



PRDM1 (BLIMP1) Polyclonal Antibody

Product Specifications	
Catalog Number:	905-697-100
Host:	Rabbit
Species Reactivity:	Human
Applications: <i>The optimal dilution for a specific application must be determined by the investigator</i>	ELISA: 1:1,000
Predicted m.w:	~88 kDa
Concentration:	See product label
Purification:	Protein G Affinity
Format:	PBS with 0.09% (W/V) sodium azide
Storage:	Store at -20°C
Immunogen:	Purified recombinant GST fusion protein corresponding to the amino terminus of human BLIMP1

Background:

This gene encodes a protein that acts as a repressor of beta-interferon gene expression. The protein binds specifically to the PRDI (positive regulatory domain I element) of the IFN β gene promoter. Transcription of this gene increases upon virus induction. Two alternatively spliced transcript variants that encode different isoforms have been reported. The protein encoded by this transcript has also been named PRDI-BF1 α .

References:

1. Gory, I., *et al.* (2003) *J Immunol.* **170**, 3125-3133.
2. Borson, N.D., *et al.* (2002) *Blood* **100**, 4629-4639.
3. Vasanwala, F.H., *et al.* (2002) *J Immunol.* **169**, 922-1929.
4. Shaffer, A.L., *et al.* (2002) *Immunity* **17**, 51-62.
5. Medina, F., *et al.* (2002) *Blood* **99**, 2154-2161.