

## FOXO1/FKHR antibody (pAb)

**Catalog Nos:** 39629, 39630

**RRID:** AB\_2614952

**Isotype:** IgG

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

**Reactivity:** Human

**Volumes:** 200  $\mu$ l, 10  $\mu$ l

**Purification:** Affinity Purified

**Host:** Rabbit

**Molecular Weight:** 90 kDa

**Background:** FOXO1/FKHR (forkhead box 01) belongs to the forkhead family of transcription factors that contain a distinct forkhead domain. Chromosomal aberrations involving FOXO1, such as translocation with PAX3, are associated with avascular rhabdomyosarcoma. The specific function of this gene is undetermined.

**Immunogen:** This FOXO1/FKHR antibody was raised against a peptide from the C-terminal region of human FOXO1/FKHR.

**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:

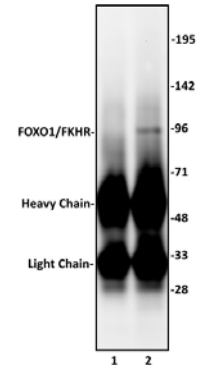
Applications Validated by Active Motif:

IP: 10  $\mu$ l per IP

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

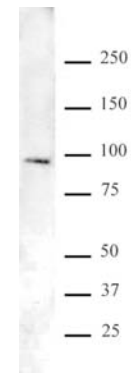
**Storage and Guarantee:** Some products may be shipped at room temperature. This will not affect their stability or performance. Avoid repeated freeze/thaw cycles by aliquoting items into single-use fractions for storage at  $-20^{\circ}\text{C}$  for up to 2 years. Keep all reagents on ice when not 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.



### FOXO1/FKHR antibody (pAb) tested by Immunoprecipitation.

10  $\mu$ l of FOXO1/FKHR antibody was used to immunoprecipitate FOXO1/FKHR from 500  $\mu$ g of K562 nuclear cell extract (lane 2). 10  $\mu$ l of rabbit IgG was used as a negative control (lane 1). The immunoprecipitated protein was detected by Western blotting using the FOXO1/FKHR antibody at a dilution of 1:500.



### FOXO1/FKHR antibody (pAb) tested by Western blot.

U2OS nuclear extract (20  $\mu$ g per lane) was probed with 39629 at a 1:500 dilution.