

FOXG1 antibody (pAb)

Catalog Nos: 61211, 61212

RRID: AB_2793554

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

Reactivity: Human, Mouse

Volumes: 100 µl, 10 µl

Purification: Affinity Purified

Host: Rabbit

Isotype: IgG

Molecular Weight: 55 kDa

Background: FOXG1 (Forkhead box protein G1, FKH2, FKHL1) is a transcriptional repressor that serves an important role in the development of the brain. FOXG1 is expressed in the neural plate early in development and is involved in the specification, proliferation and differentiation of the telencephalon. FOXG1 is also essential for proper patterning of the eye and development of the olfactory epithelium.

Immunogen: This FOXG1 antibody was raised against a peptide within the C-terminal region of human FOXG1.

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: 5 µl per ChIP

ChIP-Seq: 5 µl each

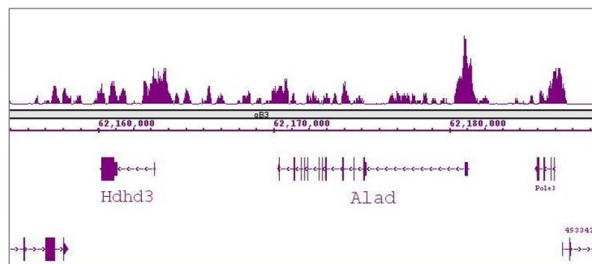
WB: 1:500 - 1:1,000 dilution

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](#).

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

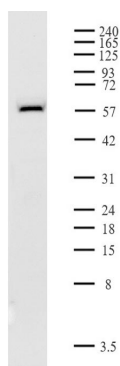
Storage and Guarantee: Some products may be shipped at room temperature. This will not affect their stability or performance. Avoid repeated freeze/thaw cycles by aliquoting items into single-use fractions for storage at -20°C for up to 2 years. Keep all reagents on ice when not in storage. This product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

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



FOXG1 antibody (pAb) tested by ChIP-Seq.

ChIP was performed using the ChIP-IT® High Sensitivity Kit (Cat. No. 53040) with 30 ug of mouse whole brain chromatin and 2 µl of antibody. ChIP DNA was sequenced on the Illumina HiSeq and 20 million sequence tags were mapped to identify FOXG1 binding sites. The image shows binding across a region of chromosome 4. You can view the complete data set in the UCSC Genome Browser, starting at this specific location, [here](#).



FOXG1 antibody (pAb) tested by Western blot.

Detection of Goosecoid by Western blot analysis. Cytosolic extract of Jurkat cells (40 µg). Both were probed with FOXG1 antibody (pAb) at a 1:500 dilution.