

Cantharidin

PP2A inhibitor

Potent protein phosphatase 2A (PP2A) inhibitor. At higher concentrations, also inhibits protein phosphatase 1 (PP1).

Citations: 5

[View Online »](#)

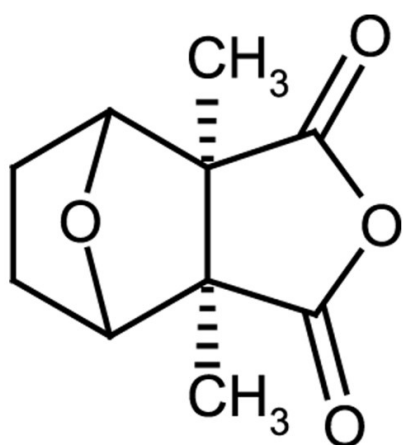
Ordering Information

[Order Online »](#)

ALX-270-063-M025	25mg
ALX-270-063-M100	100mg

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended.
Handling	Protect from light.
Long Term Storage	Ambient
Shipping	Ambient Temperature

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	Hexahydro-3a,7a-dimethyl-4,7-epoxyisobenzo-furan-1,3-dione
Appearance	White to off-white solid.
CAS	56-25-7
Couple Target	Serine/threonine-protein phosphatase
Couple Type	Inhibitor
Formula	$C_{10}H_{12}O_4$
MI	14: 1752
MW	196.2
Purity	≥98% (GC)
Solubility	Soluble in DMSO or 100% ethanol.

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-
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)