

## HNF-3 $\beta$ / FOXA2 antibody (pAb)

**Catalog Nos:** 39827, 39828

**RRID:** AB\_2793357

**Isotype:** Serum

**Application(s):** IHC, WB

**Reactivity:** Human, Rat

**Volumes:** 100  $\mu$ l, 10  $\mu$ l

**Purification:** None

**Host:** Rabbit

**Molecular Weight:** 50 kDa

**Background:** HNF-3 $\beta$  (Hepatocyte nuclear factor 3 beta, forkhead box A2, FoxA2) is a homeodomain containing transcription factor and is a member of the Hepatocyte Nuclear Factor family (HNF-1 $\alpha$ , HNF-1 $\beta$ , HNF-3 $\alpha$ , HNF-3 $\beta$ , HNF-3 $\gamma$ , HNF-4 $\alpha$ , HNF-4 $\gamma$ , HNF-6 $\alpha$  and HNF-6 $\beta$ ), 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 $\alpha$  and HNF-3 $\gamma$  have also been shown to be involved in gastrulation events such as body axis formation.

**Immunogen:** This HNF-3 $\beta$  / FOXA2 antibody was raised against a recombinant protein corresponding to amino acids 7-86 of rat HNF-3 $\beta$  / FOXA2.

**Buffer:** Rabbit serum containing 30% glycerol and 0.035% sodium azide. Sodium azide is highly toxic.

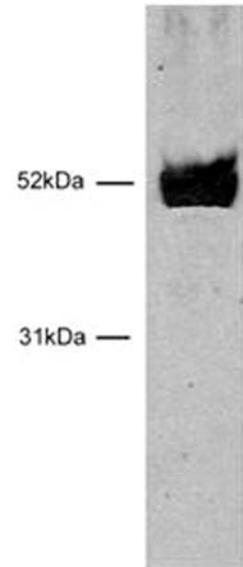
### Application Notes:

Applications Validated by Active Motif:

WB: 1:2,000 - 1:5,000 dilution

**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 $\beta$ / FOXA2 antibody (pAb) tested by Western blot.

Extract of HeLa cells expressing HNF-3 $\beta$  / FOXA2 (50  $\mu$ g per lane) probed with HNF-3 $\beta$  / FOXA2 antibody (pAb) (1:5,000 dilution).