

Polymyxin B . sulfate

PKC inhibitor

Antibiotic. Inhibitor of protein kinase C (PKC). Breaks the bacterial membrane by incorporating into the phospholipid of the outer membrane and activating phospholipase.

Citations: 10

[View Online »](#)

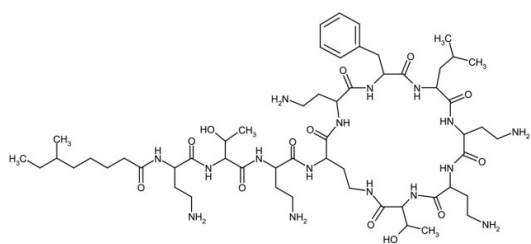
Ordering Information

[Order Online »](#)

ALX-380-040-G001	1g
ALX-380-040-G005	5g

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

Appearance	White to off-white powder.
CAS	1405-20-5
Couple Target	Phospholipase, PKC
Couple Type	Activator, Inhibitor
MI	14: 7573
Purity	≥80% (HPLC) (sum of polymyxins B1, B2, B3, B1-I)
Purity Detail	Mixture of Polymyxin B1, B2, B3 and B1-I.
RTECS	TR1150000
Solubility	Soluble in water (50mg/ml).
Source	Synthetic.

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)