

NBQX

AMPA/kainate receptor antagonist

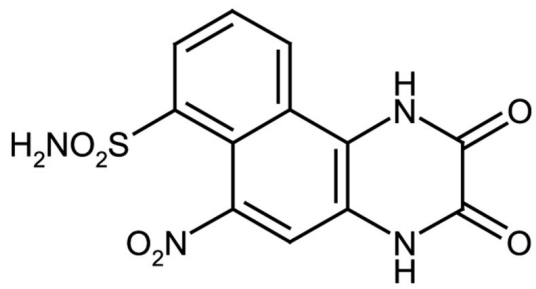
Most potent and selective AMPA/kainate receptor antagonist. For the more soluble form see NBQX . disodium salt (Prod. No. ALX-550-272).

Citations: 10

[View Online »](#)

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	+4°C
Shipping	Ambient Temperature

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	2,3-Dihydroxy-6-nitro-7-sulphamoylbenzo(f)-quinoxaline, 6-Nitro-7-sulphamoylbenzo(f)-quinoxaline-2,3-dione
Appearance	Yellow crystalline powder.
CAS	118876-58-7
Couple Target	AMPA receptor, Kainate receptor
Couple Type	Inhibitor, Ligand
Formula	$C_{12}H_8N_4O_6S$
MW	336.3
Purity	≥99% (HPLC)
Solubility	Soluble in DMSO. A 1mM aqueous solution may be prepared by taking up in >2.5 eq. of NaOH and backtitrating to pH 8.0-8.5 with conc. HCl.

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)