

# Phospho-VIM-S39 Polyclonal Antibody

(Catalog # A72599)

## Background

This gene encodes a member of the intermediate filament family. Intermediate filaments, along with microtubules and actin microfilaments, make up the cytoskeleton. The protein encoded by this gene is responsible for maintaining cell shape, integrity of the cytoplasm, and stabilizing cytoskeletal interactions. It is also involved in the immune response, and controls the transport of low-density lipoprotein (LDL)-derived cholesterol from a lysosome to the site of esterification. It functions as an organizer of a number of critical proteins involved in attachment, migration, and cell signaling. Mutations in this gene causes a dominant, pulverulent cataract.

## Description

Phospho-VIM-S39 Polyclonal Antibody. Unconjugated. Raised in: Rabbit.

## Formulation

Buffer: PBS with 0.02% sodium azide, 50% glycerol, pH7.3.

## Specificity

Human, Mouse, Rat

## Isotype

IgG

## Uniprot ID

P08670

## Purification

Affinity Purification

## Immunogen

A synthetic phosphorylated peptide around S39 of human VIM (NP\_003371.2).

## Storage

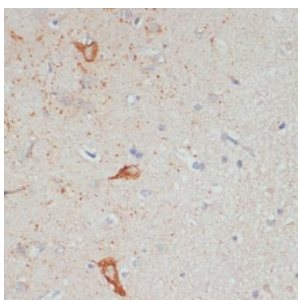
Shipped at 4°C. Upon receipt, store at -20°C. Avoid freeze / thaw cycles

## Alternative Names

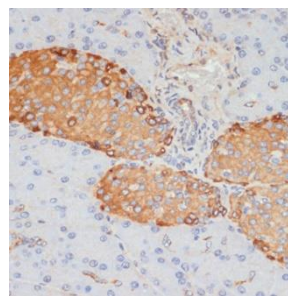
VIM; CTRCT30; HEL113; vimentin

## Application

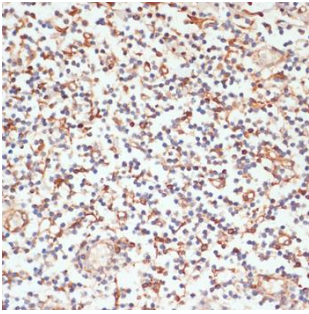
WB, IHC; Recommended dilution: WB 1:500 - 1:2000, IHC 1:50 - 1:100



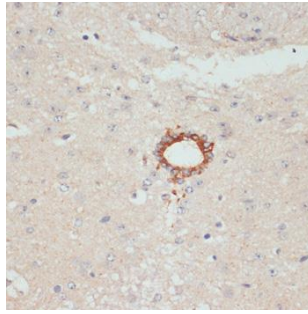
Immunohistochemistry of paraffin-embedded rat brain using Phospho-VIM-S39 Polyclonal Antibody at dilution of 1:100 (40x lens).



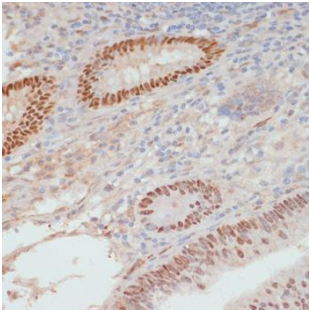
Immunohistochemistry of paraffin-embedded rat pancreas using Phospho-VIM-S39 Polyclonal Antibody at dilution of 1:100 (40x lens).



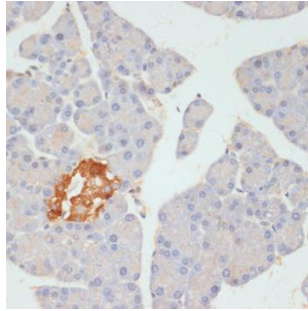
Immunohistochemistry of paraffin-embedded human tonsil using Phospho-VIM-S39 Polyclonal Antibody at dilution of 1:100 (40x lens).



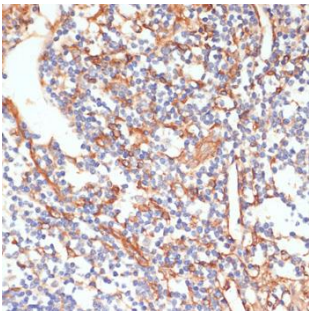
Immunohistochemistry of paraffin-embedded mouse spinal cord using Phospho-VIM-S39 Polyclonal Antibody at dilution of 1:100 (40x lens).



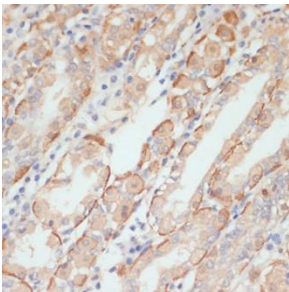
Immunohistochemistry of paraffin-embedded human colon carcinoma using Phospho-VIM-S39 Polyclonal Antibody at dilution of 1:100 (40x lens).



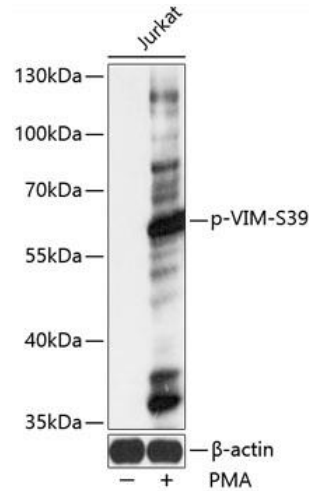
Immunohistochemistry of paraffin-embedded mouse pancreas using Phospho-VIM-S39 Polyclonal Antibody at dilution of 1:100 (40x lens).



Immunohistochemistry of paraffin-embedded human appendix using Phospho-VIM-S39 Polyclonal Antibody at dilution of 1:100 (40x lens).



Immunohistochemistry of paraffin-embedded human stomach using Phospho-VIM-S39 Polyclonal Antibody at dilution of 1:100 (40x lens).



Western blot analysis of extracts of Jurkat cells, using Phospho-VIM-S39 antibody at 1:2000 dilution. Jurkat cells were treated by PMA/TPA (200nM) for 10 minutes.

Secondary antibody: HRP Goat Anti-Rabbit IgG (H+L) at 1:10000 dilution.

Lysates/proteins: 25ug per lane.

Blocking buffer: 3% BSA.

Exposure time: 1s.