

Recombinant PRDM11 protein

Catalog No: 81278, 81978

Expressed In: Baculovirus

Quantity: 20, 1000 µg

Concentration: 0.5 µg/µl

Source: Human

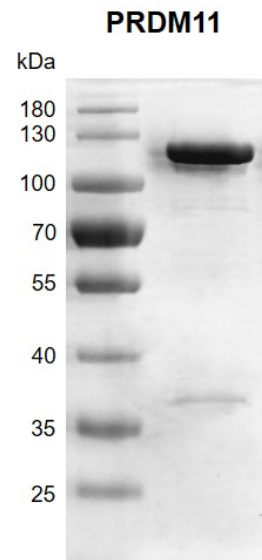
Buffer Contents: Recombinant PRDM11 protein is supplied in 25 mM HEPES-NaOH pH 7.5, 300 mM NaCl, 10% glycerol, 0.04% Triton X-100, 0.5 mM TCEP.

Background: PRDM11 (PR/SET Domain 11), also called as PFM8, is a member of the PR-domain family and acts as a transcription regulator. Some reported that PRDM11 is a putative tumor suppressor that controls the expression of key oncogenes. It's been demonstrated that overexpression of PRDM11 inhibits cell proliferation and induces apoptosis. PRDM11 may also regulate the expression of cell cycle genes.

Protein Details: Recombinant PRDM11 was expressed in a baculovirus expression system as the full length protein accession number XP_011518530.1) with an N-terminal FLAG tag. The molecular weight of PRDM11 is 112.4 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 this product is shown.

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 PRDM11 protein gel
10% SDS-PAGE gel with Coomassie blue staining

MW: 112.4 kDa

Purity: >90%