

PRDM6 Polyclonal Antibody

(Catalog # A-2006)

Background

PRDM6 is a member of PR domain family of proteins associated with gene regulation and chromatin remodeling. It functions as a transcriptional repressor through interactions with Histone H4 deacetylases and methyltransferases. Expressed in vascular precursor cells, PRDM6 is believed to play a role in the development and differentiation of vascular tissue.

Description

PRDM6 Polyclonal Antibody. Unconjugated. Raised in: Rabbit.

Formulation

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

Specificity

Human, Mouse, Rat

Isotype

IgG

Uniprot ID

Q9NQX0

Purification

Affinity Purified

Immunogen

Recombinant Protein of Human PRDM6

Storage

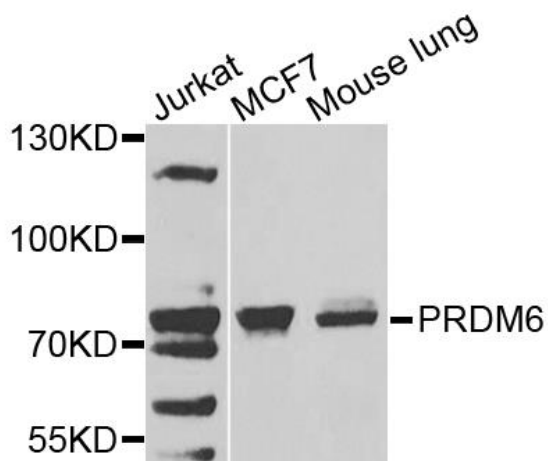
Shipped at 4°C. Store at -20°C. Avoid multiple freeze/thaw cycles.

Alternative Names

Gm92 antibody, PFM3 antibody, PR domain containing 6 antibody, PR domain zinc finger protein 6 antibody, Prdm6 antibody, PRISM antibody, Putative histone-lysine N-methyltransferase PRDM6 antibody, RGD1559594 antibody

Application

WB; Recommended dilution: WB: 1:500 - 1:2000



Western blot analysis of extracts of various cells, using PRDM6 Polyclonal Antibody.