
RHOH Polyclonal Antibody

(Catalog # A72661)

Background

The protein encoded by this gene is a member of the Ras superfamily of guanosine triphosphate (GTP)-metabolizing enzymes. The encoded protein is expressed in hematopoietic cells, where it functions as a negative regulator of cell growth and survival. This gene may be hypermutated or misexpressed in leukemias and lymphomas. Chromosomal translocations in non-Hodgkin's lymphoma occur between this locus and B-cell CLL/lymphoma 6 (BCL6) on chromosome 3, leading to the production of fusion transcripts. Alternative splicing in the 5' untranslated region results in multiple transcript variants that encode the same protein.

Description

RHOH Polyclonal Antibody. Unconjugated. Raised in: Rabbit.

Formulation

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

Specificity

Human, Mouse

Isotype

IgG

Uniprot ID

Q15669

Purification

Affinity Purification

Immunogen

Recombinant fusion protein containing a sequence corresponding to amino acids 1-191 of human RHOH (NP_004301.1).

Storage

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

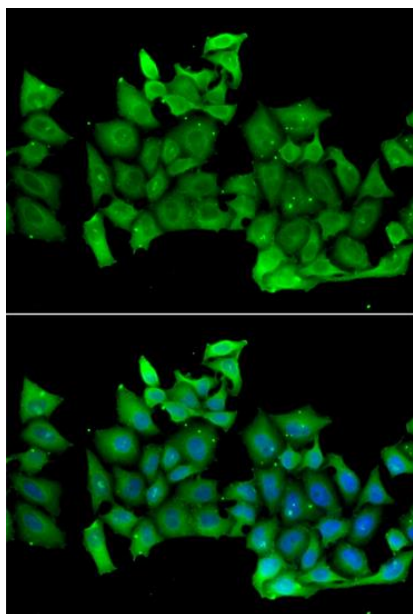
Alternative Names

RHOH; ARHH; TTF; ras homolog family member H

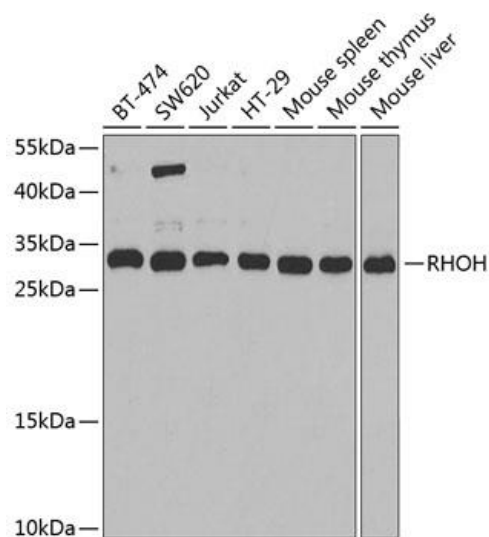
Application

WB, IF; Recommended dilution: WB: 1:500-1:2000, IF:1:10-1:100

Exposure time: 90s.



Immunofluorescence analysis of MCF-7 cells using RHOH Polyclonal Antibody. Blue: DAPI for nuclear staining.



Western blot analysis of extracts of various cell lines, using RHOH 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.