

Oligomycin A

ATPase inhibitor

Inhibits mitochondrial ATPases (F_1F_0). Stimulates lysosome acidification due to shunting of membrane potential. Induces apoptosis in a variety of cell types. Induces autophagy in the IPLB-LdFB insect cell line.

Citations: 11

[View Online »](#)

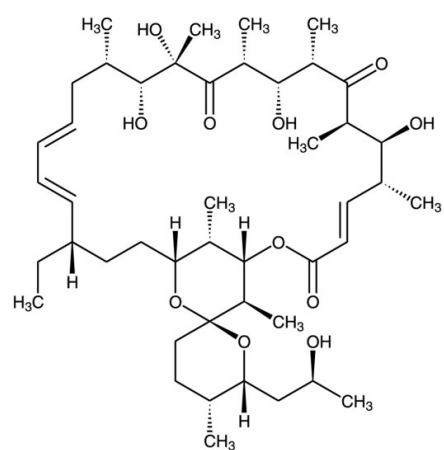
Ordering Information

[Order Online »](#)

BML-CM111-0005	5mg
BML-CM111-0025	25mg

Manuals, SDS & CofA

[View Online »](#)



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, as supplied, at -20°C. Stock solutions are stable for up to 3 months at -20°C.
Long Term Storage	-20°C
Shipping	Ambient Temperature

Regulatory Status

RUO - Research Use Only

Product Details

Appearance	White solid.
CAS	579-13-5
Couple Target	ATPase
Couple Type	Inhibitor
Formula	$C_{45}H_{74}O_{11}$
MI	14: 6833
MW	791.1
Purity	≥97% (HPLC, TLC)
RTECS	RK3328000
Solubility	Soluble in DMSO (300mg/ml), 100% ethanol (200mg/ml) or acetone (800mg/ml).

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