

Cypermethrin

Calcineurin inhibitor

Synthetic type II pyrethroid. Potent calcineurin (protein phosphatase 2B) inhibitor $IC_{50} \sim 0.03$ nM. May also have varied effects on ion channels. Serious doubts have been raised about the efficacy of this inhibitor.

Citations: 4

[View Online »](#)

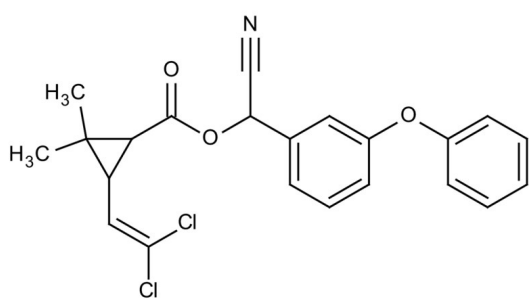
Ordering Information

[Order Online »](#)

BML-PR100-0050	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	Alphamethrine, Pestanal, (R,S)- α -Cyano-3-phenoxybenzyl(1RS)-cis,trans-3-(2,2-dichlorovinyl)-2,2-dimethylcyclopropane carboxylate
Appearance	White to off-white solid.
CAS	52315-07-8
Couple Target	Serine/threonine-protein phosphatase
Couple Type	Inhibitor
Formula	$C_{22}H_{19}Cl_2NO_3$
MI	14: 2767
MW	416.3
Purity	$\geq 96\%$ (HPLC) (mixture of isomers:racemic and cis/trans)
RTECS	GZ1250000
Solubility	Soluble in ethanol (25mg/ml). Insoluble in water.



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
info-usa@enzolifesciences.com

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
info-eu@enzolifesciences.com

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
info-fr@enzolifesciences.com

France
Phone: +33 472 440 655
info-fr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info-de@enzolifesciences.com

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