

## Recombinant PRDM10 protein

**Catalog No:** 31396, 31796

**Expressed In:** Baculovirus

**Quantity:** 20, 1000 µg

**Concentration:** 0.4 µg/µl

**Source:** Human

**Buffer Contents:** Full length recombinant PRDM10 protein expressed in Sf9 cells at a concentration of 0.4 µg/µl in 25 mM HEPES, pH 7.5, 300 mM NaCl, 0.2 mg/ml 3X FLAG peptide and 5% glycerol.

**Background: PRDM10 (PR Domain Containing 10)** is a member of the PRDI-BF1 and RIZ homology domain containing (PRDM) family of transcriptional regulators. PRDM10 is a transcription factor that contains C2H2-type zinc-fingers and that possesses histone methyltransferase activity. PRDM10 also contains a positive regulatory domain which has been found in several other zinc-finger transcription factors including those involved in B cell differentiation and tumor suppression. Studies of the mouse counterpart suggest that this protein may be involved in the development of the central nerve system (CNS), as well as in the pathogenesis of neuronal storage disease. Multiple alternatively spliced transcript variants encoding distinct isoforms have been observed.

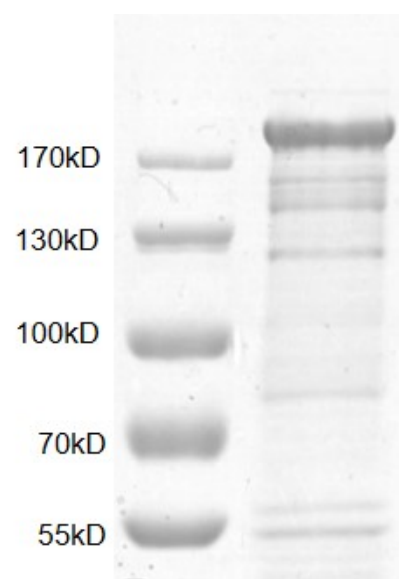
**Protein Details:** Recombinant PRDM10 (accession number NP\_004982.2) was expressed in Sf9 cells and contains an N-terminal FLAG tag with an observed molecular weight of 134 kDa (appears larger in SDS-PAGE gel image). The recombinant protein is >70% pure by SDS-PAGE.

**Application Notes:** Recombinant PRDM10 is suitable for use in the study of enzyme kinetics, inhibitor screening, and selectivity profiling.

**Specific Activity:** Unknown.

**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 guaranteed for 6 months from date of receipt.

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



**Recombinant PRDM10 protein gel.** PRDM10 protein was run on an SDS-PAGE gel and stained with Coomassie Blue.