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 $_{50}$ =1-10µM) due to binding to inositol phospholipids. Neomycin inhibits phosphatidylcholine-phospholipase D activity (IC $_{50}$ =65µM). 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

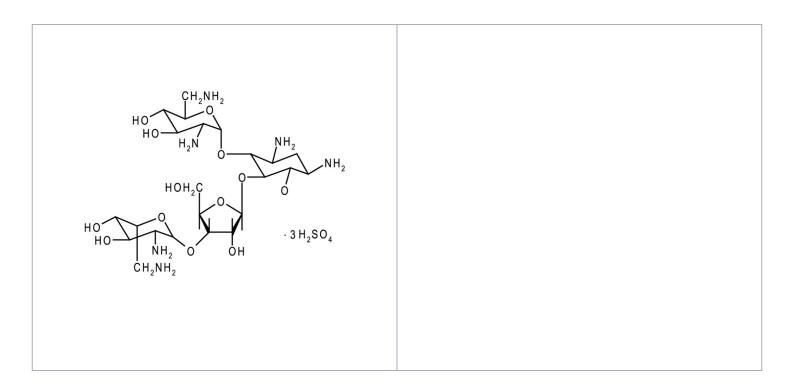
Order Online »

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 $C_{23}H_{46}N_6O_{13}$. $3H_2SO_4$

MI 14: 6454

MW 614.7 . 294.2

Purity ≥90% (TLC)

RTECS QP4375000

Solubility Soluble in water (10mg/ml).

Last modified: May 29, 2024

