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, Ca2+/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, Ca2+/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

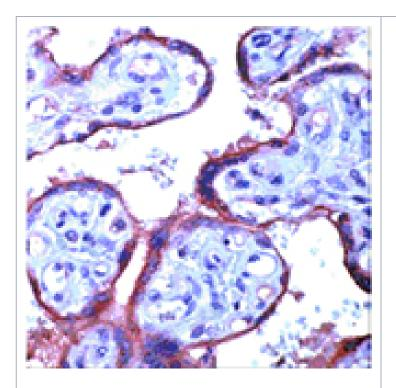
Order Online »

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 GenelD)

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 ReactivityBovine, 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 infousa@enzolifesciences.com 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