

Lysophosphatidic acid

LPA receptor ligand

Lysophosphatidic acid (LPA) is a multifunctional intercellular phospholipid messenger. LPA stimulates the growth of a variety of cells including fibroblasts, vascular smooth muscle cells, endothelial cells and keratinocytes. It acts as a proliferative and anti-apoptotic factor. Ligand for LPA1 (EDG-2), LPA2 (EDG-4) and LPA3 (EDG-7) receptors. It also inhibits differentiation of neural stem cells into neurons.

Citations: 12

[View Online »](#)

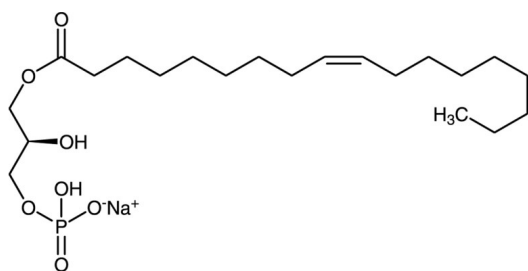
Ordering Information

[Order Online »](#)

BML-LP100-0005	5mg
BML-LP100-0025	25mg

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended. Store as supplied at -20°C for up to 1 year. Store solutions at -20°C for up to 1 month.
Long Term Storage	-20°C
Shipping	Ambient Temperature

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	1-Oleoyl-lysophosphatidic acid . sodium salt
Appearance	Waxy solid.
CAS	325465-93-8
Couple Target	LPA receptor
Couple Type	Ligand
Formula	$C_{21}H_{40}O_7P \cdot Na$
MW	458.5
Purity	≥98% (TLC)
Solubility	Soluble in water (5 mg/mL with sonication) or DMSO (0.5 mg/mL warm).

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)