
PTPN13 Polyclonal Antibody

(Catalog # A72996)

Background

The protein encoded by this gene is a member of the protein tyrosine phosphatase (PTP) family. PTPs are signaling molecules that regulate a variety of cellular processes including cell growth, differentiation, mitotic cycle, and oncogenic transformation. This PTP is a large intracellular protein. It has a catalytic PTP domain at its C-terminus and two major structural domains: a region with five PDZ domains and a FERM domain that binds to plasma membrane and cytoskeletal elements. This PTP was found to interact with, and dephosphorylate, Fas receptor and I κ B α through the PDZ domains. This suggests it has a role in Fas mediated programmed cell death. This PTP was also shown to interact with GTPase-activating protein, and thus may function as a regulator of Rho signaling pathways. Four alternatively spliced transcript variants, which encode distinct proteins, have been reported.

Description

PTPN13 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

Q12923

Purification

Affinity Purification

Immunogen

Recombinant fusion protein containing a sequence corresponding to amino acids 250-500 of human PTPN13 (NP_542416.1).

Storage

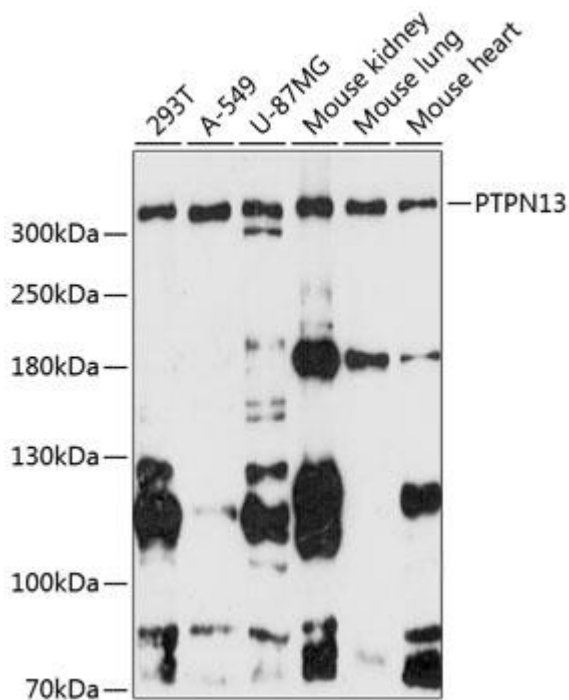
Shipped at 4°C. Upon receipt, store at -20°C. Avoid freeze / thaw cycles

Alternative Names

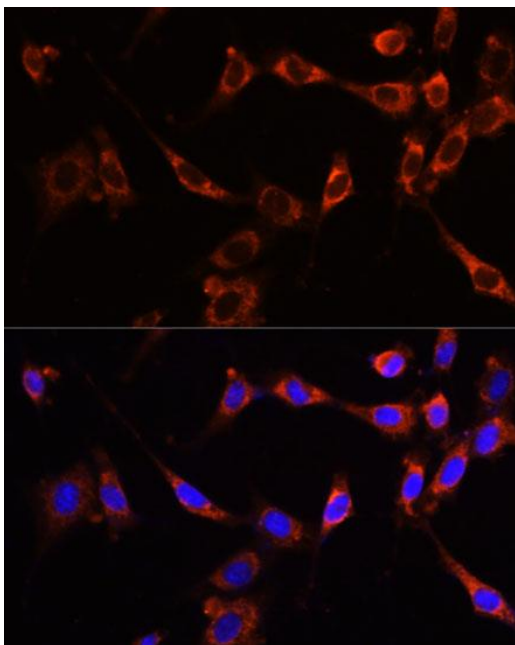
PTPN13; FAP-1; PNP1; PTP-BAS; PTP-BL; PTP1E; PTPL1; PTPLE; hPTP1E; tyrosine-protein phosphatase non-receptor type 13

Application

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



Western blot analysis of extracts of various cell lines, using PTPN13 Polyclonal Antibody at 1:3000 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.



Immunofluorescence analysis of NIH-3T3 cells using PTPN13 Polyclonal Antibody at dilution of 1:100 (40x lens). Blue: DAPI for nuclear staining.