemother.* **12(4)**, 474-478 (1977).
3. Liu, M., Kato, M., and Hashimoto, Y. Neuromuscular blocking effects of the aminoglycoside antibiotics arbekacin, astromicin, isepamicin and netilmicin on the diaphragm and limb muscles in the rabbit. *Pharmacology* **63(3)**, 142-146 (2001).

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, 10/21/2022

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