

# 1-O-Hexadecyl-2-O-docosa-hexaenoyl-sn-glycero-3-phosphorylcholine

**Ultra-pure** A biosynthetic precursor of PAF *via* the lipid remodeling pathway.

Citations: 1

[View Online »](#)

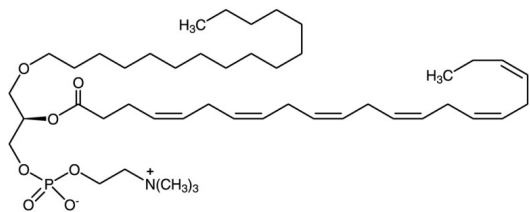
Ordering Information

[Order Online »](#)

BML-L117-0005	5mg
---------------	-----

Manuals, SDS & CofA

[View Online »](#)



## Handling & Storage

<b>Use/Stability</b>	As indicated on product label or CoA when stored as recommended. Store, as supplied, at -80°C for up to 1 year. Subject to oxidation – Protect from air. Open ampule and aliquot into suitable vials. Store at -20°C or colder. For experiments with cells or tissues, aliquots of ethanol solutions may be added directly to aqueous media.
<b>Long Term Storage</b>	-80°C
<b>Shipping</b>	Dry Ice

**Regulatory Status** RUO - Research Use Only

## Product Details

<b>Appearance</b>	Pale yellow solid dissolved in liquid.
<b>CAS</b>	132213-85-5
<b>Formula</b>	$C_{46}H_{86}NO_7P$
<b>Formulation</b>	Supplied as solid dissolved in ethanol.
<b>MW</b>	792.2
<b>Purity</b>	≥98%

**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-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)