

Fenobam

mGluR5 antagonist

Potent and selective non-competitive mGluR5 antagonist that displays inverse agonist properties; blocks mGluR5 constitutive activity *in vitro* (IC_{50} =87nM). Acts at an allosteric modulatory site shared with MPEP (Prod. No. ALX-550-368) and binds mGluR5 with K_d values of 54nM and 31nM for rat and human receptors respectively. Displays anxiolytic activity following oral administration *in vivo*.

Citations: 4

[View Online »](#)

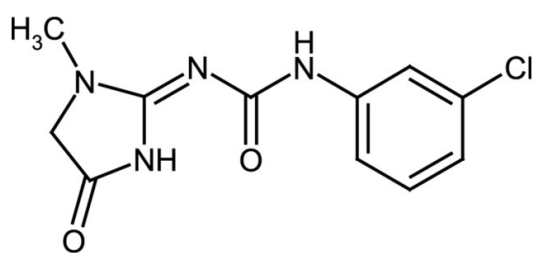
Ordering Information

[Order Online »](#)

ALX-550-413-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 +4°C

Shipping Ambient Temperature

Regulatory Status RUO - Research Use Only

Product Details

Appearance Off-white to pale orange solid.

CAS 57653-26-6

Couple Target Metabotropic glutamate receptor

Couple Type Inhibitor

Formula $C_{11}H_{11}ClN_4O_2$

Identity Determined by NMR.

MW 266.7

Purity ≥98% (1H-NMR)

Solubility Soluble in DMSO or methanol.

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)