

Cyclothiazide

AMPA receptor activator

Potently inhibits the desensitization of AMPA receptors. Antihypertensive.

Citations: 8

[View Online »](#)

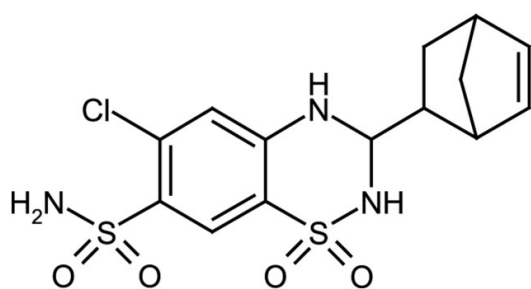
Ordering Information

[Order Online »](#)

ALX-550-338-M050	50mg
------------------	------

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 3-Bicyclo[2.2.1]hept-5-en-2-yl-6-chloro-3,4-dihydro-2H-1,2,4-benzothiadiazine-7-sulfonamide 1,1-dioxide, 6-Chloro-3,4-dihydro-3[2-norbornen-5-yl]-2H-1,2,4-benzothiadiazine-7-sulfonamide 1,1-dioxide

Appearance White to off-white solid.

CAS 2259-96-3

Couple Target AMPA receptor

Couple Type Activator, Ligand

Formula $C_{14}H_{16}ClN_3O_4S_2$

MI 14: 2754

MW 389.9

Purity ≥98% (HPLC (UV))

Solubility Soluble in DMSO, methanol (36mg/ml) or dilute aqueous base (36mg/ml); insoluble in water.

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