

Recombinant PRDM6 protein

Catalog No: 31495, 31899

Expressed In: Baculovirus

Quantity: 20, 1000 µg

Concentration: 0.2 µg/µl

Source: Human

Buffer Contents: Full length recombinant PRDM6 protein was expressed in Sf9 cells and is supplied at a concentration of 0.2 µg/µl in 25 mM Hepes pH 7.5, 300 mM NaCl, 5% Glycerol, 0.04% Triton X-100, 0.2 mM TCEP.

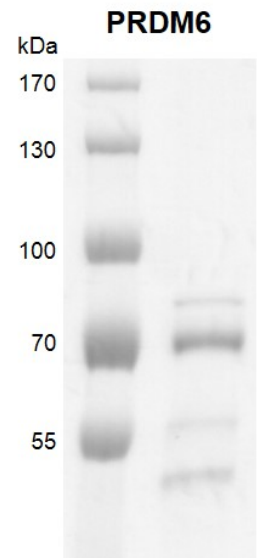
Background: PRDM6 (PR Domain Containing 6) is a putative histone methyltransferase that acts as a transcriptional repressor of smooth muscle gene expression. PRDM6 can regulate the phenotypic plasticity of smooth muscle cells by suppressing differentiation and maintaining the proliferative potential of vascular SMCs. It also plays a role in endothelial cells by inhibiting endothelial cell proliferation, survival and differentiation.

Protein Details: Recombinant PRDM6 protein (accession number NP_001129711.1) was expressed in Sf9 cells and contains an N-terminal FLAG tag with a molecular weight of 64.4 kDa. The recombinant protein is >75% pure by SDS-PAGE.

Application Notes: Recombinant PRDM6 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 PRDM6 protein gel. Recombinant PRDM6 run on an 8% SDS-PAGE gel and stained with Coomassie blue.