

L-AP3

mGluR antagonist

Selective antagonist of the PI-linked metabotropic glutamate response.

Citations: 5

[View Online »](#)

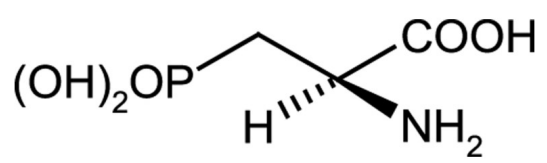
Ordering Information

[Order Online »](#)

ALX-550-038-M005	5mg
------------------	-----

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended.
Handling	Protect from light.
Long Term Storage	+4°C
Shipping	Ambient Temperature

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name	L(+)-2-Amino-3-phosphonopropanoic acid
Appearance	White hygroscopic solid.
CAS	23052-80-4
Couple Target	Metabotropic glutamate receptor
Couple Type	Inhibitor, Ligand
Formula	$C_3H_8NO_5P$
MW	169.1
Purity	≥97%
Solubility	Soluble in water.

Last modified: May 29, 2024



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
info-usa@enzolifesciences.com

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
info-eu@enzolifesciences.com

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
info-be@enzolifesciences.com

France
Phone: +33 472 440 655
info-fr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info-de@enzolifesciences.com

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
info-uk@enzolifesciences.com