

Vancomycin . hydrochloride

Antibiotic

Glycopeptide antibiotic. Interferes with cell wall synthesis. Effective against Gram-positive bacteria.

Citations: 3

[View Online »](#)

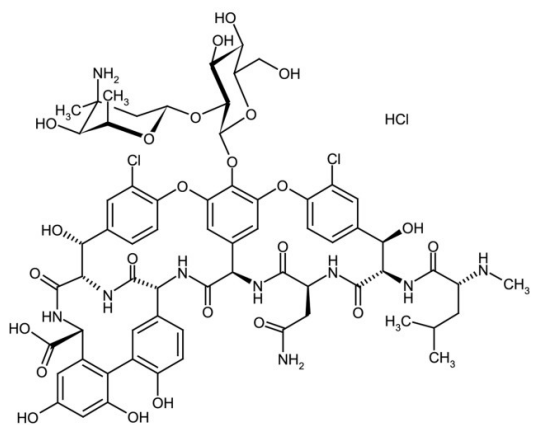
Ordering Information

[Order Online »](#)

ALX-380-279-M250	250mg
ALX-380-279-G001	1g

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

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

Long Term Storage -20°C

Shipping Ambient Temperature

Regulatory Status RUO - Research Use Only

Product Details

Activity Potency is >900µg/mg (on dry basis).

Appearance White to light brown powder.

CAS 1404-93-9

Formula $C_{66}H_{75}Cl_2N_9O_{24} \cdot HCl$

MI **14:** 9929

MW 1449.3 . 36.5

RTECS YW4380000

Solubility Soluble in water (50mg/ml).

Source Isolated from *Streptomyces lysosup*.

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)