

Kanamycin . sulfate

Antibiotic

Aminoglycoside antibiotic with broad antibacterial spectrum, including mycobacteria, many Gram-positive and most Gram-negative bacteria. Inhibits bacterial protein synthesis. Kanamycin resistance is widely used as a selection marker for obtaining transgenic plants.

Citations: 4

[View Online »](#)

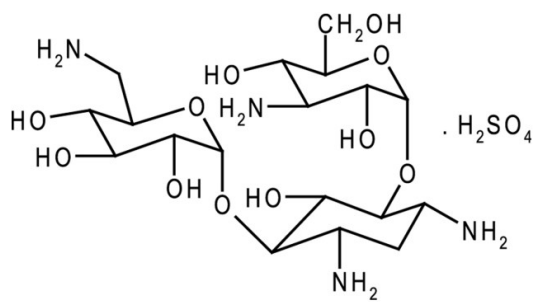
Ordering Information

[Order Online »](#)

ALX-380-049-G005	5g
------------------	----

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability As indicated on product label or CoA when stored as recommended.

Long Term Storage +4°C

Shipping Ambient Temperature

Regulatory Status RUO - Research Use Only

Product Details

Activity ≥700 µg/mg

Appearance White to off-white crystalline powder.

CAS 25389-94-0

Couple Type Inhibitor

Formula $C_{18}H_{36}N_4O_{11} \cdot H_2SO_4$

MI 14: 5281

MW 484.5 . 98.1

Solubility Soluble in water.

Source Isolated from *Streptomyces kanamyceticus*.

Last modified: May 29, 2024



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
[info-
usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
[info-
eu@enzolifesciences.com](mailto:info-eu@enzolifesciences.com)

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
[info-
be@enzolifesciences.com](mailto:info-be@enzolifesciences.com)

France
Phone: +33 472 440 655
[info-
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany
Phone: +49 7621 5500 526
[info-
de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
[info-
uk@enzolifesciences.com](mailto:info-uk@enzolifesciences.com)