

1-O-Hexadecyl-2-O-acetyl-sn-glycero-3-phosphorylcholine

PAF receptor ligand

Platelet activating factor (PAF) with a C₁₆ alkyl group. Initiates receptor-mediated responses in platelets and a variety of other cells. Rabbit platelet PAF receptor K_D=0.1 nM. EC₅₀ for stimulation of phosphoinositide metabolism in hippocampal cell line: 1.2-1.5 nM.

Citations: 3

[View Online »](#)

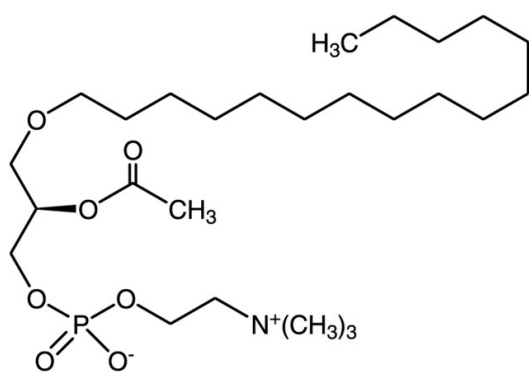
Ordering Information

[Order Online »](#)

BML-L100-0005	5mg
---------------	-----

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended. Stable for at least 1 year after receipt when stored, as supplied, at -20°C. Stock solutions are stable for up to 2 months at -20°C.
Long Term Storage	-20°C
Shipping	Ambient Temperature

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	PAF C-16
Appearance	White solid.
CAS	74389-68-7
Couple Target	PAF receptor
Couple Type	Ligand
Formula	$C_{26}H_{54}NO_7P$
MI	14: 7524
MW	523.7
Purity	≥98% (TLC)
Solubility	Soluble in 100% ethanol (>100 mg/ml). Soluble in water (>100mg/ml) after sonication for 1 minute at 40°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-
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)