

# CNQX

## AMPA/kainate receptor antagonist

Potent, competitive kainate/quisqualate (non-NMDA) receptor antagonist. For the more soluble form see CNQX . disodium salt (Prod. No. ALX-550-283).

Citations: 4

[View Online »](#)

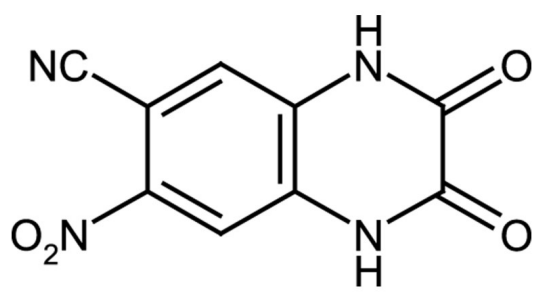
### Ordering Information

[Order Online »](#)

ALX-550-042-M005	5mg
ALX-550-042-M010	10mg
ALX-550-042-M050	50mg

### Manuals, SDS & CofA

[View Online »](#)



## Handling & Storage

<b>Use/Stability</b>	As indicated on product label or CoA when stored as recommended. Stock solutions are stable for several months when stored at -20°C.
<b>Handling</b>	Protect from light.
<b>Long Term Storage</b>	Ambient
<b>Shipping</b>	Ambient Temperature

## Regulatory Status

RUO - Research Use Only

## Product Details

<b>Alternative Name</b>	6-Cyano-7-nitroquinoxaline-2,3-dione, FG-9065
<b>Appearance</b>	Off-white to yellow powder.
<b>CAS</b>	115066-14-3
<b>Couple Target</b>	AMPA receptor, Kainate receptor
<b>Couple Type</b>	Inhibitor, Ligand
<b>Formula</b>	$C_9H_4N_4O_4$
<b>MW</b>	232.2
<b>Purity</b>	≥98% (TLC)
<b>Solubility</b>	Soluble in DMSO (10mg/ml) 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)