

Recombinant PLZF protein

Catalog No: 31545, 31945

Lot No: 31515001

Expressed In: Baculovirus

Quantity: 20 µg

Concentration: 0.35 µg/µl

Source: Human

Buffer Contents: Recombinant full length PLZF protein is supplied at a concentration of 0.35 µg/µl in 25 mM HEPES pH 7.5, 300 mM NaCl, 5% glycerol, 0.04% Triton X-100, 0.2 mM TCEP.

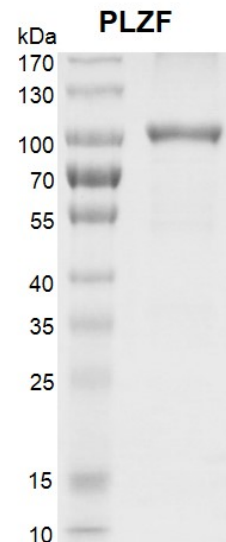
Background: PLZF (Promyelocytic Leukemia Zinc-Finger protein, also known as ZBTB16) is a member of the POK (POZ and Krüppel) family of transcriptional repressors. PLZF is expressed in early, but not differentiated, hematopoietic cells and is required for spermatogonial stem cell maintenance. In addition, PLZF regulates genes involved in cellular proliferation and differentiation (cyclin A2, Myc and Hox genes) and targeted deletion of PLZF in mice results in limb and axial skeleton and limb deformations.

Protein Details: Recombinant human PLZF was expressed in a baculovirus expression system as the full length protein (accession number NP_005997.2) with an N-terminal FLAG tag. The molecular weight of the protein is of 77.3 kDa. The IRF3 protein is >85% pure by SDS-PAGE.

Application Notes: This protein is useful for the study of protein-protein interaction, screening inhibitors, 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 PLZF protein gel.
PLZF protein was run on a 10% SDS-PAGE gel and stained with Coomassie Blue.