

Aiolos antibody (mAb)

Catalog Nos: 39657, 39658

RRID: AB_2615065

Clone: 8B2

Isotype: IgG

Application(s): ChIP, WB

Reactivity: Mouse

Quantities: 100 µg, 10 µg

Purification: Protein G Chromatography

Host: Mouse

Concentration: 1 µg/µl

Molecular Weight: 70 kDa

Background: Aiolos (also known as IKZF3, Ikaros family zinc finger 3 and ZNFN1A3) belongs to a family of zinc-finger proteins that also include Ikaros and Helios. Aiolos is a transcription factor that is important in the regulation of lymphocyte proliferation and differentiation, which is preferentially expressed in peripheral blood leukocytes, spleen and thymus tissue. Both Ikaros and Aiolos can also participate in chromatin remodeling.

Immunogen: This Aiolos antibody was raised against the N-terminal half of mouse Aiolos protein.

Buffer: Purified IgG in 70 mM Tris (pH 8), 105 mM NaCl, 31 mM glycine, 0.07 mM EDTA, 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.