

Bafilomycin A1

Macrolide antibiotic; vacuolar H⁺ 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-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

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

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

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

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

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