Nitric oxide synthase (neuronal) polyclonal antibody

This antibody is covered by our Worry-Free Guarantee.

Citations: 1

View Online »

Ordering Information

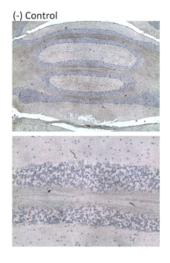
Order Online »

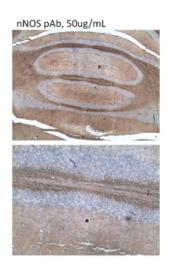
ALX-210-529-C100

100µg

Manuals, SDS & CofA

View Online »





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 NotesDetects 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

(CN1411RLRSESIAFIEESK1425) 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



info-

eu@enzolifesciences.com