

Elaiophylin

Inhibitor of testosterone 5-reductase and P-ATPase

Macrolide antibiotic inhibiting testosterone 5-reductase and plasma-proton ATPase (P-ATPase). Shows antifungal, antiprotozoal, antitumor, antihelminthic and immunosuppressive activity.

Citations: 9

[View Online »](#)

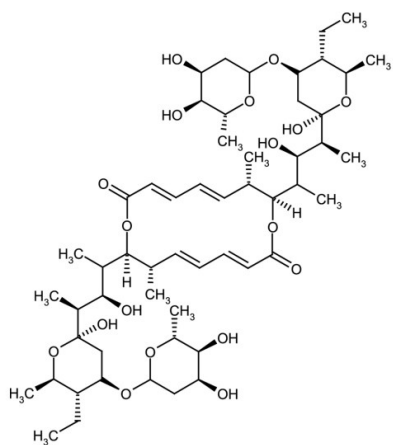
Ordering Information

[Order Online »](#)

ALX-380-212-M005	5mg
------------------	-----

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended. Stable for at least 1 year after receipt when stored at -20°C.
Handling	Protect from light when in solution.
Short Term Storage	+4°C
Long Term Storage	-20°C
Shipping	Ambient Temperature

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	Antibiotic 255-E, Antibiotic 5001B, Antibiotic 56-62, Antibiotic 846 I, SNA 4606-3, Azalomycin B, Gopalamicin, Salbomycin
Appearance	White to off-white solid.
CAS	37318-06-2
Couple Target	ATPase, Testosterone 5-reductase
Couple Type	Inhibitor
Formula	C ₅₄ H ₈₈ O ₁₈
Identity	Determined by 1H-NMR.
MW	1025.3
Purity	≥95% (HPLC)
RTECS	CL8380000
Solubility	Soluble in 100% ethanol, methanol, dimethyl formamide or DMSO.
Source	Isolated from <i>Streptomyces</i> sp. MT-1.



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)