

## NOX2-gp91phox Polyclonal Antibody

(Catalog #A70477)

### Background

Cytochrome b (-245) is composed of cytochrome b alpha (CYBA) and beta (CYBB) chain. It has been proposed as a primary component of the microbicidal oxidase system of phagocytes. CYBB deficiency is one of five described biochemical defects associated with chronic granulomatous disease (CGD). In this disorder, there is decreased activity of phagocyte NADPH oxidase; neutrophils are able to phagocytize bacteria but cannot kill them in the phagocytic vacuoles. The cause of the killing defect is an inability to increase the cell's respiration and consequent failure to deliver activated oxygen into the phagocytic vacuole.

### Description

NOX2-gp91phox 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

P04839

### Purification

Affinity Purification

### Immunogen

A synthetic peptide corresponding to a sequence within amino acids 385-570 of human NOX2/gp91phox (NP\_000388.2)

### Storage

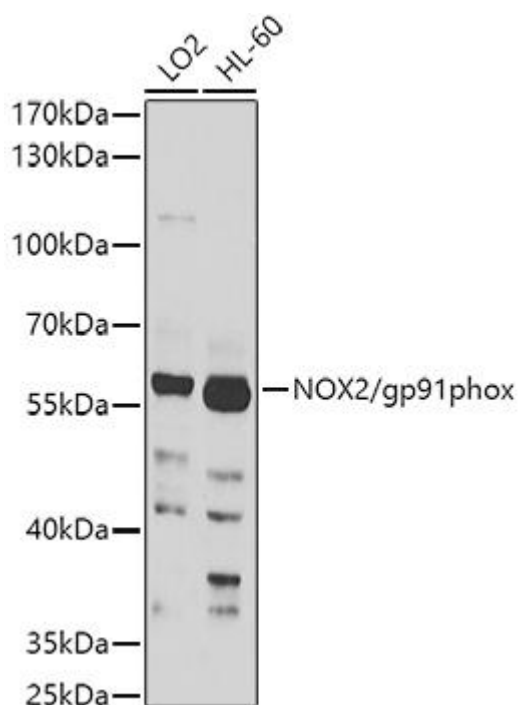
Shipped at 4°C. Upon receipt, store at -20°C. Avoid freeze / thaw cycles

### Alternative Names

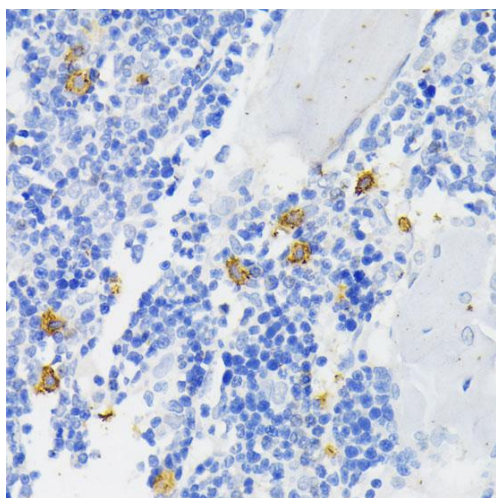
CYBB; AMCBX2; CGD; GP91-1; GP91-PHOX; GP91PHOX; IMD34; NOX2; p91-PHOX; cytochrome b-245 beta chain

### Application

WB, IHC, IF, ELISA; Recommended dilution: WB 1:500 - 1:1000, IHC 1:50 - 1:200, IF 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 various lysates using NOX2-gp91phox Polyclonal Antibody at 1:1000 dilution.  
Secondary antibody: HRP-conjugated 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.



Immunohistochemistry analysis of paraffin-embedded Rat bone marrow using NOX2-gp91phox Polyclonal Antibody at dilution of 1:200 (40x lens). Microwave antigen retrieval performed with 0.01M PBS Buffer (pH 7.2) prior to IHC staining.