

2-Methoxy-antimycin

A

Bcl-2 inhibitor

An analog of antimycin A (mixture of antimycin A1, A2, A3 and A4) which binds to the hydrophobic groove of Bcl-2 and Bcl-X_L thereby inducing apoptosis. It displays selective cytotoxicity for Bcl-X_L-expressing cells (10µg/ml). Unlike antimycin A it does not exhibit general mitochondrial toxicity.

Citations: 3

[View Online »](#)

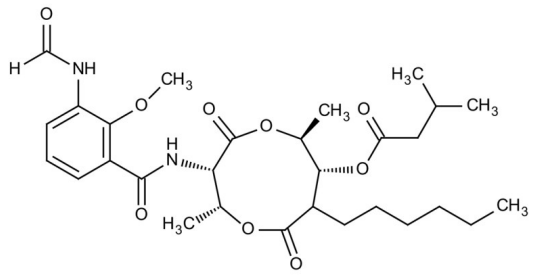
Ordering Information

[Order Online »](#)

BML-CM123-0001	1mg
----------------	-----

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended. Store, as supplied, at -20°C for up to 1 year. Store solutions at -20°C for up to 3 months.
Long Term Storage	-20°C
Shipping	Ambient Temperature

Regulatory Status

RUO - Research Use Only

Product Details

Appearance	White solid.
Couple Target	Bcl-2
Couple Type	Inhibitor
Formula	$C_{27}H_{38}N_2O_9$
MW	534.6
MeltingPoint	63.5-65°C
Purity	≥85%
Solubility	Soluble in DMSO (12 mg/ml) or ethanol (12 mg/ml)

Last modified: May 29, 2024



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