

# Bafilomycin A1

Macrolide antibiotic; vacuolar H<sup>+</sup> ATPase inhibitor.

Macrolide isolated from *Streptomyces sp.* that has been shown to decrease multi-drug resistance. Specific inhibitor of vacuolar (V-type) proton ATPases. It also shows antibiotic properties and induces apoptosis.

## Ordering Information

[Order Online »](#)

LKT-B0025-MC01	0.1mg
LKT-B0025-MC05	0.5mg
LKT-B0025-M001	1mg
LKT-B0025-M005	5mg

## Manuals, SDS & CofA

[View Online »](#)

## Handling & Storage

Handling	Light sensitive
Long Term Storage	-20°C
Shipping	Ambient Temperature

**Regulatory Status** RUO - Research Use Only

## Product Details

CAS	88899-55-2
Couple Type	Inhibitor
Formula	$C_{35}H_{58}O_9$
MW	622.83
Purity	≥90%
Solubility	Soluble in DMSO, ethanol and methanol.



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)