

# bis(Methylthio)gliotoxin

## PAF receptor ligand

Inhibits PAF-induced platelet aggregation ( $IC_{50}=8.4\text{ }\mu\text{M}$ ) and collagen-induced aggregation ( $IC_{50}=84.2\text{ }\mu\text{M}$ ) with no effect on arachidonic acid or ADP induced aggregation. Inhibits PAF-induced bronchoconstriction *in vivo* in the guinea-pig (69% at 1.0 mg/kg).

Citations: 5

[View Online »](#)

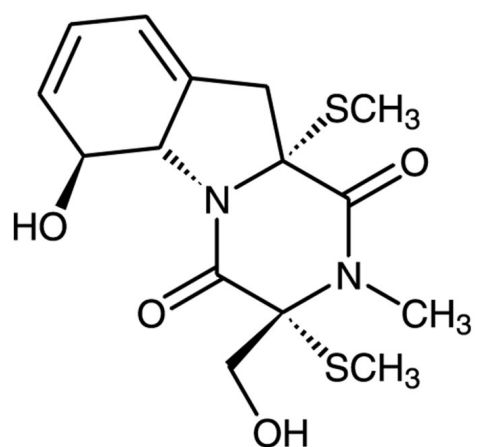
## Ordering Information

[Order Online »](#)

BML-L122-0001	1mg
BML-L122-0005	5mg

## Manuals, SDS & CofA

[View Online »](#)



## Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended. Store, as supplied, at -20°C for up to 1 year. Store solutions at -20°C for up to 3 months.
Long Term Storage	-20°C
Shipping	Ambient Temperature

**Regulatory Status** RUO - Research Use Only

## Product Details

Alternative Name	FR-49175
Appearance	White solid.
CAS	74149-38-5
Couple Target	PAF receptor
Couple Type	Ligand
Formula	$C_{15}H_{20}N_2O_4S_2$
MW	356.5
Purity	≥98% (TLC)
Solubility	Soluble in DMSO (25mg/ml) and ethanol (25mg/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-  
fr@enzolifesciences.com](mailto:info-fr@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)