

Carbamyl-PAF

PAF receptor activator

Biologically potent, non-metabolizable analog of platelet-activating factor (PAF). Induces half-maximal PMNL degranulation response at 16nM and binds stereospecifically to high and low affinity PAF receptors. Completely resists inactivation by acetylhydrolase.

Citations: 4

[View Online »](#)

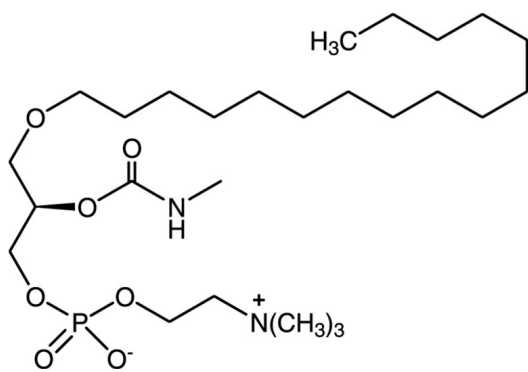
Ordering Information

[Order Online »](#)

BML-L120-0005	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. Ethanol stock solutions are stable for up to 3 months at -20°C. Aqueous solutions should be prepared fresh daily.
Long Term Storage	-20°C
Shipping	Ambient Temperature

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	1-O-Hexadecyl-2-N-methylcarbamyl-sn-glycero-3-phosphocholine
Appearance	White solid.
CAS	91575-58-5
Couple Target	PAF receptor
Couple Type	Activator, Ligand
Formula	$C_{26}H_{55}N_2O_7P$
MW	538.7
Purity	≥98% (TLC)
Solubility	Soluble in 100% ethanol.

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)