

(2R,4R)-APDC

mGluR agonist

Selective agonist for group II mGluRs. Inhibits the NADPH-oxidase dependent ROS system.

Citations: 14

[View Online »](#)

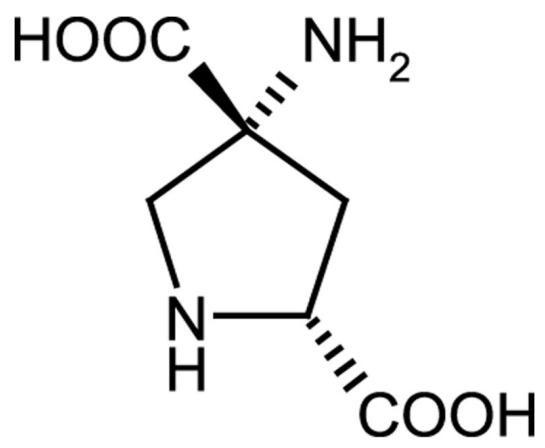
Ordering Information

[Order Online »](#)

ALX-550-178-M001	1mg
ALX-550-178-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 Ambient

Shipping Ambient Temperature

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name LY-314,593, (2R,4R)-4-Aminopyrrolidine-2,4-dicarboxylic acid

Appearance Off-white powder.

CAS 169209-63-6

Couple Target Metabotropic glutamate receptor

Couple Type Activator, Ligand

Formula $C_6H_{10}N_2O_4$

MW 174.2

Purity ≥98% (HPLC)

Solubility Soluble in water.



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-
de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

Germany
Phone: +49 7621 5500 526
[info-
uk@enzolifesciences.com](mailto:info-uk@enzolifesciences.com)

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
[info-
uk@enzolifesciences.com](mailto:info-uk@enzolifesciences.com)