

AIDA

mGluR agonist

Potent and selective ligand for group I mGluRs. Shows no effect on group II or group III mGluRs. Has no effect on ionotropic glutamate receptors.

Citations: 9

[View Online »](#)

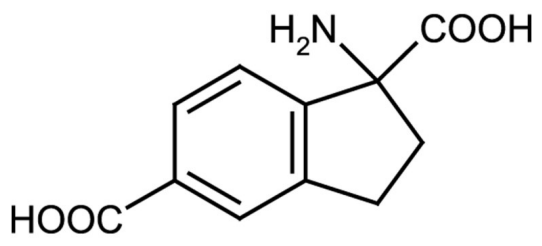
Ordering Information

[Order Online »](#)

ALX-550-101-M025	25mg
------------------	------

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	(±)-1-Aminoindane-1,5-dicarboxylic acid, UPF 523
CAS	168560-79-0
Couple Target	Metabotropic glutamate receptor
Couple Type	Activator, Ligand
Formula	$C_{11}H_{11}NO_4$
MW	221.2
Purity	≥96%
Solubility	Soluble in water.



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