

# 2-(Phosphonomethyl)-pentanedioic acid

NAALADase inhibitor

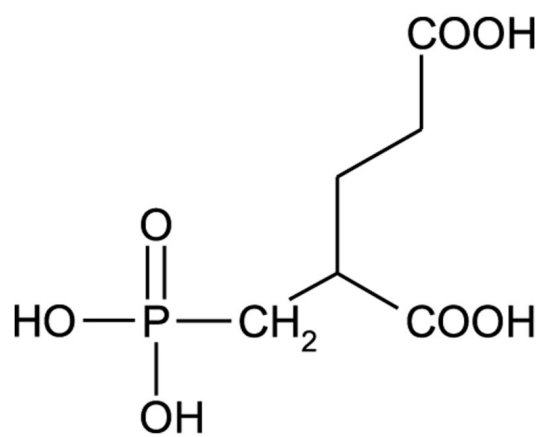
Potent inhibitor of the neuropeptidase N-acetylated  $\alpha$ -linked acidic dipeptidase (NAALADase), which is a membrane-bound peptidase that hydrolyses N-acetyl-L-aspartate-L-glutamate (NAAG) (Prod. No. ALX-151-020), a major brain peptide.

Citations: 3

[View Online »](#)

Manuals, SDS & CofA

[View Online »](#)



## Handling & Storage

**Use/Stability** As indicated on product label or CoA when stored as recommended.

**Long Term Storage** -20°C

**Shipping** Ambient Temperature

**Regulatory Status** RUO - Research Use Only

## Product Details

<b>Alternative Name</b>	PMPA
<b>Appearance</b>	Viscous, colorless oil.
<b>CAS</b>	173039-10-6
<b>Couple Target</b>	NAALADase
<b>Couple Type</b>	Inhibitor
<b>Formula</b>	$C_6H_{11}O_7P$
<b>Identity</b>	Determined by NMR.
<b>MW</b>	226.1
<b>Purity</b>	≥97% (HPLC)
<b>Solubility</b>	Soluble in water, methanol or DMSO.

**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-  
be@enzolifesciences.com](mailto:info-be@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)