

# (±)-Lisofylline

Potent inhibitor of phosphatidic acid generation

A potent inhibitor of phosphatidic acid generation ( $IC_{50}=0.6 \mu M$ ). It protects mice from endotoxic shock and attenuates sepsis-induced acute lung injury in the pig. Anti-inflammatory agent. Enhances glucose-stimulated insulin secretion. Inhibits IL-12-mediated STAT4 activation.

Citations: 1

[View Online »](#)

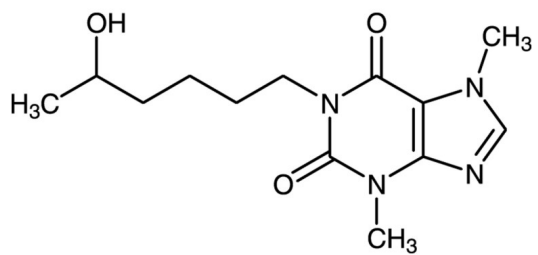
Ordering Information

[Order Online »](#)

BML-LP102-0010	10mg
----------------	------

Manuals, SDS & CofA

[View Online »](#)



## Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended.
Long Term Storage	Ambient
Shipping	Ambient Temperature

**Regulatory Status** RUO - Research Use Only

## Product Details

Appearance	White to off-white solid.
CAS	6493-06-7
Formula	$C_{13}H_{20}N_4O_3$
MW	280.3
Purity	≥98% (HPLC)

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)