

PRDM12 Polyclonal Antibody

(Catalog #A-2012)

Background

PRDM12 belongs to PR-domain family of proteins associated with cell differentiation and cancer. It is believed to play a role in transcription regulation and act as a tumor suppressor. Inactivation of this gene can happen through the loss of mRNA expression, frame shift mutation, missense mutation, and chromosomal deletion.

Description

PRDM12 Polyclonal Antibody. Unconjugated. Raised in: Rabbit.

Formulation

Liquid in PBS containing 50% glycerol, 0.5% BSA* and 0.02% sodium azide.

Specificity

Human, Mouse

Isotype

IgG

Uniprot ID

Q9H4Q4

Purification

Affinity Purified

Immunogen

Synthesized peptide derived from the Internal region of Human PRDM12

Storage

Shipped at 4°C (Blue Ice). Upon receipt, store at -20°C or -80°C. Avoid repeated freeze.

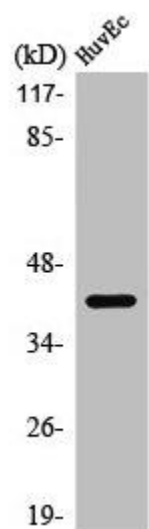
Alternative Names

PRDM12; PFM9; PR domain zinc finger protein 12; PR domain-containing protein 12

Application

WB, IF, ELISA; Recommended dilution: WB:1/500-1/2000, IF:1/200-1/1000, ELISA:1/20000

**The BSA is derived from animal or animal-derived material of negligible amounts. The animal-derived material is subject to heat treatment at a temperature higher than 65°C for at least three hours and acid treatment with pH value of less than 5 for at least three hours, thereby being free of Bovine Spongiform Encephalopathy (BSE) and Transmissible Spongiform Encephalopathy (TSE). Additionally, as all BSA-containing products are strictly intended for research purposes and not for diagnostic or therapeutic use, they are not subject to certification authority oversight.*



Western Blot analysis of HuvEC cells using PRDM12 Polyclonal Antibody