

eNOS polyclonal antibody

Nitric oxide (NO) is synthesized from L-arginine, oxygen, and NADPH by three known isoforms of heme-containing flavoproteins termed NO synthases (NOS I-III). Analysis of cDNA clones has identified three distinct NOS genes in mammals: neuronal (nNOS/NOS-I), endothelial (eNOS/NOS-III), and inducible (iNOS/NOS-II). eNOS and nNOS are constitutively expressed, agonist-triggered, Ca^{2+} /calmodulin-dependent, and are inhibited by L-arginine analogues (L-NNA, L-NMMA). They are expressed in endothelium, adrenal glands, brain, and platelets. iNOS is inducible, Ca^{2+} /calmodulin-independent, and also inhibited by NMMA and L-NNA. It is expressed in macrophages, hepatocytes, tumor cells, vascular smooth muscle and endothelial cells.

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

Citations: 5

[View Online »](#)

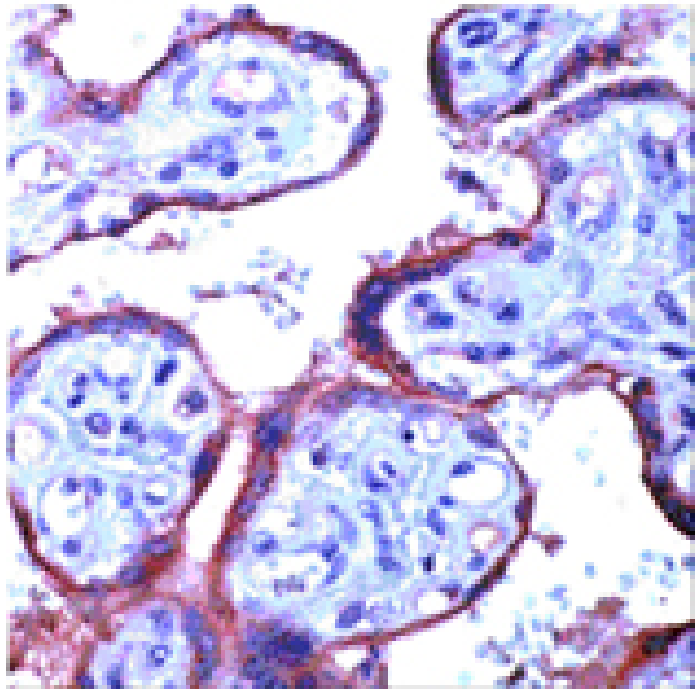
Ordering Information

[Order Online »](#)

ADI-905-386-1	1ml
---------------	-----

Manuals, SDS & CofA

[View Online »](#)



immunohistochemistry analysis of human placenta
stained with eNOS, pAb.

Handling & Storage

Long Term Storage +4°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name	NOS3, cNOS, NOS III, Nitric oxide synthase (endothelial)
Application	IHC
Formulation	Liquid. In PBS, pH 7.4, containing BSA and sodium azide.
Gene/Protein Identifier	4846 (Entrez GeneID)
Host	Rabbit
Immunogen	Synthetic peptide corresponding to the sequence near the C-terminus of human eNOS. The sequence is conserved in mouse, rat, bovine, dog and pig.
Purity Detail	Affinity purified.
Recommendation Dilutions/Conditions	Immunohistochemistry (1:70)Suggested dilutions/conditions may not be available for all applications.Optimal conditions must be determined individually for each application.
Species Reactivity	Bovine, Dog, Human, Mouse, Porcine, Rat
UniProt ID	P29474
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-
fr@enzolifesciences.com](mailto:info-fr@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)