

# Bafilomycin A<sub>1</sub>

A selective V-ATPase inhibitor that disrupts lysosomal acidification, essential for autophagy flux analysis and lysosomal function studies.

Bafilomycin A1 is a macrolide antibiotic produced by *Streptomyces* with diverse biological activities. Bafilomycin A1 acts as a potent and selective inhibitor of vacuolar-type H<sup>+</sup>-ATPase (V-type) (IC<sub>50</sub>=0.04-0.4nmol/mg). Bafilomycin A1 is a valuable tool for distinguishing among different types of ATPases. Bafilomycin A1 inhibits endosomal acidification, lysosomal cholesterol trafficking, as well as autophagosome maturation and protein degradation. Bafilomycin A1 also blocks pH regulation in brain cells.

Citations: 193

[View Online »](#)

## Ordering Information

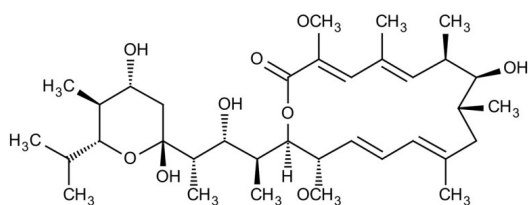
[Order Online »](#)

BML-CM110-0100	100µg
----------------	-------

## Manuals, SDS & CofA

[View Online »](#)

- Potent and selective inhibitor of vacuolar H<sup>+</sup>ATPases
- Compound used to study autophagy in a variety of cells
- Highly cited



## 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. Stock solutions are stable for up to 3 months at -20°C.
Handling	Protect from light.
Long Term Storage	-20°C
Shipping	Blue Ice

## Regulatory Status

RUO - Research Use Only

## Product Details

Alternative Name	NSC 381866
Appearance	White powder to colorless film.
CAS	88899-55-2
Couple Target	ATPase
Couple Type	Inhibitor
Formula	$C_{35}H_{58}O_9$
MW	622.8
Purity	≥95% (HPLC, TLC)
Solubility	Soluble in DMSO (5mg/ml), 100% ethanol, or 100% methanol (5mg/ml).
Source	<i>Streptomyces griseus</i>

## Technical Info / Product Notes

Typical concentrations for use in cell culture are 50-100nM.

Note: Product is not sterile.



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-  
fr@enzolifesciences.com](mailto:info-fr@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)

Last modified: May 29, 2024