

Concanamycin A

H⁺ ATPase inhibitor

Antibiotic. More potent and specific H⁺-ATPase inhibitor than bafilomycin A₁ (Prod. No. BML-CM110). Inhibits acidification of organelles such as lysosomes and the Golgi apparatus. Blocks cell surface expression of viral glycoproteins without affecting their synthesis. Exhibits cytotoxic effects in a number of cell lines in a cell viability assay. Induces nitric oxide (NO) production. Inhibits Autophagy.

Citations: 24

[View Online »](#)

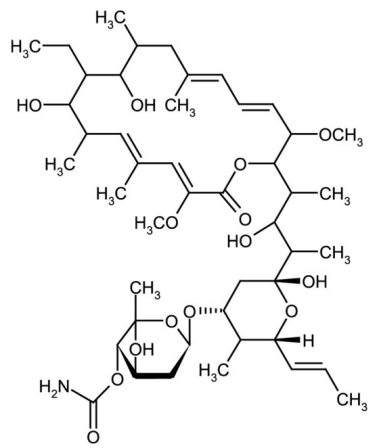
Ordering Information

[Order Online »](#)

ALX-380-034-C025	25µg
ALX-380-034-C100	100µg

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. After reconstitution protect from light . Solutions are more stable when stored at -20°C.
Short Term Storage	+4°C
Long Term Storage	-20°C
Shipping	Ambient Temperature

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name	Folimycin
Appearance	White to off-white solid (sub-milligram quantities will appear as clear film).
CAS	80890-47-7
Couple Target	ATPase
Couple Type	Inhibitor
Formula	$C_{46}H_{75}NO_{14}$
MW	866.1
Purity	≥95% (HPLC)
RTECS	CB9732000
Solubility	Soluble in methanol, DMSO or acetonitrile; insoluble in water.
Source	Isolated from <i>Streptomyces</i> sp.



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
Last modified: **May 29, 2024**
usa@enzolifesciences.com

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

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

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

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

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