

Oligomycin B

F-ATPase inhibitor

Minor component of the oligomycin complex. Macrolide antibiotic that inhibits membrane bound mitochondrial ATPase (F_1F_0), preventing phosphoryl group transfer. Induces apoptosis.

Citations: 4

[View Online »](#)

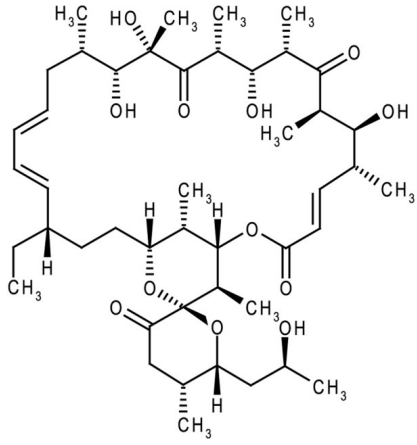
Ordering Information

[Order Online »](#)

ALX-380-038-M005	5mg
ALX-380-038-46190	5mg

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 28-Oxooligomycin A

Appearance White to off-white powder.

CAS 11050-94-5

Couple Target ATPase

Couple Type Inhibitor

Formula $C_{45}H_{72}O_{12}$

MI 14: 6833

MW 805.1

Purity ≥95% (HPLC)

RTECS RK3330000

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

Source Isolated from *Streptomyces* sp. MST-AS5339.

Last modified: May 29, 2024



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
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