

RBBP5 antibody (pAb)

Catalog Nos: 61405, 61406

RRID: AB_2793622

Isotype: IgG

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

Reactivity: Human

Volumes: 100 µl, 10 µl

Purification: Affinity Purified

Host: Rabbit

Molecular Weight: 70 kDa

Background: As part of the MLL1/MLL complex, **RBBP5** (retinoblastoma binding protein 5) is involved in methylation and dimethylation at 'Lys-4' of histone H3. H3 'Lys-4' methylation represents a specific tag for epigenetic transcriptional activation.

Immunogen: This antibody was raised against a peptide within the C-terminal region of human RBBP5.

Buffer: Purified IgG in PBS with 30% glycerol and 0.035% sodium azide. Sodium azide is highly toxic.

Application Notes:

Applications Validated by Active Motif:

ChIP: 10 µl per ChIP

ChIP-Seq: 10 µl each

WB: 1:500 - 1:5,000 dilution

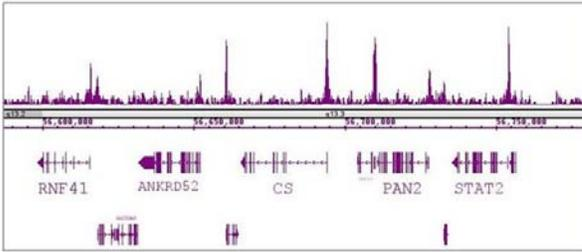
ChIP-Seq validation was performed by Active Motif's Epigenetics Services; the complete data set is available in the UCSC Genome Browser by clicking [here](#).

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.

RBBP5 antibody (pAb) tested by CHIP-Seq.

ChIP was performed using the ChIP-IT[®] High Sensitivity Kit (Cat. No. 53040) with 30 ug of human PBMC chromatin and 7 µl of antibody. ChIP DNA was sequenced on the Illumina HiSeq and 21 million sequence tags were mapped to identify RBBP5 binding sites. The image shows binding across a region of chromosome 12. You can view the complete data set in the UCSC Genome Browser, starting at this specific location, [here](#).



RBBP5 antibody (pAb) tested by Western blot.

MCF-7 whole-cell extract (20 µg) probed with RBBP5 antibody at a dilution of 1:1,000.

