

Gentamicin sulfate

Broad-spectrum antibiotic

Broad-spectrum cell culture antibiotic that is nontoxic to viruses and mammalian cells at antibacterial and antimycoplasmal concentrations. Due to its extended stability and slow development of bacterial resistance, it is a useful antibiotic in long-term virus and tissue culture studies.

Citations: 6

[View Online »](#)

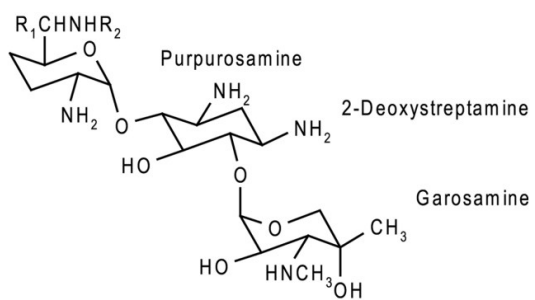
Ordering Information

[Order Online »](#)

ALX-380-003-G005	5g
ALX-380-003-G010	10g
ALX-380-003-G025	25g

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended.
Handling	Hygroscopic.
Long Term Storage	+4°C
Shipping	Ambient Temperature

Regulatory Status

RUO - Research Use Only

Product Details

Activity	Potency is ≥ 590 IU/mg (anhydrous material).
Appearance	White to off-white powder.
CAS	1405-41-0
Formula	$C_{63}H_{131}N_{15}O_{25}S$
MI	14: 4392
MW	1530.9
RTECS	LY2625000
Solubility	Soluble in water; almost insoluble in alcohol or ether.
Source	Isolated from <i>Micromonospora</i> sp.

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-
fr@enzolifesciences.com](mailto:info-fr@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)