

PLZF antibody (mAb)

Catalog Nos: 39987, 39988

RRID: AB_2793419

Clone: 2A9

Isotype: IgG2a

Application(s): ChIP, ICC, IF, IP, WB

Reactivity: Human, Mouse

Quantities: 100 µg, 10 µg

Purification: Protein G Chromatography

Host: Mouse

Concentration: 1 µg/µl

Molecular Weight: 90 kDa

Background: PLZF (Promyelocytic Leukemia Zinc-Finger protein) 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.

Immunogen: This PLZF antibody was raised against a recombinant protein corresponding to full-length mouse PLZF with a GST-Tag.

Buffer: PBS pH 7.5 containing 30% glycerol and 0.035% sodium azide. Sodium azide is highly toxic.

Application Notes:

Storage and Guarantee: Some products may be shipped at room temperature. This will not affect their stability or performance. Avoid repeated freeze/thaw cycles by aliquoting items into single-use fractions for storage at -20°C for up to 2 years. Keep all reagents on ice when not in storage. This product is guaranteed for 12 months from date of receipt.

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