

Streptothricin . sulfate

Antibiotic

Antibiotic. Active against both Gram-positive and Gram-negative bacteria and some fungi.

Citations: 2

[View Online »](#)

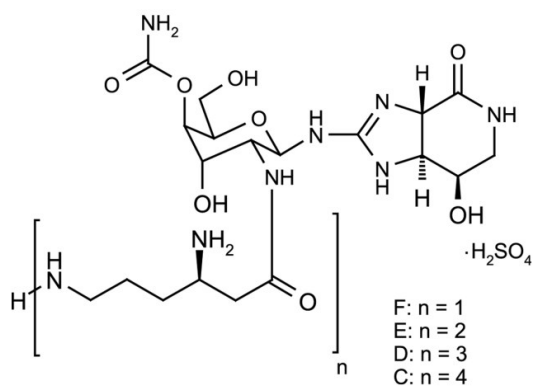
Ordering Information

[Order Online »](#)

ALX-380-025-M005	5mg
------------------	-----

Manuals, SDS & CofA

[View Online »](#)



F: $n = 1$
E: $n = 2$
D: $n = 3$
C: $n = 4$

Handling & Storage

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

Long Term Storage +4°C

Shipping Ambient Temperature

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name Nourseothricin . sulfate

Appearance White to off-white powder.

CAS 96736-11-7

Formula Streptothricin D: $C_{31}H_{58}N_{12}O_{10}$ Streptothricin F: $C_{19}H_{34}N_8O_8$

MI 14: 8829

MW Streptothricin D: 758.9 Streptothricin F: 502.5

Purity Detail Mixture of streptothricin F (n=1) and D (n=3).

RTECS RD4240000

Solubility Soluble in water.

Source Isolated from *Streptomyces noursei*.

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-be@enzolifesciences.com)

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
[info-
be@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)