

N-Methyl-D-aspartic acid

NMDA receptor agonist

Selective ionotropic NMDA receptor agonist. Excitatory amino acid.

Citations: 2

[View Online »](#)

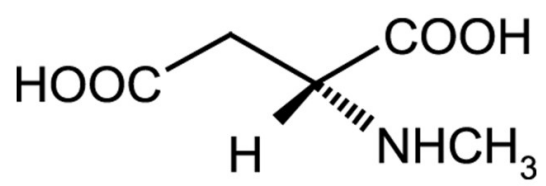
Ordering Information

[Order Online »](#)

ALX-105-009-M050	50mg
------------------	------

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 NMDA

Appearance White crystalline powder.

CAS 6384-92-5

Couple Target NMDA receptor

Couple Type Activator

Formula $C_5H_9NO_4$

Identity Determined by NMR.

MI 14: 6662

MW 147.1

Purity $\geq 98\%$ (TLC)

RTECS CI9457000

Solubility Soluble in dilute aqueous base; slightly soluble in water or 100% ethanol.



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