

## ASH2L antibody (pAb)

---

**Catalog Nos:** 39099, 39100

**RRID:** AB\_2615057

**Isotype:** Serum

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

**Reactivity:** Human, Wide Range Predicted

**Volumes:** 100 µl, 10 µl

**Purification:** None

**Host:** Rabbit

**Molecular Weight:** 77 kDa

---

**Background:** ASH2L (absent, small or homeotic discs 2) is a trithorax protein and is a component of different protein complexes, including *i.e.* mixed-lineage leukemia (MLL) oncoproteins, which contain histone methyltransferase activity. These complexes trimethylate the lysine 4 of histone H3. Histone H3 trimethyl Lys4 is one of histone modifications that are part of elaborate mechanisms that evolved to control the structure of the chromatin.

**Immunogen:** This ASH2L antibody was raised against full-length recombinant human ASH2L protein.

**Buffer:** Rabbit serum containing 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 - 1:2,000 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](#).

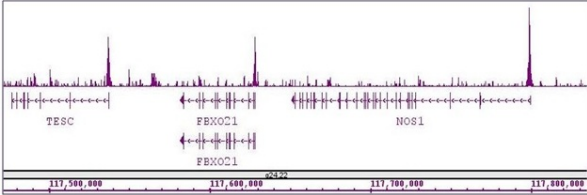
\*Note: many chromatin-bound proteins are not soluble in a low salt nuclear extract and fractionate to the pellet. Therefore, we recommend a High Salt / Sonication Protocol when preparing nuclear extracts for Western blot.

**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.

**ASH2L antibody (pAb) 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 undifferentiated hESC cells and 10 µl of antibody. ChIP DNA was sequenced on the Illumina HiSeq and 20 million sequence tags were mapped to identify ASH2L 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](#).



**ASH2L pAb tested by Western blot.**

ASH2L detection by Western blot. The analysis of ASH2L was performed using HeLa nuclear extract and ASH2L pAb at a 1:2,000 dilution.

