

# Cypermethrin

## Calcineurin inhibitor

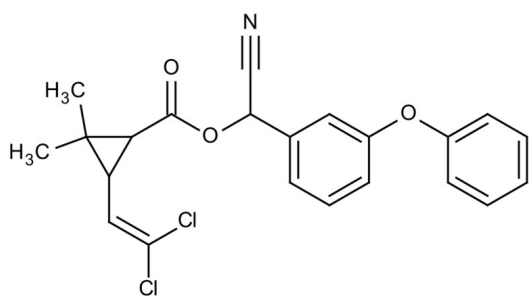
Synthetic type II pyrethroid. Potent calcineurin (protein phosphatase 2B) inhibitor  $IC_{50} \sim 0.03$  nM. May also have varied effects on ion channels. Serious doubts have been raised about the efficacy of this inhibitor.

Citations: 4

[View Online »](#)

Manuals, SDS & CofA

[View Online »](#)



## Handling & Storage

<b>Use/Stability</b>	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.
<b>Long Term Storage</b>	Ambient
<b>Shipping</b>	Ambient Temperature

**Regulatory Status** RUO - Research Use Only

## Product Details

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