

Neomycin . sulfate

Inhibitor of PLC and PLD

Aminoglycoside antibiotic that inhibits translation by binding to the small subunit of prokaryotic ribosomes. Inhibitor of phospholipase C (PLC; IC_{50} =1-10 μ M) due to binding to inositol phospholipids. Neomycin inhibits phosphatidylcholine-phospholipase D activity (IC_{50} =65 μ M). Blocks voltage sensitive Ca^{2+} channels without affecting Na^+/Ca^{2+} antiporters in nerve cells. Induces PLA2 activation in human platelets.

Citations: 16

[View Online »](#)

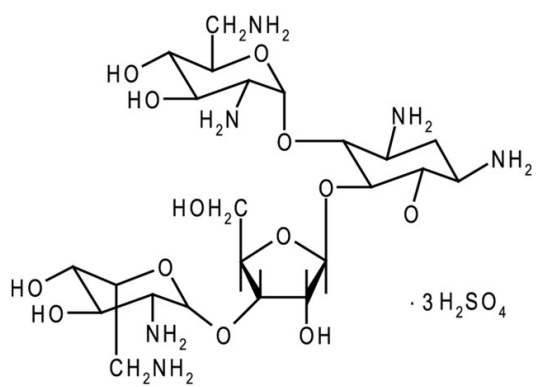
Ordering Information

[Order Online »](#)

ALX-380-035-G025	25g
------------------	-----

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended.
Handling	Protect from light and moisture.
Long Term Storage	-20°C
Shipping	Ambient Temperature

Regulatory Status

RUO - Research Use Only

Product Details

Appearance	White to off-white solid.
CAS	1405-10-3
Couple Target	Calcium channel, PLC, PLD
Couple Type	Blocker, Inhibitor
Formula	$C_{23}H_{46}N_6O_{13} \cdot 3H_2SO_4$
MI	14: 6454
MW	614.7 . 294.2
Purity	≥90% (TLC)
RTECS	QP4375000
Solubility	Soluble in water (10mg/ml).

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)