

## Aiolos antibody (pAb)

**Catalog Nos:** 39293, 39294

**RRID:** AB\_2615037

**Isotype:** Serum

**Application(s):** IP, WB

**Reactivity:** Mouse

**Volumes:** 200 µl, 10 µl

**Purification:** None

**Host:** Rabbit

**Molecular Weight:** 58 kDa

**Background:** Aiolos transcription factor is a member of the Ikaros family of zinc-finger proteins. Aiolos is a hemopoietic-specific transcription factor that regulates B lymphocyte proliferation and differentiation. It can also participate in chromatin remodeling.

**Immunogen:** This Aiolos antibody was raised against a recombinant GST-Tagged protein corresponding to the C-terminal half of mouse Aiolos.

**Buffer:** Rabbit antisera with 30% glycerol containing 0.035% sodium azide. Sodium azide is highly toxic.

**Application Notes:**

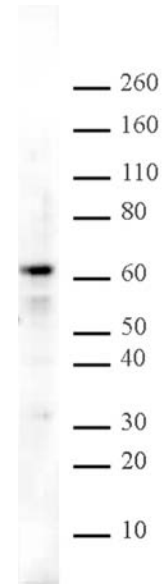
Applications Validated by Active Motif:

WB: 1:500 dilution

The protein migrates around 70 kDa, and multiple isoforms may be detected due to alternative splicing as seen in the example Western blot image.

**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.



**Aiolos pAb tested by Western blot.**

Detection of Aiolos by Western blot analysis. Aiolos was detected using a 1:1,000 dilution of Aiolos pAb with 20 µg mouse spleen tissue extract.