

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 L-arginine, 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 exists in its latent form until it is activated by the binding of calmodulin following 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](#).

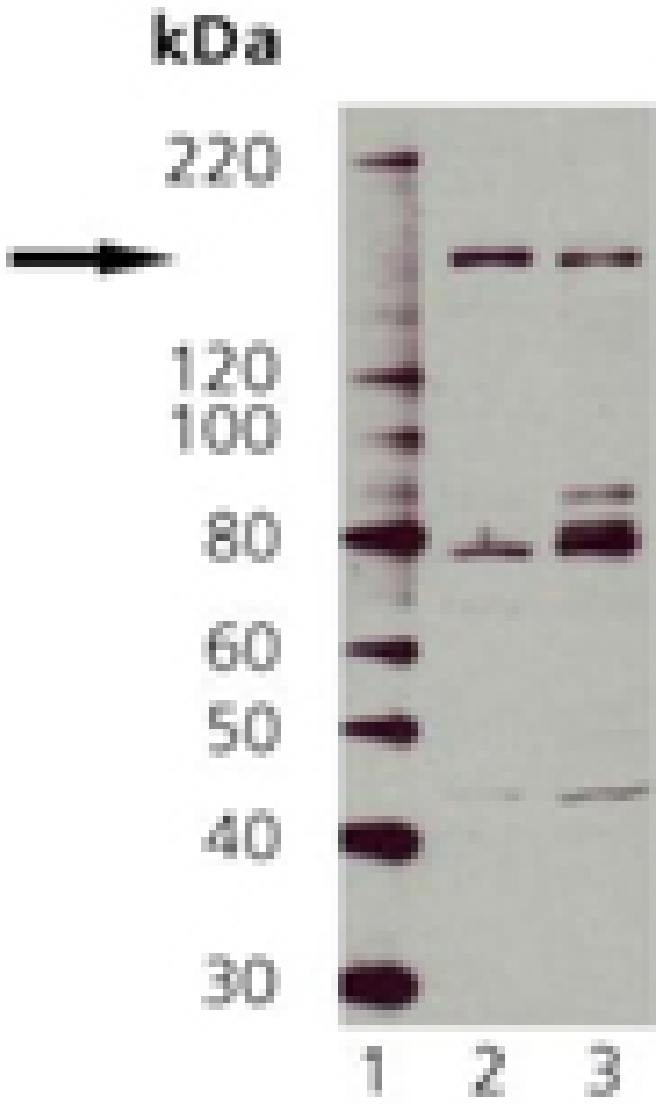
Ordering Information

[Order Online »](#)

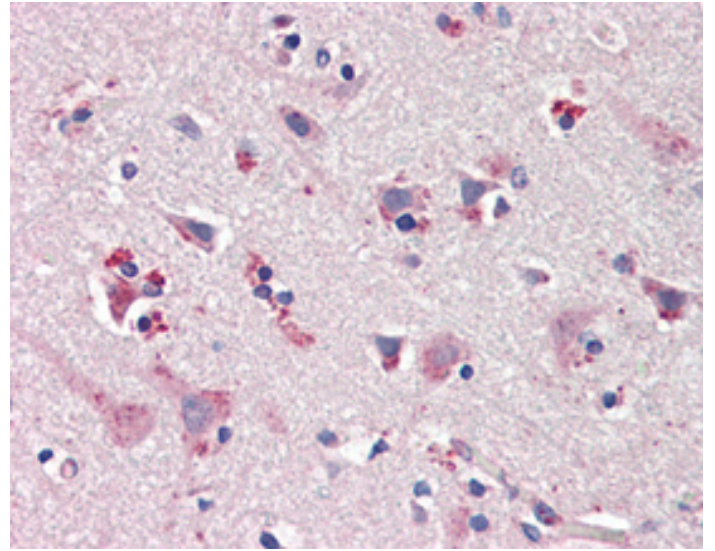
ADI-KAP-NO030-D	50µg
-----------------	------

Manuals, SDS & CofA

[View Online »](#)



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



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

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 Notes	Detects 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/Conditions	Western 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](#)



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)

Last modified: May 29, 2024