

Permethrin

Calcineurin inhibitor

Synthetic type I pyrethroid. Structurally similar to the protein phosphatase 2B (calcineurin) inhibitors cypermethrin and deltamethrin but weakly active or inactive depending on assay conditions. May be used as a negative control for cypermethrin (Prod. No. BML-PR100) and deltamethrin (BML-PR101).

Citations: 1

[View Online »](#)

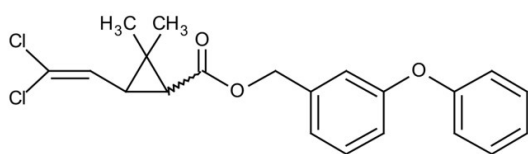
Ordering Information

[Order Online »](#)

BML-PR103-0050	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	3-Phenoxybenzyl (1R,S)-cis,trans-3-(2,2-dichlorovinyl)-2,2-dimethylcyclo-propanecarboxylate
Appearance	Beige solid.
CAS	52645-53-1
Couple Target	Serine/threonine-protein phosphatase
Couple Type	Inhibitor
Formula	$C_{21}H_{20}Cl_2O_3$
MI	14: 7180
MW	391.3
MeltingPoint	35-34°C
Purity	≥96% (HPLC)
RTECS	GZ1255000
Solubility	Soluble in DMSO (40mg/ml) or ethanol (25mg/ml). Insoluble in water.



Last modified: May 29, 2024

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

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

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

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

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

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