

## Recombinant PRDM14 protein

**Catalog No:** 31397, 31797

**Expressed In:** Baculovirus

**Quantity:** 20 µg

**Concentration:** 0.15 µg/µl

**Source:** Human

**Buffer Contents:** Recombinant PRDM14 protein is supplied at a concentration of 0.15 µg/µl in 25 mM HEPES pH 7.5, 300 mM NaCl, 5% glycerol, 0.04% Triton X-100, 0.2 mM TCEP.

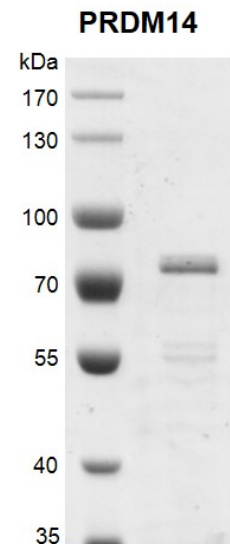
**Background: PRDM14 (PR Domain Containing 14)** is a member of the PRDI-BF1 and RIZ homology domain containing (PRDM) family of transcriptional regulators. PRDM14 possesses histone methyltransferase activity. PRDM14 is a transcription factor that has both positive and negative roles on transcription. It is required for the maintenance of embryonic stem cell identity and the reacquisition of pluripotency in somatic cells. PRDM14 may play an essential role in germ cell development at two levels: the reacquisition of potential pluripotency, including Sox2 up-regulation, and successful epigenetic reprogramming, characterized by EHMT1 repression. PRDM14 directly up-regulates the expression of pluripotency gene Oct4 through its proximal enhancer. PRDM14 binds to the DNA consensus sequence 5'-GGTC[TC]CTAA-3.'

**Protein Details:** Recombinant PRDM14 was expressed in a baculovirus expression system as the full length protein (accession number NP\_004982.2) with an N-terminal FLAG tag. The molecular weight of PRDM14 is 67 kDa. The purity of the protein is > 80% by SDS-PAGE.

**Application Notes:** Recombinant PRDM14 is suitable for use in the study of enzyme kinetics, inhibitor screening, and selectivity profiling.

**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 PRDM14 protein gel.** PRDM14 protein was run on an 8% SDS-PAGE gel and stained with Coomassie Blue.