

Phosphatidic acid, dioleoyl

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. Activates PKC ζ .

Citations: 4

[View Online »](#)

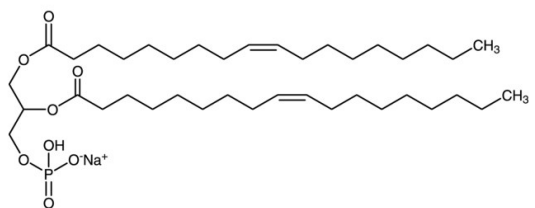
Ordering Information

[Order Online »](#)

BML-LP104-0010	10mg
----------------	------

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability As indicated on product label or CoA when stored as recommended.

Long Term Storage -20°C

Shipping Ambient Temperature

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name 1,2-Dioleoyl-sn-glycero-3-phosphate . Na

Appearance White powder.

CAS 108392-02-5

Couple Target PKC

Couple Type Activator

Formula $C_{39}H_{73}O_8P \cdot Na$

Identity Determined by IR.

MW 724

Purity ≥98% (GC)

Solubility Soluble in chloroform.

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)