
NOSTRIN Polyclonal Antibody

(Catalog # A72096)

Background

Nitric oxide (NO) is a potent mediator in biologic processes such as neurotransmission, inflammatory response, and vascular homeostasis. NOSTRIN binds the enzyme responsible for NO production, endothelial NO synthase (ENOS; MIM 163729), and triggers the translocation of ENOS from the plasma membrane to vesicle-like subcellular structures, thereby attenuating ENOS-dependent NO production.

Description

NOSTRIN Polyclonal Antibody. Unconjugated. Raised in: Rabbit.

Formulation

Buffer: PBS with 0.01% thimerosal, 50% glycerol, pH7.3.

Specificity

Human, Mouse, Rat

Isotype

IgG

Uniprot ID

Q8IIV9

Purification

Affinity Purification

Immunogen

Recombinant fusion protein containing a sequence corresponding to amino acids 207-506 of human NOSTRIN (NP_001034813.2).

Storage

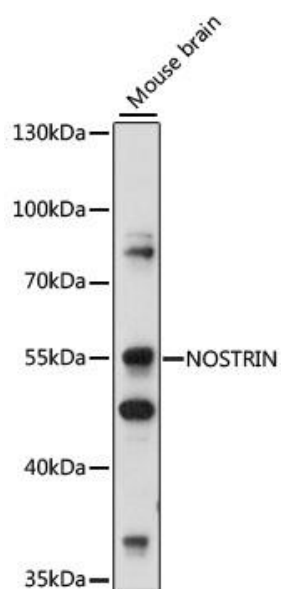
Shipped at 4°C. Upon receipt, store at -20°C. Avoid freeze / thaw cycles

Alternative Names

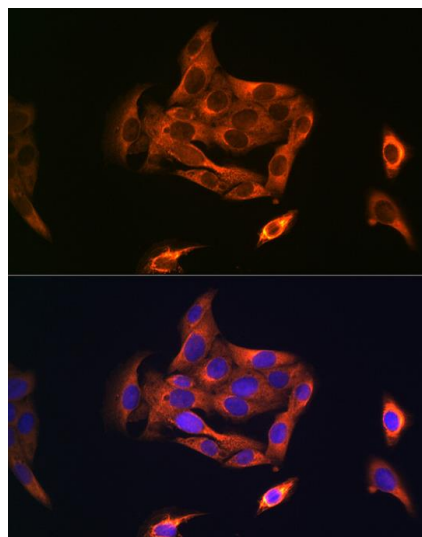
NOSTRIN; DaIP2; nostrin

Application

WB, IF; Recommended dilution: WB 1:500 - 1:1000, IF/ICC 1:50 - 1:200



Western blot analysis of extracts of various cell lines, using NOSTRIN antibody at 1:1000 dilution.
Secondary antibody: HRP Goat Anti-Rabbit IgG (H+L) (AS014) at 1:10000 dilution.
Lysates/proteins: 25µg per lane.
Blocking buffer: 3% nonfat dry milk in TBST.



Immunofluorescence analysis of U-2 OS cells using NOSTRIN Rabbit pAb at dilution of 1:100 (40x lens). Blue: DAPI for nuclear staining.