

MPEP . hydrochloride

mGluR5 antagonist

Potent and selective antagonist for mGluR5 receptor. Systemically active *in vivo*.

Citations: 13

[View Online »](#)

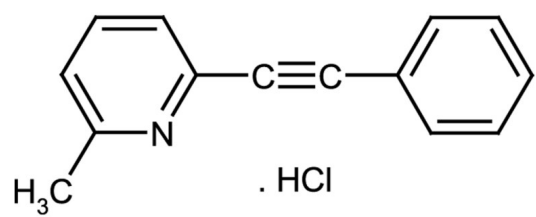
Ordering Information

[Order Online »](#)

ALX-550-368-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

Alternative Name 6-Methyl-2-(phenylethynyl)-pyridine . HCl

Appearance Off-white powder.

CAS 96206-92-7

Couple Target Metabotropic glutamate receptor

Couple Type Inhibitor, Ligand

Formula $C_{14}H_{11}N \cdot HCl$

Identity Determined by NMR.

MW 193.3 . 36.5

Purity ≥99%

Solubility Soluble in water, DMSO or methanol.

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