

AML-1/Runx1 antibody (pAb)

Catalog No: 39000

RRID: AB_2793147

Isotype: IgG

Application(s): ChIP, ICC, IF, IHC, IP, WB

Reactivity: Human, Mouse

Volume: 100 µl

Purification: Affinity Purified

Host: Rabbit

Concentration: 0.278 µg/µl

Molecular Weight: 53 kDa

Background: AML-1 (acute myeloid leukemia 1, Runx1, CBFA2, PEBP2αB) is a transcription factor (activator) for various hematopoietic-specific genes (for example, GM-CSF, CSF1R and TCRβ). AML-1 forms a heterodimeric DNA binding complex with its partner protein CBFβ and regulates the expression of cellular genes by binding to promoter and enhancer elements. AML-1 is believed to be involved in the development of normal hematopoiesis and is commonly translocated in hematopoietic cancers.

Immunogen: This AML-1/Runx1 antibody was raised against a synthetic peptide corresponding to amino acid residues 231-245 of human AML-1/Runx1.

Buffer: PBS containing 0.2% gelatin, 0.01% thimersol and 0.1% sodium azide. Thimersol and sodium azide are highly toxic.

Application Notes:

Validated Applications:

IHC (FFPE): 1:100 - 1:500

WB*: 1:500 - 1:1000

Published Applications:

ChIP

IP

WB

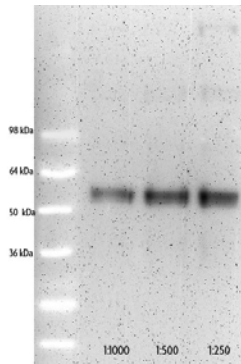
ICC/IF

See references for more information. Individual optimization may be required.

*For optimal results, primary antibody incubations should be performed overnight at 4°C. The addition of 0.1% Tween 20 to all Blotto solutions may also reduce background. Individual optimization may be required.

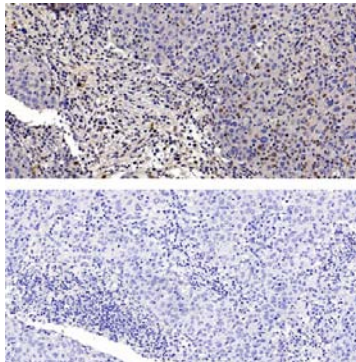
Storage and Guarantee: Some products may be shipped at room temperature. This will not affect their stability or performance. Avoid repeated freeze/thaw cycles by aliquoting items into single-use fractions for storage at -20°C for up to 2 years. Keep all reagents on ice when not in storage. This product is guaranteed for 12 months from date of receipt.

This product is for research use only and is not for use in diagnostic procedures.



AML-1/Runx1 antibody tested by Western blot.

Detection of AML-1/Runx1 by Western blot analysis. Positive detection of AML-1/Runx1 in nuclear extracts derived from Jurkat cells at 1:1,000 (lane 1), 1:500 (lane 2), and 1:250 (lane 3) dilutions.



AML-1 / Runx1 antibody (pAb) tested by Immunohistochemistry

Nuclear staining pattern is detected in Formalin-fixed, paraffin-embedded tissue sections from human breast carcinoma. Top Panel: CTCF antibody at 1:100 dilution. Bottom Panel: No primary antibody (2nd step antibody alone)