

(2S,4R)-4-Methylglutamic acid

Kainate receptor agonist

Potent and highly selective kainate receptor agonist, with an IC_{50} for inhibition of [3H]-kainate binding of 35 nM and almost 3,000- and 200-fold selectivity for kainate receptors over AMPA and NMDA receptors respectively. Also selectively inhibits the cloned excitatory amino acid transporter EAAT2 at higher concentrations.

Citations: 4

[View Online »](#)

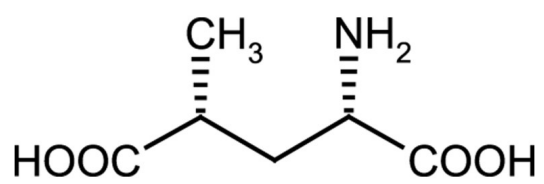
Ordering Information

[Order Online »](#)

ALX-550-186-M005	5mg
------------------	-----

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

Alternative Name SYM 2081

Appearance White powder.

CAS 31137-74-3

Couple Target Kainate receptor

Couple Type Activator

Formula $C_6H_{11}NO_4$

MW 161.2

Purity ≥97%

Solubility Soluble in dilute aqueous acid.



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)