nNOS polyclonal antibody

The diffusible free radical gas nitric oxide (NO) affects a variety of physiological functions, and is a key regulator of the cardiovascular, nervous, and immune systems. NO is synthesized in many tissues from Larginine, oxygen, and NADPH by three known isoforms of a heme-containing flavoprotein termed NO synthase (nNOS/NOS-I, iNOS/NOS-II, and eNOS/NOS-III). nNOS is a constitutively expressed neuronal NOS isoform that exits in its latent form until it is activated by the binding of calmodulin follo wing elevation of intracellular calcium levels. The C-terminus of nNOS contains a conserved serine residue, Ser1417, analogous to Ser1177 of the constitutively expressed endothelial NOS isoform (eNOS). Phosphorylation of Ser1417 is believed to regulate nNOS activation, particularly in glucose-sensing neurons, where inhibition of AMPK pathways by glucose and leptin serve to suppress nNOS activity, whereas activation of AMPK by insulin leads to nNOS activation.

This antibody is covered by our Worry-Free Guarantee.

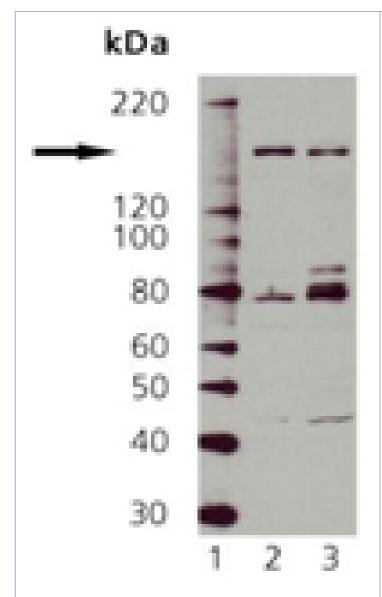
Ordering Information

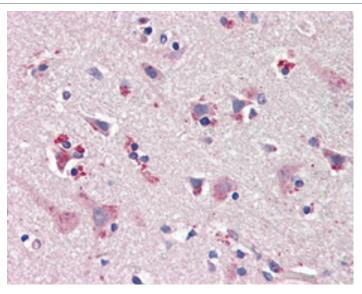
Order Online »

ADI-KAP-NO030-D	50µg
ADI-KAP-NO030-F	200μg

Manuals, SDS & CofA

View Online »





IHC Analysis: Human brain, cortex tissue stained with nNOS, pAb at 20µg/ml.

Western blot analysis: Lane 1: MW marker, Lane 2: Rat Brain, Lane 3: Mouse Brain.

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 Nitric oxide synthase (neuronal), NOS I

Application IHC (PS), WB

Application NotesDetects a band of ~155kDa by Western blot.

Formulation Liquid. In PBS, pH 7.2, containing 50% glycerol and 0.09%

sodium azide.

Host Rabbit

Immunogen Synthetic peptide corresponding to the sequence near the

C-terminus of human nNOS.

Purity Detail Protein A affinity purified.

Recommendation Dilutions/ConditionsWestern Blot (1:250, ECL)Suggested dilutions/conditions

may not be available for all applications. Optimal conditions

must be determined individually for each application.

Source Purified from rabbit serum.

Species Reactivity Mouse, Rat

UniProt ID P29475

Worry-free Guarantee This antibody is covered by our Worry-Free Guarantee



European Sales Office ENZO LIFE SCIENCES (ELS) AG Phone: +41 61 926 8989 infoeu@enzolifesciences.com Belgium, The Netherlands & Luxembourg Phone: +32 3 466 0420 infobe@enzolifesciences.com

France
Phone: +33 472 440 655
infofr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
infode@enzolifesciences.com

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
infouk@enzolifesciences.com