

PRODUCT INFORMATION

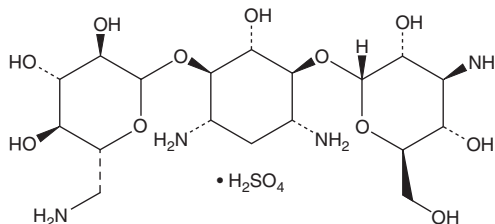


Kanamycin A (sulfate)

Item No. 15321

CAS Registry No.: 25389-94-0
Formal Name: O-3-amino-3-deoxy- α -D-glucopyranosyl-(1 \rightarrow 6)-O-[6-amino-6-deoxy- α -D-glucopyranosyl-(1 \rightarrow 4)]-2-deoxy-D-streptomine, monosulfate

Synonym: Ophthalmokalixan
MF: C₁₈H₃₆N₄O₁₁ • H₂SO₄
FW: 582.6
Purity: \geq 90%
Supplied as: A crystalline solid
Storage: -20°C
Stability: \geq 4 years
Item Origin: Synthetic



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

Laboratory Procedures

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

Kanamycin A is a broad spectrum antibiotic isolated from a soil streptomycetes (*S. kanamyceticus*). This aminoglycoside alters translation by prokaryotic ribosomes, resulting in bacterial cell death.¹ In molecular biology, kanamycin A is used as a selective agent to isolate transformants that are expressing a kanamycin resistance gene, usually neomycin phosphotransferase (NptII/neo). This mode of selection is most commonly used in plant and mammalian biology.

Reference

1. Misumi, M. and Tanaka, N. Mechanism of inhibition of translocation by kanamycin and viomycin: A comparative study with fusidic acid. *Biochem. Biophys. Res. Commun.* **92**(2), 647-654 (1980).

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, 11/11/2022

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