

Cyclopiazonic acid

ATPase inhibitor

Mycotoxin. Cell permeable. Induces the release of intracellular stored Ca^{2+} without increasing IP3 levels via inhibition of endoplasmic reticulum Ca^{2+} -ATPase. Highly specific inhibitor of Ca^{2+} -ATPase of sarcoplasmic reticulum, completely inhibiting the enzyme at 6-8 nmol/mg protein (at 0.5-2.0 μM ATP).

Citations: 20

[View Online »](#)

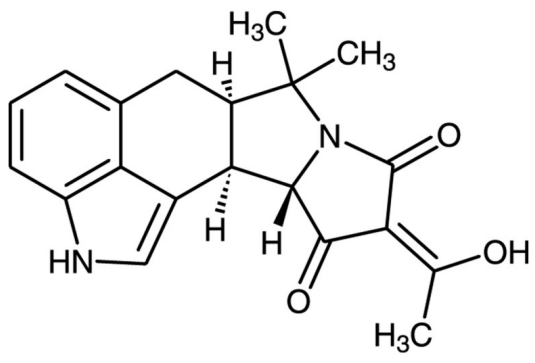
Ordering Information

[Order Online »](#)

BML-CA415-0010	10mg
----------------	------

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability As indicated on product label or CoA when stored as recommended.

Long Term Storage -20°C

Shipping Ambient Temperature

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name	CPA
Appearance	Off-white to faint yellow solid.
CAS	18172-33-3
Couple Target	ATPase
Couple Type	Inhibitor
Formula	$C_{20}H_{20}N_2O_3$
MW	336.4
Purity	≥98% (TLC)
RTECS	UY8587000
Solubility	Soluble in DMSO (20mg/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-
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-
de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

Germany
Phone: +49 7621 5500 526
[info-
uk@enzolifesciences.com](mailto:info-uk@enzolifesciences.com)

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