

PRDM12 Antibody [#C11694]

Catalog Number:	C11694
Concentration:	1mg/ml
Swiss-Prot No.:	Q9H4Q4
Other Names:	PR domain-containing protein 12, PR domain zinc finger protein 12, PRDM12, PFM9
Storage/Stability:	Store at -20°C/1 year
Form of Antibody:	Rabbit IgG in phosphate buffered saline (without Mg ²⁺ and Ca ²⁺), pH 7.4, 150mM NaCl, 0.02% sodium azide and 50% glycerol.
Immunogen:	The antiserum was produced against synthesized peptide derived from internal of human PRDM12.
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	PRDM12 antibody detects endogenous levels of total PRDM12 protein.
Reactivity:	Human, Mouse
Applications:	WB: 1:500~1:1000 ELISA: 1:20000
References:	Yang X.-H., Submitted (JUL-2001) to the EMBL/GenBank/DDBJ databases.