

HNF-3 α / FOXA1 antibody (mAb)

Catalog Nos: 39837, 39838

RRID: AB_2793362

Clone: 2F83

Isotype: IgG

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

Reactivity: Human, Mouse

Quantities: 100 μ g, 10 μ g

Purification: Protein G Chromatography

Host: Mouse

Concentration: 1 μ g/ μ l

Molecular Weight: 50 kDa

Background: HNF-3 α (Hepatocyte nuclear factor 3 alpha, forkhead box A1, FoxA1) is a homeodomain containing transcription factor and is a member of the Hepatocyte Nuclear Factor family (HNF-1 α , HNF-1 β , HNF-3 α , HNF-3 β , HNF-3 γ , HNF-4 α , HNF-4 γ , HNF-6 α and HNF-6 β), whose members regulate the transcription of numerous hepatocyte genes in adult liver. HNF-3 members bind to DNA as monomers using a winged helix DNA-binding domain. HNF-3 α and HNF-3 γ have also been shown to be involved in gastrulation events such as body axis formation.

Immunogen: This HNF-3 α / FOXA1 antibody was raised against a recombinant protein corresponding to amino acids 7-103 of rat HNF-3 α /FOXA1.

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:

ChIP: 5 - 10 μ g per ChIP

ChIP-Seq: 5 - 10 μ g each

WB: 1 - 5 μ g/ml dilution

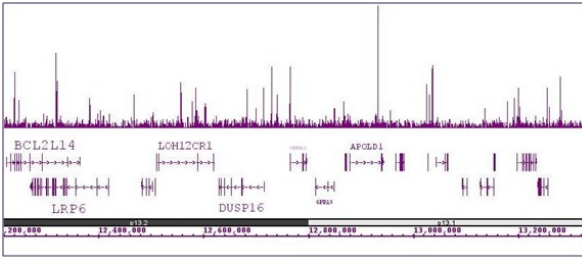
ChIP-Seq validation was performed by Active Motif's Epigenetics Services; the complete data set is available in the UCSC Genome Browser by clicking [here](#).

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.

HNF-3 α / FOXA1 antibody (mAb) tested by ChIP-Seq.

ChIP was performed using the ChIP-IT[®] High Sensitivity Kit (Cat. No. 53040) with 30 μ g of chromatin from LnCaP cells and 4 μ g of antibody. ChIP DNA was sequenced on the Illumina HiSeq and 20 million sequence tags were mapped to identify HNF-3 α / FOXA1 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](#).



HNF-3 α / FOXA1 antibody (mAb) tested by Western blot.

Extract of HeLa cells over-expressing mouse HNF-3 α / FOXA1 (50 μ g per lane) probed with HNF-3 β / FOXA1 antibody (mAb) (1:5,000 dilution).

