

Colistin . sulfate

Antibiotic

Polymyxin antibiotic. Binds to lipids on the cell cytoplasmic membrane. Disrupts cell wall integrity. Effective against Gram-negative bacteria.

Citations: 4

[View Online »](#)

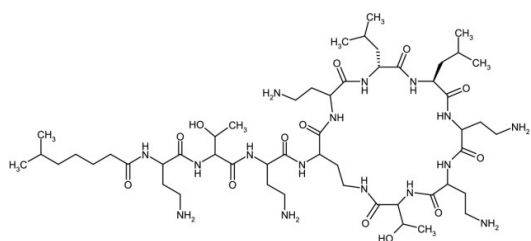
Ordering Information

[Order Online »](#)

ALX-380-272-M500	500mg
------------------	-------

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 >19'900 u/mg (on dry basis).

Alternative Name Polymyxin E . sulfate

Appearance White to off-white powder.

CAS 1264-72-8

MI **14:** 2479

Purity Detail Mixture of >30 compounds, primarily colistin A and B.

RTECS TR1500000

Solubility Soluble in water (50mg/ml).

Source Isolated from *Bacillus colistinus*.

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)