

5-Nitro-6,7-dichloro- 1,4-dihydro-2,3- quinoxalinedione

NMDA receptor antagonist

Potent and selective NMDA receptor glycine site antagonist.

Citations: 3

[View Online »](#)

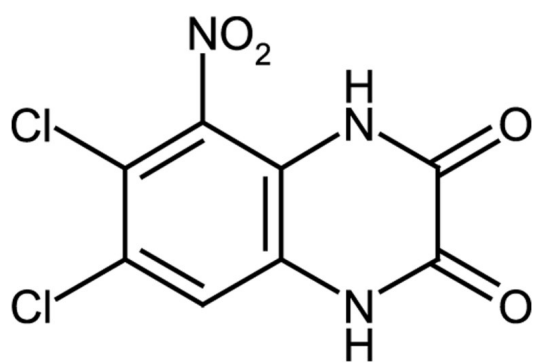
Ordering Information

[Order Online »](#)

ALX-550-228-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 ACEA-1021

Appearance Yellow powder.

CAS 153504-81-5

Couple Target NMDA receptor

Couple Type Inhibitor

Formula $C_8H_3Cl_2N_3O_4$

MW 276.0

Purity ≥98% (TLC; elemental analysis)

Solubility Soluble in DMSO or dilute aqueous base.

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)