

CNQX . disodium salt

AMPA/kainate receptor antagonist

Water soluble, potent competitive kainate/quisqualate (non-NMDA) receptor antagonist. More soluble form of CNQX (Prod. No. ALX-550-042).

Citations: 2

[View Online »](#)

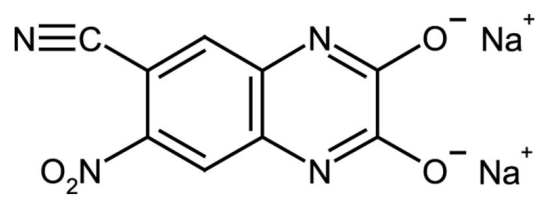
Ordering Information

[Order Online »](#)

ALX-550-283-M010	10mg
ALX-550-283-M025	25mg

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended.
Handling	Protect from light.
Long Term Storage	-20°C
Shipping	Ambient Temperature

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name	6-Cyano-7-nitroquinoxaline-2,3-dione . 2Na
Appearance	Rust red solid.
CAS	115066-14-3 (non-salt)
Couple Target	AMPA receptor, Kainate receptor
Couple Type	Inhibitor, Ligand
Formula	$C_9H_2N_4O_4 \cdot 2Na$
MW	230.1 . 46.0
Purity	≥98% (TLC)
Solubility	Soluble in water (5mg/ml).

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