

# eNOS 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.

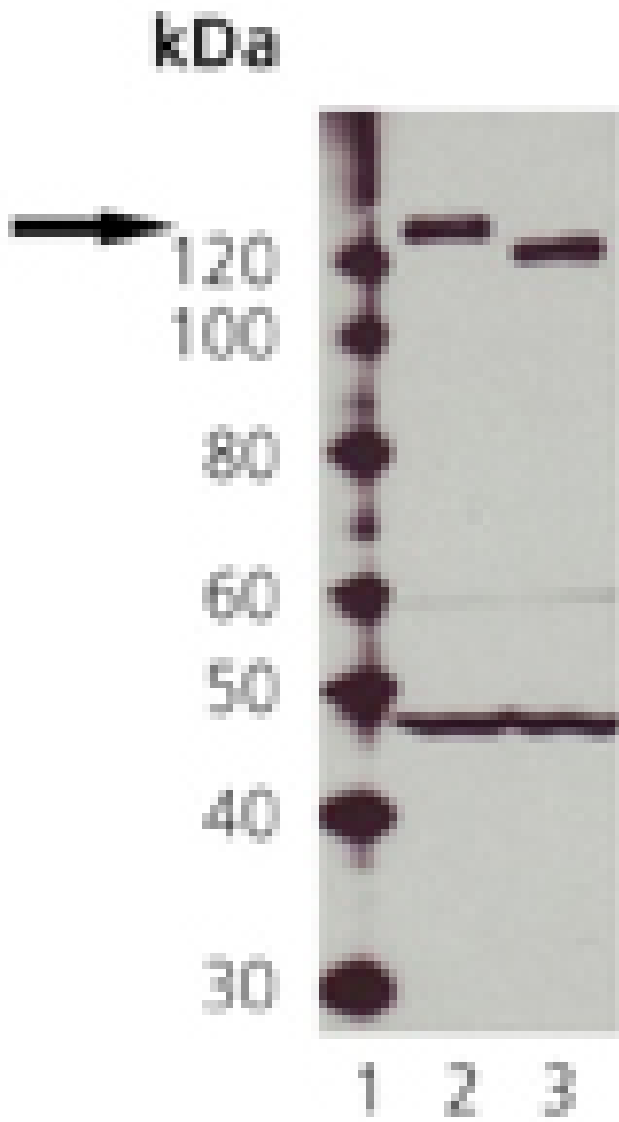
This antibody is covered by our [Worry-Free Guarantee](#).

Citations: 6

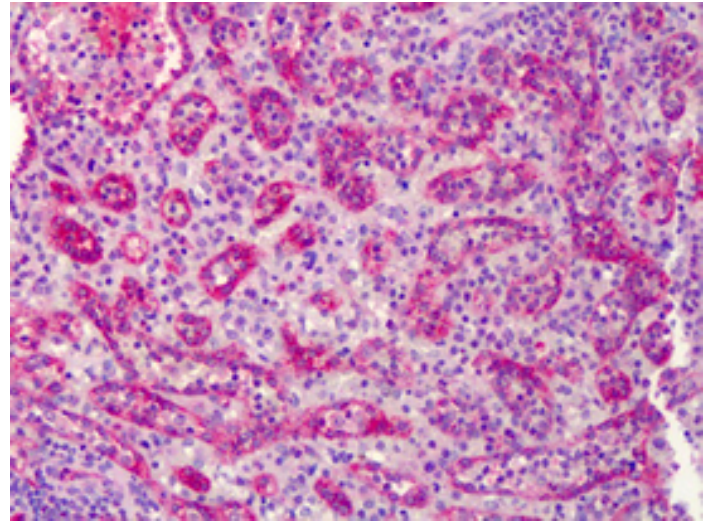
[View Online »](#)

Manuals, SDS & CofA

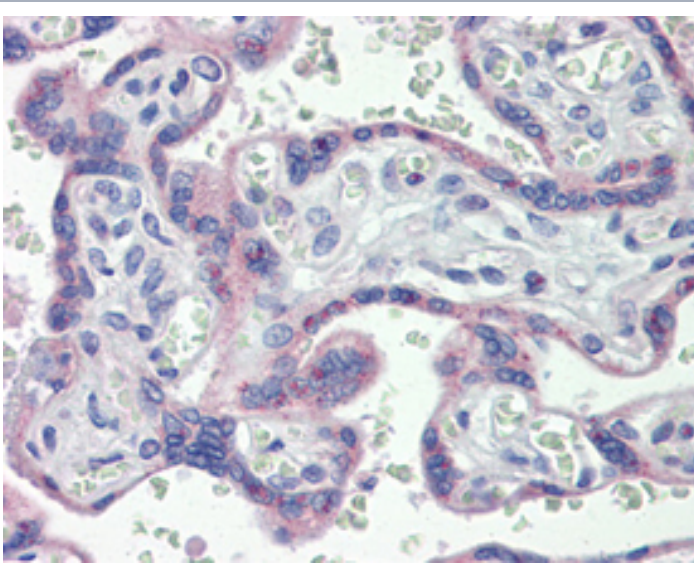
[View Online »](#)



Western blot analysis: Lane 1 MW marker, Lane 2: Rat Brain, and Lane 3: Mouse Brain probed with eNOS, pAb.



IHC Analysis: Human spleen tissue stained with eNOS, pAb at 10µg/ml.



IHC Analysis: Human placenta tissue stained with eNOS, pAb at 10 $\mu$ 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** NOS3, cNOS, NOS III, Nitric oxide synthase (endothelial)

**Application** IHC (PS), WB

**Application Notes** Detects a band of ~135kDa 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 eNOS.

**Purity Detail** Peptide affinity purified.

**Recommendation Dilutions/Conditions** Western Blot (1:1,000, 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** Human, Mouse, Rat

**UniProt ID** P29474

**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-  
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

France  
Phone: +33 472 440 655  
[info-  
de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

Germany  
Phone: +49 7621 5500 526  
[info-  
uk@enzolifesciences.com](mailto:info-uk@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