

# MTEP

## mGluR5 antagonist

Potent and highly selective antagonist for mGluR5 with a 5-fold higher anxiolytic activity than MPEP (Prod. No. ALX-550-368). Shows fewer off-target effects than MPEP.

Citations: 8

[View Online »](#)

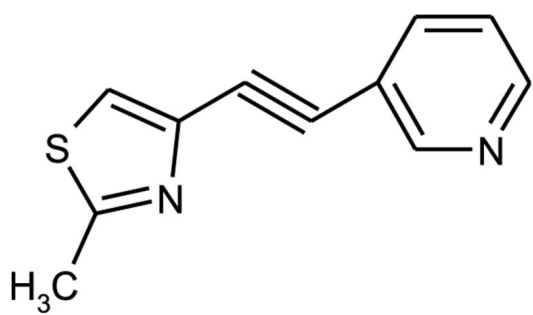
### Ordering Information

[Order Online »](#)

ALX-550-381-M005	5mg
------------------	-----

### 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	+4°C
Shipping	Ambient Temperature

## Regulatory Status

RUO - Research Use Only

## Product Details

Appearance	White to yellow solid.
CAS	329205-68-7
Couple Target	Metabotropic glutamate receptor
Couple Type	Inhibitor, Ligand
Formula	$C_{11}H_8N_2S$
MW	200.3
Purity	≥98% (HPLC)
Solubility	Soluble in DMSO, 100% ethanol 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)