

Calmidazolium chloride

Calmodulin inhibitor

Binds reversibly to calmodulin, inhibiting calmodulin-mediated enzyme activation. It inhibits the Ca^{2+} /calmodulin-induced activation of myosin light chain kinase in a concentration dependent manner ($\text{IC}_{50}=1\text{ }\mu\text{M}$). Cell permeable. Inhibits calmodulin N-methyltransferase ($K_i=65\text{ nM}$). Blocks Ca^{2+} -calmodulin binding to NOS. Inhibitor of voltage-gated Ca^{2+} channels

Citations: 13

[View Online »](#)

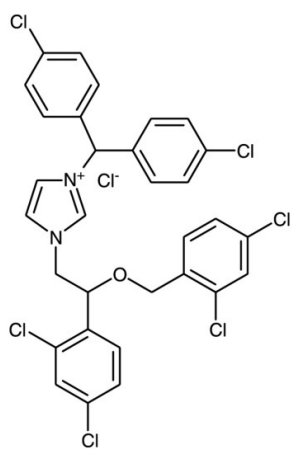
Ordering Information

[Order Online »](#)

BML-CA300-0010	10mg
BML-CA300-0050	50mg

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

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

Long Term Storage +4°C

Shipping Ambient Temperature

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name R24571

Appearance White powder.

CAS 57265-65-3

Couple Target Calcium channel, Calmodulin, CAM KMT

Couple Type Blocker, Inhibitor

Formula $C_{31}H_{23}Cl_7N_2O$

MW 687.7

Purity ≥98% (HPLC)

Solubility Soluble in 100% ethanol (>100mg/ml), DMSO (>100mg/ml), methanol, chloroform or propylene glycol. Almost insoluble in water.

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-
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)