

(R)-CPP

NMDA receptor antagonist

Highly potent NMDA receptor antagonist.

Citations: 3

[View Online »](#)

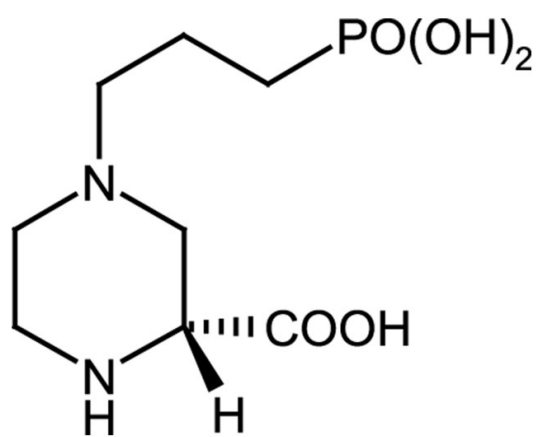
Ordering Information

[Order Online »](#)

ALX-550-231-M001	1mg
------------------	-----

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

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

Long Term Storage Ambient

Shipping Ambient Temperature

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name 3-((R)-2-Carboxypiperazin-4-yl)-propyl-1-phosphonic acid

Appearance White crystalline powder.

CAS 126453-07-4

Couple Target NMDA receptor

Couple Type Inhibitor, Ligand

Formula $C_8H_{17}N_2O_5P$

MW 252.2

Purity $\geq 97\%$ (HPLC)

Solubility Soluble in dilute aqueous base or water.

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)