

Ristocetin A . sulfate

Antibiotic. Inducer of platelet aggregation

Potent antibacterial antibiotic. Induces platelet aggregation by binding to a factor absent in people suffering from von Willebrand's disease and is therefore an important diagnostic tool. Chiral selector for the separation of enantiomers by HPLC, TLC and electrophoresis.

Citations: 5

[View Online »](#)

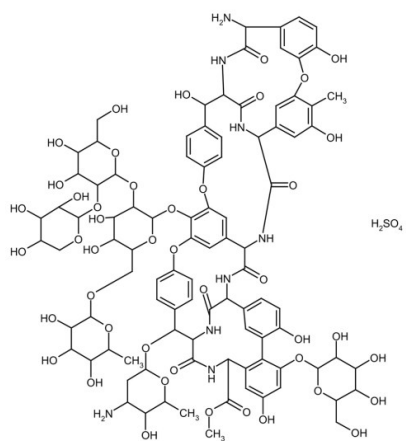
Ordering Information

[Order Online »](#)

ALX-380-237-M010	10mg
------------------	------

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 Ristocetin . sulfate, Ristomycin III . sulfate, Spontin . sulfate, Riston . sulfate

Appearance Light tan solid.

CAS 11140-99-1

Formula $C_{95}H_{110}N_8O_{44} \cdot H_2SO_4$

MI **14:** 8234

MW 2067.9 . 98.1

Purity ≥95% (HPLC)

RTECS VJ8595000

Solubility Soluble in water, methanol or DMSO.

Source Isolated from *Amycolatopsis* sp. MST-AS5924.

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)