

E₆ Berbamine

Calmodulin inhibitor

A potent, selective, cell-permeable calmodulin antagonist. It inhibits calmodulin-dependent enzymes such as MLC kinase ($IC_{50}=8\text{ }\mu\text{M}$) and phosphodiesterase I ($IC_{50}=0.53\text{ }\mu\text{M}$).

Citations: 2

[View Online »](#)

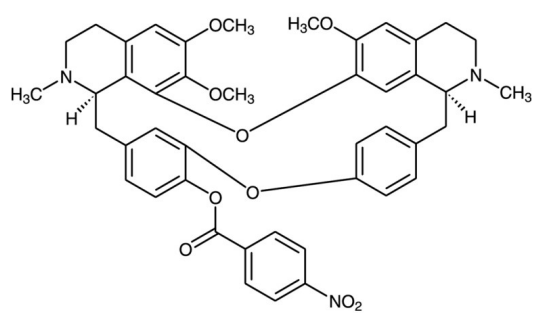
Ordering Information

[Order Online »](#)

BML-CA302-0050	50mg
----------------	------

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

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

Long Term Storage -20°C

Shipping Ambient Temperature

Regulatory Status RUO - Research Use Only

Product Details

Appearance Yellow solid.

CAS 114784-59-7

Couple Target Calmodulin, MLCK, PDE

Couple Type Inhibitor

Formula $C_{44}H_{43}N_3O_9$

MW 757.8

Purity ≥98% (TLC)

Solubility Soluble in DMSO (25 mg/ml) or ethanol (25 mg/ml).



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