

DL-AP5

NMDA receptor antagonist

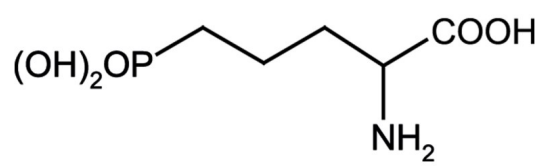
Potent and selective NMDA receptor antagonist.

Citations: 5

[View Online »](#)

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	DL-2-Amino-5-phosphonopentanoic acid, DL-2-Amino-5-phosphonovaleric acid
Appearance	White to off-white powder.
CAS	76326-31-3
Couple Target	NMDA receptor
Couple Type	Inhibitor, Ligand
Formula	$C_5H_{12}NO_5P$
MW	197.1
Purity	≥99% (HPLC, TLC)
Solubility	Soluble in 100mM NaOH or 25mM 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