

Virginiamycin M1

Streptogramin antibiotic

Antibiotic that inhibits bacterial protein synthesis.

Virginiamycin M1 inhibits bacterial protein synthesis at the level of aminoacyl-tRNA binding and peptide bond formation. It inactivates the 50S ribosome. When used in combination with Virginiamycin S1 it is synergistic. Virginiamycin is used as a performance promoter in animal husbandry. It is chemically modified to make therapeutic drugs such as quinupristin and dalfopristin.

Citations: 1

[View Online »](#)

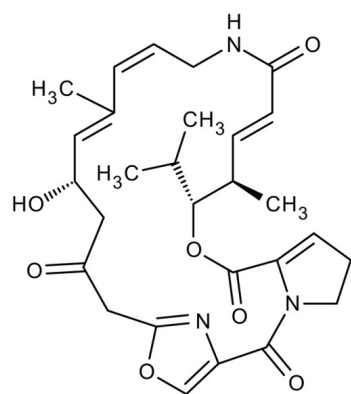
Ordering Information

[Order Online »](#)

ENZ-CHM105-0005	5mg
-----------------	-----

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 Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name Mikamycin A, Staphilomycin, Ostreogrycin A

Appearance White solid.

CAS 21411-53-0

Formula $C_{28}H_{35}N_3O_7$

MW 525.6

Purity ≥95% (HPLC)

Source Isolated from *Streptomyces virginiae*



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-
be@enzolifesciences.com](mailto:info-be@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)