

Nitric oxide synthase (neuronal) polyclonal antibody

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

Citations: 2

[View Online »](#)

Ordering Information

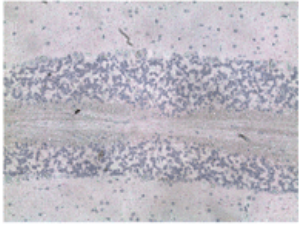
[Order Online »](#)

ALX-210-529-C100	100µg
------------------	-------

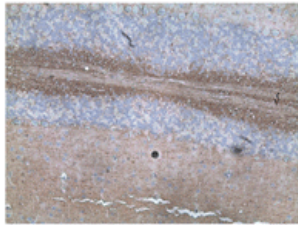
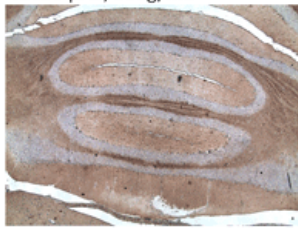
Manuals, SDS & CofA

[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)