
SNX13 Polyclonal Antibody

(Catalog #A75633)

Background

This gene encodes a PHOX domain- and RGS domain-containing protein that belongs to the sorting nexin (SNX) family and the regulator of G protein signaling (RGS) family. The PHOX domain is a phosphoinositide binding domain, and the SNX family members are involved in intracellular trafficking. The RGS family members are regulatory molecules that act as GTPase activating proteins for G alpha subunits of heterotrimeric G proteins. The RGS domain of this protein interacts with G alpha(s), accelerates its GTP hydrolysis, and attenuates G alpha(s)-mediated signaling. Overexpression of this protein delays lysosomal degradation of the epidermal growth factor receptor. Because of its bifunctional role, this protein may link heterotrimeric G protein signaling and vesicular trafficking.

Description

SNX13 Polyclonal Antibody. Unconjugated. Raised in: Rabbit.

Formulation

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

Specificity

Human, Mouse

Isotype

IgG

Uniprot ID

Q9Y5W8

Purification

Affinity Purified

Immunogen

Recombinant protein (or fragment). Please contact EpigenTek for more information.

Storage

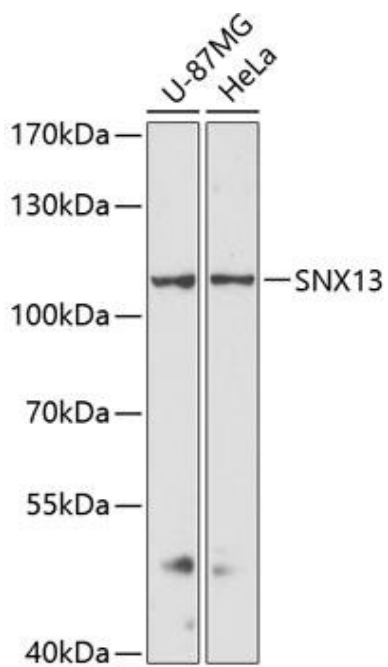
Store at -20°C. Avoid freeze / thaw cycles.

Alternative Names

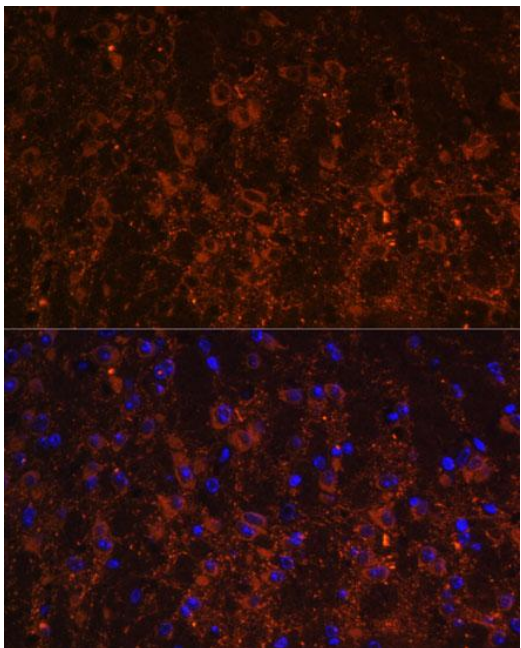
RGS-PX1; SNX13

Application

WB, IF-P, ELISA; Recommended dilution: WB: 1:500 - 1:2000; IF-P: 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 SNX13 Polyclonal Antibody at 1:3000 dilution.
Secondary antibody: HRP-conjugated Goat anti-Rabbit IgG (H+L) at 1:10000 dilution.
Lysates/proteins: 25 μ g per lane.
Blocking buffer: 3% nonfat dry milk in TBST.
Exposure time: 90s.



Immunofluorescence analysis of paraffin-embedded mouse brain using SNX13 Polyclonal Antibody at dilution of 1:100.
Secondary antibody: Cy3-conjugated Goat anti-Rabbit IgG (H+L) at 1:500 dilution. Blue: DAPI for nuclear staining.