

Spectinomycin . sulfate

Aminocyclitol antibiotic

Antibiotic. Inhibits protein synthesis (elongation) by interfering with peptidyl tRNA translocation. Effective against Gram-negative and Gram-positive bacteria (*Gonococcus* only). Mutation in *rpsE* (the gene for ribosomal protein S5) prevents binding of spectinomycin.

Citations: 2

[View Online »](#)

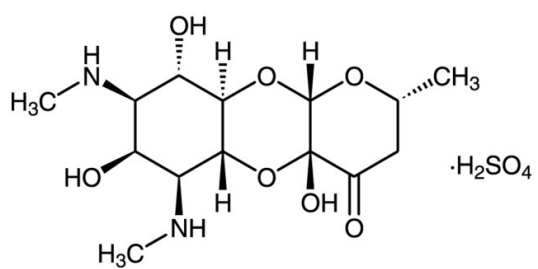
Ordering Information

[Order Online »](#)

BML-A281-0010	10g
---------------	-----

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

Alternative Name U-18409E

Appearance White solid.

CAS 64058-48-6

Formula $C_{14}H_{24}N_2O_7 \cdot H_2SO_4$

MW 430.4

Purity ≥94% (HPLC)

Solubility Soluble in water.

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-
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)