

## FUBP1 polyclonal antibody - Classic

**Cat. No.** C15410233

**Type:** Polyclonal

**Source:** Rabbit

**Lot #:** 40912

**Size:** 25 µl/100 µl

**Concentration:** 1 µg/µl

**Specificity:** Human.

**Purity:** Purified by peptide affinity chromatography.

**Storage:** Store at -20°C; for long storage, store at -80°C. Avoid multiple freeze-thaw cycles.

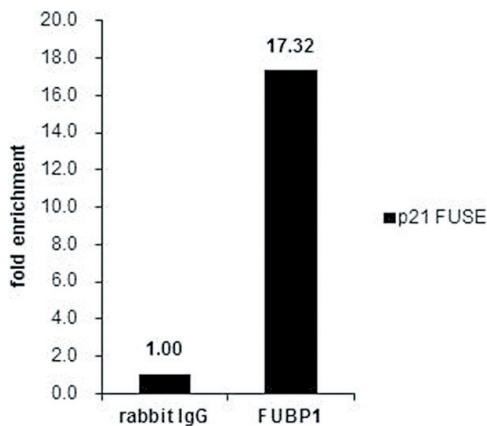
**Precautions:** This product is for research use only. Not for use in diagnostic or therapeutic procedures.

### Applications

	Suggested dilution*	Results
ChIP assay	5 µg/ChIP	Figure 1
ICC/IF	1:100-1:1,000	Figure 5
IHC	1:100-1:1,000	Figure 3
Immunoprecipitation	1:100-1:500	Figure 2
WB	1:100-1:1,000	Figure 4

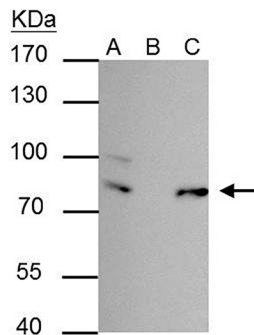
\*Optimal dilutions/concentrations should be determined by the researcher.

### Results



**Figure 1. FUBP1 ChIP results**

ChIP was performed with HeLa chromatin extract and 5 µg of either control rabbit IgG or FUBP1 antibody. The precipitated DNA was detected by PCR with primer set targeting to p21 FUSE.

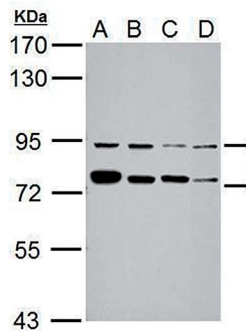
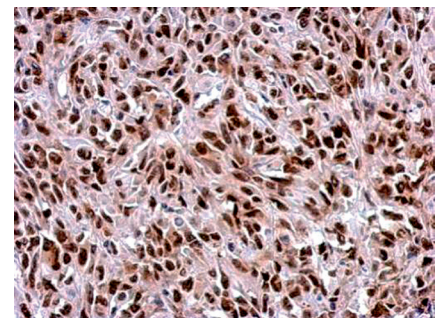


**Figure 2. FUBP1 IP results**

FUBP1 antibody immunoprecipitates FUBP1 protein in IP experiments. IP Sample: 293T whole cell extract A: 30 µg whole cell extract of FUBP1 protein expressing 293T cells B: Control with 2.5 µg of pre-immune rabbit IgG C : Immunoprecipitation of FUBP1 by 2.5 µg of FUBP1 antibody. The immunoprecipitated FUBP1 protein was detected by western blot with the FUBP1 antibody diluted 1:1,000.

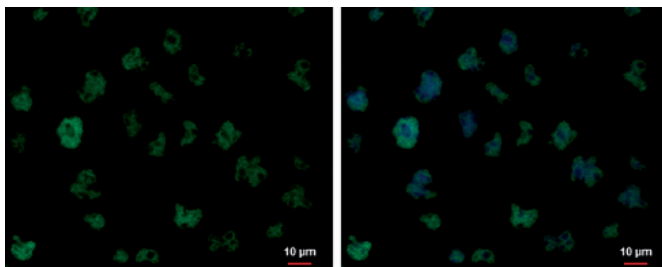
**Figure 3. IHC**

Immunohistochemical analysis of paraffin-embedded HeLa xenograft, using FUBP1 antibody at a 1:500 dilution.



**Figure 4. Western blot**

Sample: 30 µg of whole cell lysate  
 A: Jurkat  
 B: K562  
 C: THP-1  
 D: NCI-H929  
 7.5% SDS PAGE  
 FUBP1 antibody diluted 1:5,000



**Figure 5. FUBP1 IF results**

FUBP1 antibody detects FUBP1 protein at nucleus by immunofluorescent analysis.  
 Sample: Jurkat cells were fixed in 4% paraformaldehyde at RT for 15 min.  
 Green: FUBP1 protein stained by FUBP1 antibody diluted 1:500.  
 Blue: Hoechst 33342 staining.