

Venturicidin A

Inhibitor of mitochondrial ATP synthase

Macrolide antibiotic. Potent inhibitor of mitochondrial ATP synthase complex, acting on the F_0 membrane sector. Shows antifungal activity.

Citations: 5

[View Online »](#)

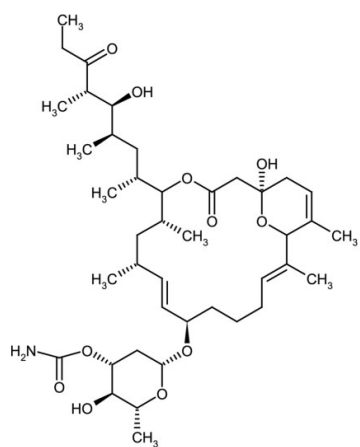
Ordering Information

[Order Online »](#)

ALX-380-211-M001	1mg
------------------	-----

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 Aabomycin A1

Appearance White to off-white solid.

CAS 33538-71-5

Couple Target ATPase

Couple Type Inhibitor

Formula $C_{41}H_{67}NO_{11}$

Formulation Lyophilized.

MI 14: 9947

MW 750.0

Purity ≥95% (HPLC)

RTECS YX4556000

Solubility Soluble in 100% ethanol, methanol, dimethyl formamide or DMSO.

Source Isolated from *Streptomyces* sp. MST-117594.



ENZO LIFE SCIENCES,
INC.
Phone: +1 909 424 1100
info-usa@enzolifesciences.com

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
info-eu@enzolifesciences.com

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
info-be@enzolifesciences.com

France
Phone: +33 472 440 655
info-fr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info-de@enzolifesciences.com

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
info-uk@enzolifesciences.com