

L-AP4

mGluR agonist

Selective agonist for group III mGluRs. NMDA agonist. Potent synaptic depressant.

Citations: 9

[View Online »](#)

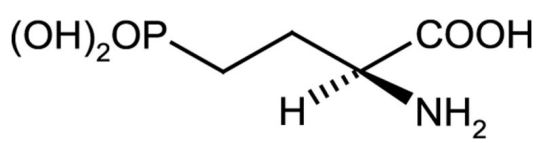
Ordering Information

[Order Online »](#)

ALX-550-026-M025	25mg
------------------	------

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	Ambient
Shipping	Ambient Temperature

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	L(+)-2-Amino-4-phosphonobutanoic acid
Appearance	White crystalline powder.
CAS	23052-81-5
Couple Target	Metabotropic glutamate receptor, NMDA receptor
Couple Type	Activator, Ligand
Formula	$C_4H_{10}NO_5P$
MW	183.1
Purity	≥97% (HPLC)
Solubility	Soluble in dilute aqueous base.

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-be@enzolifesciences.com)

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
[info-
be@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)