

# Nitric oxide synthase (neuronal) polyclonal antibody

This antibody is covered by our [Worry-Free Guarantee](#).

Citations: 2

[View Online »](#)

Ordering Information

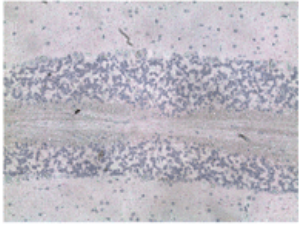
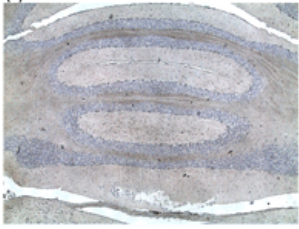
[Order Online »](#)

ALX-210-5290	100µg
--------------	-------

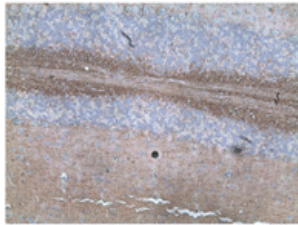
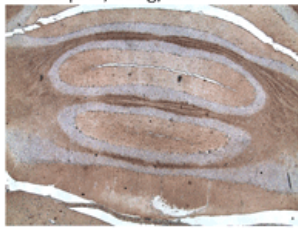
Manuals, SDS & CofA

[View Online »](#)

(-) Control



nNOS pAb, 50ug/mL



Immunohistochemistry analysis of mouse and rat brain tissue (paraffin sections) with Nitric oxide synthase (neuronal) (human), pAb (prod. no. ALX-210-529).

## Handling & Storage

**Handling** Avoid freeze/thaw cycles.

**Long Term Storage** -20°C

**Shipping** Blue Ice

**Regulatory Status** RUO - Research Use Only

## Product Details

**Alternative Name** nNOS, NOS I

**Application** ICC, IF, IHC, WB

**Application Notes** Detects a band of ~140kDa by Western blot.

**Formulation** Liquid. In PBS containing 1mg/ml BSA and 0.05% sodium azide.

**Host** Rabbit

**Immunogen** Synthetic peptide corresponding to aa 1411-1425 (CN1411RLRSESI AFIEESK1425) of human neuronal nitric oxide synthase (nNOS; NOS I). This sequence is completely conserved in human, mouse, rat, rabbit and frog nNOS.

**Purity Detail** Epitope affinity-purified.

**Recommendation Dilutions/Conditions** Immunocytochemistry (1:50-1:500)Immunofluorescence (1:50-1:500)Immunohistochemistry (1:20-1:200)Western blot (1:100-1:1000)Suggested dilutions/conditions may not be available for all applications.Optimal conditions must be determined individually for each application.

**Species Reactivity** Human, Mouse, Rat

**UniProt ID** P29475

**Worry-free Guarantee** This antibody is covered by our [Worry-Free Guarantee](#)

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)