

# N-Palmitoyl phosphatidylethanolamine

Reduces food intake

Endogenous lipid which significantly reduced food intake in a dose-dependent fashion when administered to rats and mice (100 mg/kg, i.p.), without causing taste aversion. Has been found to enter the brain (concentrating in the hypothalamus) when administered systemically and also reduced food intake when infused intracerebroventricularly (nanomolar concentrations). This suggests that mediation of effects is *via* the CNS. Chronic infusion resulted in a reduction of both food intake and body weight.

Citations: 1

[View Online »](#)

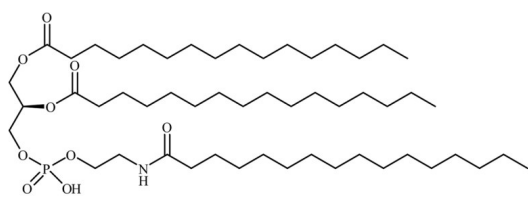
## Ordering Information

[Order Online »](#)

BML-LP108-0001	1mg
BML-LP108-0005	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. MAKE SOLUTIONS FRESH DAILY.
Long Term Storage	-20°C
Shipping	Blue Ice

## Regulatory Status

RUO - Research Use Only

## Product Details

Alternative Name	C16:0 NAPE, NPPE
Appearance	White solid.
CAS	108321-22-8
Formula	$C_{53}H_{104}NO_9P$
MW	930.4
Purity	≥98% (TLC)
Solubility	Soluble in 100% ethanol (2 mg/ml with ~30 minutes sonication at 37°C) or dimethylacetamide (1 mg/ml with ~30 minutes sonication at 37°C).

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)