
NOX4 Polyclonal Antibody

(Catalog # A72493)

Background

This gene encodes a member of the NOX family of enzymes that functions as the catalytic subunit of the NADPH oxidase complex. The encoded protein is localized to non-phagocytic cells, where it acts as an oxygen sensor and catalyzes the reduction of molecular oxygen to various reactive oxygen species (ROS). The ROS generated by this protein has been implicated in numerous biological functions, including signal transduction, cell differentiation, and tumor cell growth. A pseudogene has been identified on the other arm of chromosome 11. Alternative splicing results in multiple transcript variants.

Description

NOX4 Polyclonal Antibody. Unconjugated. Raised in: Rabbit.

Formulation

Buffer: PBS with 0.09% Sodium azide, 50% glycerol, pH7.3.

Specificity

Human, Mouse, Rat

Isotype

IgG

Uniprot ID

Q9NPH5

Purification

Affinity Purified

Immunogen

A synthetic peptide corresponding to a sequence within amino acids 328-578 of human NADPH oxidase 4 (NOX4) (NP_058627.1).

Storage

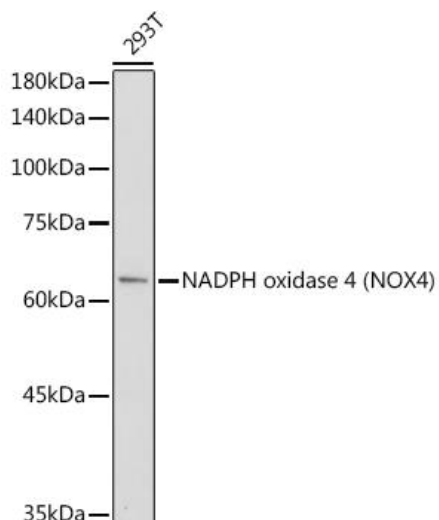
Shipped at 4°C. Upon receipt, store at -20°C. Avoid freeze / thaw cycles.

Alternative Names

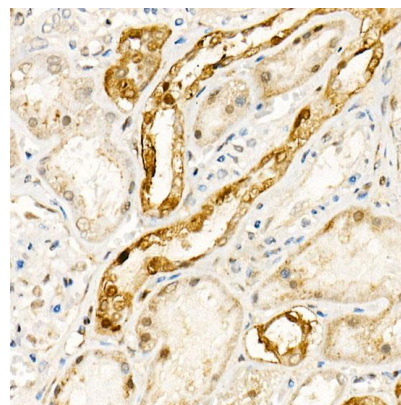
NOX4; KOX; KOX-1; RENOX; NADPH oxidase 4

Application

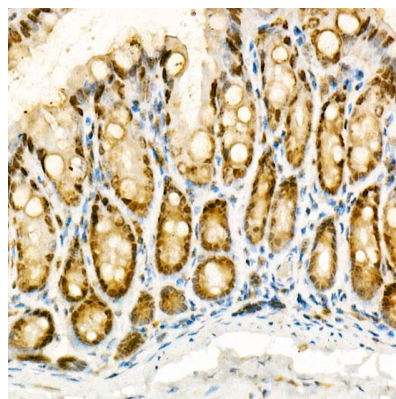
WB, IHC, ELISA; Recommended dilution: WB 1:500 - 1:1000, IHC 1:50 - 1:200, ELISA - Recommended starting concentration is 1 µg/mL. Please optimize the concentration based on your specific assay requirements.



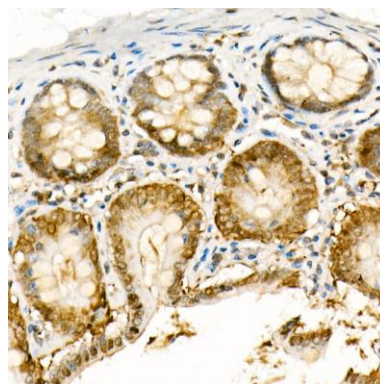
Western blot analysis of extracts of 293T cells, using NOX4 Polyclonal Antibody at 1:1000 dilution.
 Secondary antibody: HRP Goat Anti-Rabbit IgG (H+L) at 1:10000 dilution.
 Lysates/proteins: 25ug per lane.
 Blocking buffer: 3% nonfat dry milk in TBST.
 Exposure time: 90s.



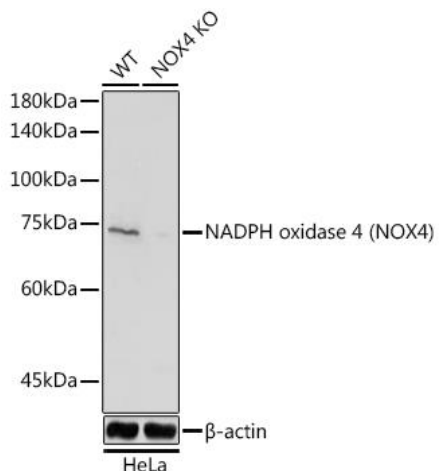
Immunohistochemistry analysis of paraffin-embedded Human kidney using NOX4 Polyclonal Antibody at dilution of 1:100 (40x lens). High pressure antigen retrieval performed with 0.01M Citrate Bufferr (pH 6.0) prior to IHC staining.



Immunohistochemistry analysis of paraffin-embedded Mouse colon using NOX4 Polyclonal Antibody at dilution of 1:100 (40x lens). High pressure antigen retrieval performed with 0.01M Citrate Bufferr (pH 6.0) prior to IHC staining.



Immunohistochemistry analysis of paraffin-embedded Rat colon using NOX4 Polyclonal Antibody at dilution of 1:100 (40x lens). High pressure antigen retrieval performed with 0.01M Citrate Bufferr (pH 6.0) prior to IHC staining.



Western blot analysis of extracts from normal (control) and NADPH oxidase 4 (NOX4) Rabbit pAb knockout (KO) HeLa cells, using NOX4 Polyclonal Antibody at 1:1000 dilution.
 Secondary antibody: HRP Goat Anti-Rabbit IgG (H+L) at 1:10000 dilution.
 Lysates/proteins: 25µg per lane.
 Blocking buffer: 3% nonfat dry milk in TBST.
 Exposure time: 60s.