

Recombinant PRDM4 protein

Catalog No: 81395, 81695**Expressed In:** Baculovirus**Quantity:** 20, 1000 µg**Concentration:** 0.2 µg/µl**Source:** Human

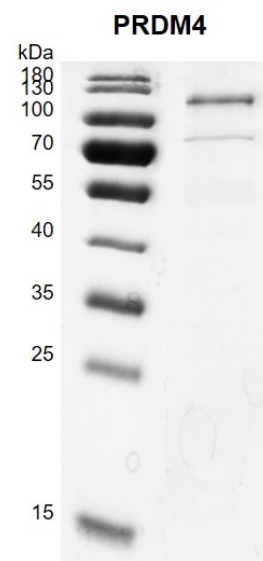
Buffer Contents: Recombinant PRDM4 protein is supplied in 25 mM HEPES-NaOH pH 7.5, 300 mM NaCl, 10% glycerol, 0.04% Triton X-100, 0.5 mM TCEP.

Background: PRDM4 (PR/SET Domain 4) is a transcription factor of the PR-domain protein family. It contains a PRdomain and multiple zinc finger motifs. Transcription factors of the PR-domain family are known to be involved in cell differentiation and tumorigenesis. PRDM4 recruits the chromatin modifier PRMT5, an arginine methyltransferase that catalyzes symmetric dimethylation of histone H4 arginine 3 (H4R3me2s) to exerts histone methyltransferase activity. An elevated expression level of this gene has been observed in PC12 cells treated with nerve growth factor, beta polypeptide (NGF). Diseases associated with PRDM4 include Pseudo-Von Willebrand Disease and Porokeratosis.

Protein Details: Recombinant PRDM4 protein was expressed in baculovirus expression system as the full length protein (accession number NP_036538.3) with a N-terminal FLAG tag. The molecular weight of the protein is 89.2 kDa.

Application Notes: This product was manufactured as described in Protein Details. Where possible, Active Motif has developed functional or activity assays for recombinant proteins. Additional characterization such as enzyme kinetic activity assays, inhibitor screening or other biological activity assays may not have been performed for every product. All available data for a given product is shown on the lot-specific Technical Data Sheet.

Storage and Guarantee: Recombinant proteins in solution are temperature sensitive and must be stored at -80°C to prevent degradation. Avoid repeated freeze/thaw cycles and keep on ice when not in storage. This product is for research use only and is not for use in diagnostic procedures. This product is guaranteed for 6 months from date of arrival.

**Recombinant PRDM4 protein**

10% SDS-PAGE Coomassie staining

MW: 89.2 kDa

Purity: ≥ 70%