

## Ikaros antibody (mAb)

---

**Catalog Nos:** 39355, 39356

**RRID:** AB\_2614961

**Clone:** 2A9

**Isotype:** IgG

**Application(s):** ChIP, ChIP-Seq, WB

**Reactivity:** Mouse

**Quantities:** 200 µg, 10 µg

**Purification:** Protein G Chromatography

**Host:** Mouse

**Concentration:** 1 µg/µl

**Molecular Weight:** 62 kDa

---

**Background:** Ikaros is a transcription factor that plays a critical role in the control of lymphohematopoiesis and immune regulation. All Ikaros isoforms share a common C-terminal domain that contains a transcription activation motif and two zinc finger motifs required for hetero- and homo-dimerization among the Ikaros isoforms, and for interactions with other proteins. Ikaros is abundantly expressed in pituitary mammosomatotrophs, where it deacetylates Histone H3 sites on the proximal growth hormone (GH) promoter to silence gene expression.

**Immunogen:** This Ikaros antibody was raised against a recombinant protein corresponding to the N-terminal half of mouse Ikaros.

**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:**

Applications Validated by Active Motif:

ChIP: 4 µg per ChIP

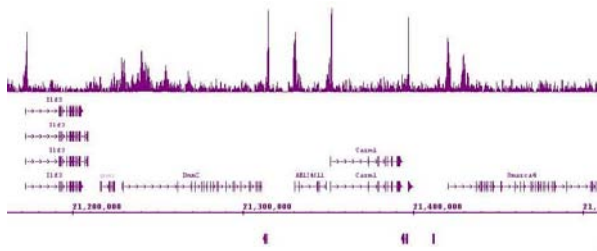
ChIP-Seq: 4 µg each

WB: 0.5 - 2 µg/ml dilution

ChIP-Seq validation was performed by Active Motif's Epigenetics Services; the complete data set is available in the UCSC Genome Browser by clicking [here](#).

**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.



**Ikaros antibody (mAb) (Clone 2A9) tested by ChIP-Seq.**

ChIP was performed using the ChIP-IT<sup>®</sup> High Sensitivity Kit (Cat. No. 53040) with 30 ug of chromatin from mouse spleen and 4 µg of antibody. ChIP DNA was sequenced on the Illumina HiSeq and 15 million sequence tags were mapped to identify Ikaros binding sites. The image shows binding across a region of chromosome 9. You can view the complete data set in the UCSC Genome Browser, starting at this specific location, [here](#).

**Ikaros antibody (mAb) (Clone 2A9) tested by Western blot.**

Detection of Ikaros by Western blot analysis. Ikaros detected in 20 µg mouse spleen whole-cell extract using Ikaros mAb (Clone 2A9) at a 2 µg/ml dilution.

