

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

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

Ultra-pure

Citations: 1

[View Online »](#)

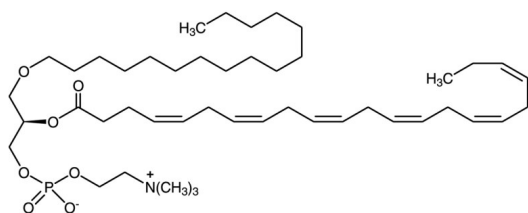
Ordering Information

[Order Online »](#)

BML-L117-0005	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 -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.
Long Term Storage	-80°C
Shipping	Dry Ice

Regulatory Status

RUO - Research Use Only

Product Details

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



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)