

N-Acetyl-Asp-Glu-OH

Glutamate receptor ligand

Endogenous neurotransmitter with high affinity for the brain glutamate receptor. N-acetylated α -linked acidic dipeptidase (NAALADase), a membrane-bound peptidase, hydrolyzes α -NAAG, a major brain peptide, to N-acetylaspartate and glutamate. Discriminates between mGluR2 and mGluR3. mGluR3 activator. NMDA receptor antagonist. Neuroprotective.

Citations: 13

[View Online »](#)

Ordering Information

[Order Online »](#)

| | |
|------------------|------|
| ALX-151-020-M010 | 10mg |
| ALX-151-020-M050 | 50mg |

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

| | |
|-------------------|--|
| Use/Stability | As indicated on product label or CoA when stored as recommended. |
| Long Term Storage | +4°C |
| Shipping | Ambient Temperature |

Regulatory Status

RUO - Research Use Only

Product Details

| | |
|------------------|--|
| Alternative Name | α -NAAG, Spaglumatic acid |
| Appearance | White to off-white powder. |
| CAS | 3106-85-2 |
| Formula | $C_{11}H_{16}N_2O_8$ |
| MW | 304.3 |
| Purity | ≥98% (TLC) |
| Solubility | Soluble in water and methanol (50mg/ml). |

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