

Deltamethrin

Calcineurin inhibitor

Synthetic type II pyrethroid. Potent calcineurin (protein phosphatase 2B) inhibitor $IC_{50} \sim 0.03$ nM. Causes an increase in neurotransmitter release at synapses and also increases intrasynaptosomal Ca^{2+} levels. May also have varied effects on ion channels.

Citations: 4

[View Online »](#)

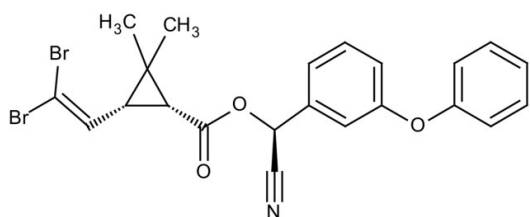
Ordering Information

[Order Online »](#)

BML-PR101-0010	10mg
BML-PR101-0050	50mg

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

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	(S)- α -Cyano-(3-phenoxybenzyl (1R)-cis-3-(2,2-dibromovinyl)-2,2-dimethylcyclopropanecarboxylate
Appearance	Beige solid.
CAS	52918-63-5
Couple Target	Serine/threonine-protein phosphatase
Couple Type	Inhibitor
Formula	$C_{22}H_{19}Br_2NO_3$
MI	14: 2883
MW	505.2
Purity	$\geq 98\%$ (External Standard Method)
Solubility	Soluble in DMSO (40mg/ml) or ethanol (15mg/ml, warm); 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-
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)