

HOXA9 antibody (pAb)

Catalog Nos: 39825, 39826

RRID: AB_2614933

Isotype: IgG

Application(s): WB

Reactivity: Human

Volumes: 100 µl, 10 µl

Purification: Affinity Purified

Host: Rabbit

Molecular Weight: 48 kDa

Background: HOXA9 (Homeobox protein A9, HOX1G) is a sequence-specific DNA binding transcription factor with a central role in both hematopoiesis and leukemia. HOXA9 is required for normal hematopoiesis and contributes to myeloid blood cell differentiation. High levels of HOXA9 expression in hematopoietic cells is a characteristic feature of acute myeloid leukemia (AML), and may contribute to the development of this disease. Overexpression of HOXA9 markedly expands hematopoietic stem cells.

Immunogen: This HOXA9 antibody was raised against a peptide within the N-terminal region of human HOXA9.

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:

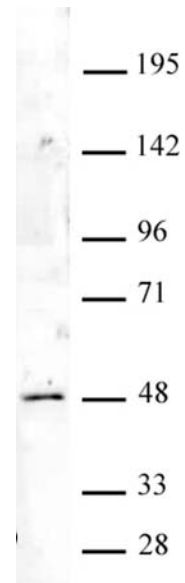
WB*: 1:500 - 1:2,000 dilution

The addition of 0.05% Tween 20 in the blocking buffer and primary antibody incubation buffer is recommended to aid in detection by Western blot. Individual optimization may be required.

*Note: many chromatin-bound proteins are not soluble in a low salt nuclear extract and fractionate to the pellet. Therefore, we recommend a High Salt / Sonication Protocol when preparing nuclear extracts for Western blot.

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.



HOXA9 antibody (pAb) tested by Western blot.

K-562 nuclear extract (20 µg per lane) probed with HOXA9 antibody at a dilution of 1:1,000.