

# Concanamycin A

## H<sup>+</sup> ATPase inhibitor

Antibiotic. More potent and specific H<sup>+</sup>-ATPase inhibitor than bafilomycin A<sub>1</sub> (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 »](#)

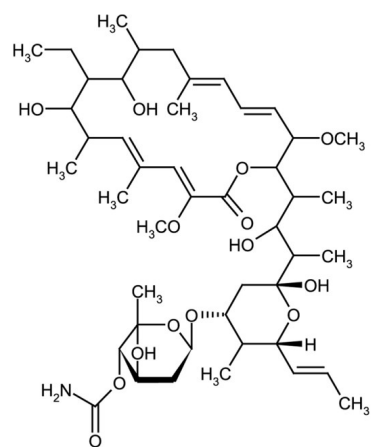
## Ordering Information

[Order Online »](#)

ALX-380-034-M001	1mg
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.



Last modified: May 29, 2024

ENZO LIFE SCIENCES,  
INC.  
Phone: 800.942.0430  
[usa@enzolifesciences.com](mailto: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-  
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

France  
Phone: +33 472 440 655  
[info-  
de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

Germany  
Phone: +49 7621 5500 526  
[info-  
uk@enzolifesciences.com](mailto:info-uk@enzolifesciences.com)

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