

# Manoalide

## Phospholipase inhibitor

Marine sesterterpenoid. Irreversibly inhibits human synovial phospholipase A<sub>2</sub> (PLA<sub>2</sub>, IC<sub>50</sub>=0.02-0.2μM depending on substrate) as well as bee (IC<sub>50</sub>=0.05μM) and cobra venom (IC<sub>50</sub>=2μM) PLA<sub>2</sub> by covalently modifying lysine residues. Inhibits human recombinant sPLA<sub>2</sub> (IC<sub>50</sub>=4.7μM) selectively over cPLA<sub>2</sub>. Also inhibits phospholipase C (PLC, IC<sub>50</sub>=1.5μM) and Ca<sup>2+</sup> channels. Inhibits superoxide generation in polymorphonuclear leukocytes. Inhibits the release of arachidonate in cultured mouse peritoneal macrophages and in a mouse *in vivo* model. Has anti-inflammatory and analgesic properties. Shows strong quorum sensing antagonism.

Citations: 14

[View Online »](#)

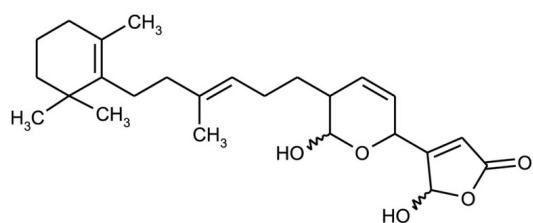
## Ordering Information

[Order Online »](#)

BML-EI177-0001	1mg
----------------	-----

## 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 at -20°C. Stock solutions are stable for up to 3 months at -20°C.
<b>Long Term Storage</b>	-20°C
<b>Shipping</b>	Ambient Temperature

## Regulatory Status

RUO - Research Use Only

## Product Details

<b>Appearance</b>	Colorless to pale waxy or filmy residue.
<b>CAS</b>	75088-80-1
<b>Couple Target</b>	Calcium channel, Phospholipase, PLA, PLC
<b>Couple Type</b>	Inhibitor
<b>Formula</b>	$C_{25}H_{36}O_5$
<b>MW</b>	416.6
<b>Purity</b>	≥98% (TLC)
<b>Solubility</b>	Soluble in DMSO (25mg/ml), 100% ethanol (25mg/ml) or methanol.



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)