

# Bafilomycin B<sub>1</sub>

V-ATPase inhibitor

Macrolide antibiotic that acts as a specific inhibitor of vacuolar-type H<sup>+</sup>-ATPase.

Citations: 7

[View Online »](#)

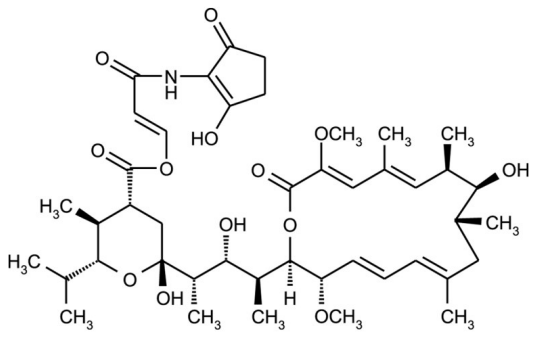
Ordering Information

[Order Online »](#)

ALX-380-063-M001	1mg
------------------	-----

Manuals, SDS & CofA

[View Online »](#)



## Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended.
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

Appearance	Yellow solid.
CAS	88899-56-3; 80112-34-1
Couple Target	ATPase
Couple Type	Inhibitor
Formula	$C_{44}H_{65}NO_{13}$
Identity	Determined by 1H-NMR.
MW	816.0
Purity	≥95% (HPLC)
Solubility	Slightly soluble in methanol or DMSO; insoluble in water.
Source	Isolated from <i>Streptomyces griseus</i> .

Last modified: May 29, 2024



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)