

Ikaros antibody (pAb)

Catalog Nos: 39291, 39292

RRID: AB_2793219

Isotype: Serum

Application(s): ChIP, EMSA, WB

Reactivity: Mouse

Volumes: 200 µl, 10 µl

Purification: None

Host: Rabbit

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: Rabbit serum containing 30% glycerol and 0.035% sodium azide. Sodium azide is highly toxic.

Application Notes:

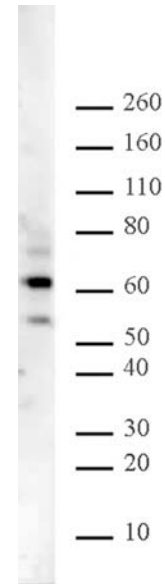
Applications Validated by Active Motif:

WB: 1:500 dilution

Ikaros may be present as multiple protein isoforms due to alternative splicing. Commonly, the two largest isoforms are detected 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.



Ikaros pAb tested by Western blot.

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