

L-NASPA

LPA receptor ligand

L-NASPA is an antagonist of LPA-induced platelet aggregation and chloride conductance in *Xenopus* oocytes. In human breast carcinoma MDA-MB231 cells it is a potent LPA agonist.

Citations: 3

[View Online »](#)

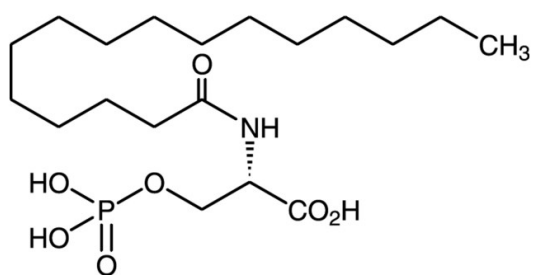
Ordering Information

[Order Online »](#)

BML-LP101-0001	1mg
----------------	-----

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended. Store, as supplied, at -80°C for up to 1 year. Store solutions at -20°C for up to 1 year.
Long Term Storage	-80°C
Shipping	Dry Ice

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	N-Palmitoyl-L-serine phosphoric acid
Appearance	White to off-white solid.
CAS	155915-46-1
Couple Target	LPA receptor
Couple Type	Ligand
Formula	$C_{19}H_{38}NO_7P \cdot H_2O$
MW	440.5
Purity	≥98% (TLC)
Solubility	Soluble in DMSO (10 mg/ml) or ethanol (10 mg/ml).

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)