

BLIMP1 / PRDM1 antibody (pAb)

Catalog Nos: 61053, 61054

RRID: AB_2793489

Isotype: IgG

Application(s): IF, WB

Reactivity: Human

Volumes: 100 μ l, 10 μ l

Purification: Affinity Purified

Host: Rabbit

Molecular Weight: 110 kDa

Background: **BLIMP1 / PRDM1** (Beta-interferon gene positive regulatory domain I-binding factor, PR domain-containing protein 1 Positive regulatory domain I-binding factor 1) is a DNA sequence-specific transcriptional repressor that binds specifically to the PRDI element in the promoter of the beta-interferon gene, as well as other genes. It is required for maturation of B-lymphocytes into immunoglobulin secreting plasma cells through direct repression of transcription factors such as Pax5, Bcl6, and Myc, which function in the lineage commitment and proliferation of B cells.

BLIMP1 / PRDM1 contains an N-terminal SET domain and four C-terminal C2H2-type zinc-finger motifs. It is also required for effector T-cell function and for the specification of primordial germ cell lineages.

Immunogen: This BLIMP1 / PRDM1 antibody was raised against a peptide within the N-terminal half of human BLIMP1 / PRDM1.

Buffer: Purified IgG in 70 mM Tris (pH 8), 105 mM NaCl, 31 mM glycine, 0.07 mM EDTA, 30% glycerol and 0.035% sodium azide. Sodium azide is highly toxic.

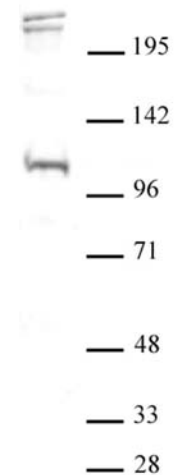
Application Notes:

Applications Validated by Active Motif:

WB: 1:500 - 1:1,000 dilution

Storage and Guarantee: Some products may be shipped at room temperature. This will not affect their stability or performance. Avoid repeated freeze/thaw cycles by aliquoting items into single-use fractions for storage at -20°C for up to 2 years. Keep all reagents on ice when not in storage. This product is guaranteed for 12 months from date of receipt.

This product is for research use only and is not for use in diagnostic procedures.



Western blot of BLIMP1 / PRDM1 antibody.
Nuclear extract of Saos-2 cells (20 μ g per lane) probed with BLIMP1 / PRDM1 antibody (pAb) at a dilution of 1:1,000.