

Phosphatidic acid, dipalmitoyl

PKC activator

A mitogenic lipid phosphate produced *via* the action of phospholipase D. Mediates agonist-dependent Raf-1 translocation and activation of the MAP kinase cascade.

Citations: 4

[View Online »](#)

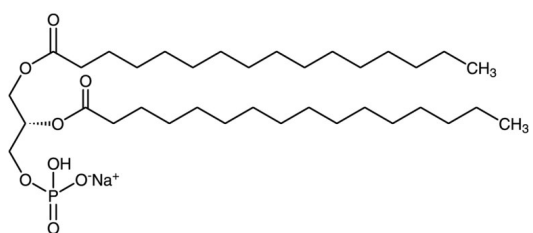
Ordering Information

[Order Online »](#)

BML-LP103-0100	100mg
----------------	-------

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 3 months at -20°C.
Long Term Storage	-20°C
Shipping	Ambient Temperature

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	1,2-Dipalmitoyl-sn-glycero-3-phosphate . sodium salt, DPPA
Appearance	White solid.
CAS	169051-60-9
Couple Target	PKC
Couple Type	Activator
Formula	$C_{35}H_{68}O_8P \cdot Na$
MW	670.9
Purity	≥98% (TLC)
Solubility	Fine suspension in ethanol (5mg/ml) with 15 minutes of sonication at room temperature.



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)