

# NBQX . disodium salt

AMPA/kainate receptor antagonist

Water soluble, potent and selective AMPA/kainate receptor antagonist. For literature references, see NBQX (Prod. No. ALX-550-044).

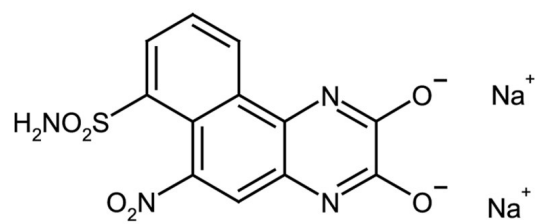
## Ordering Information

[Order Online »](#)

ALX-550-272-M005	5mg
ALX-550-272-46190	5mg
ALX-550-272-M025	25mg

## Manuals, SDS & CofA

[View Online »](#)



## Handling & Storage

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

**Regulatory Status** RUO - Research Use Only

## Product Details

<b>Alternative Name</b>	2,3-Dihydroxy-6-nitro-7-sulphamoylbenzo(f)-quinoxaline . 2Na, 6-Nitro-7-sulphamoylbenzo(f)-quinoxaline-2,3-dione . 2Na
<b>Appearance</b>	Red to brown solid.
<b>CAS</b>	118876-58-7 (non-salt)
<b>Couple Target</b>	AMPA receptor, Kainate receptor
<b>Couple Type</b>	Inhibitor, Ligand
<b>Formula</b>	$C_{12}H_6N_4O_6S \cdot 2Na$
<b>MW</b>	334.3 . 46.0
<b>Purity</b>	≥98% (HPLC)
<b>Solubility</b>	Soluble in water (5mg/ml) or DMSO.

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-  
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)