

PRODUCT INFORMATION



Apramycin (sulfate hydrate)

Item No. 19688

Formal Name: O-4-amino-4-deoxy- α -D-glucopyranosyl-(1 \rightarrow 8)-O-(8R)-2-amino-2,3,7-trideoxy-7-(methylamino)-D-glycero- α -D-allo-octodialdo-1,5:8,4-dipyranosyl-(1 \rightarrow 4)-2-deoxy-D-streptamine, sulfate, hydrate

Synonym: Nebramycin II

MF: C₂₁H₄₁N₅O₁₁ • XH₂SO₄ [XH₂O]

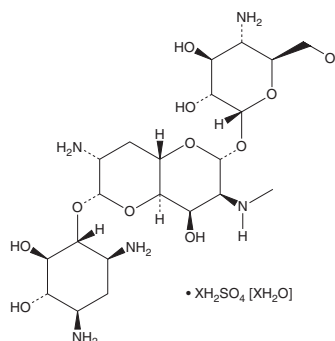
FW: 539.6

Purity: \geq 95%

Supplied as: A powder

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

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

Description

Apramycin is an aminoglycoside antibiotic first isolated from *S. tenebrarius*.¹ It is active against a variety of Gram-positive and Gram-negative bacteria, including strains which produce aminoglycoside-modifying enzymes.² Apramycin inhibits protein synthesis in bacteria both *in vivo* and *in vitro*.¹

References

1. Abe, Y., Nakagawa, S., Naito, T., *et al.* Aminoglycoside antibiotics. XIV. Synthesis and activity of 6-O-(3-amino-3-deoxy- α -D-glucopyranosyl)-and 5-O-(β -D-ribofuranosyl)apramycins. *J. Antibiot. (Tokyo)* **34(11)**, 1434-1446 (1981).
2. Perzynski, S., Cannon, M., Cundliffe, E., *et al.* Effects of apramycin, a novel aminoglycoside antibiotic on bacterial protein synthesis. *Eur. J. Biochem.* **99(3)**, 623-628 (1979).

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, 11/29/2022

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