

FUBP1 Antibody - ChIP Grade

Cat. No. C15410233-100

Type: Polyclonal	Specificity: Human
Size: 100 µl	Isotype: NA
Concentration: 1 µg/µl	Host: Rabbit
Lot No.: 40912	Purity: Affinity purified
Storage buffer: NA	Storage conditions: NA
Precautions: This product is for research use only. Not for use in diagnostic or therapeutic procedures.	

Description

Polyclonal antibody raised in rabbit against **FUBP1** (far upstream element-binding protein 1), using a recombinant protein.

Applications

Applications	Suggested dilution *	References
ChIP assay	5 µg/ChIP	Fig 1
ICC/IF	1:100 - 1:1,000	Fig 5
IHC	1:100 - 1:1,000	Fig 3
Immunoprecipitation	1:100 - 1:500	Fig 2
Western Blotting	1:100 - 1:1,000	Fig 4

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

Validation Data

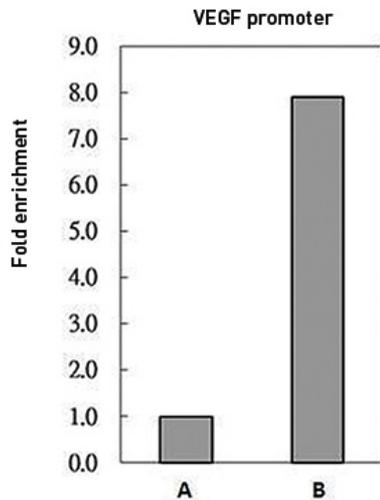


Figure 1. FUBP1 ChIP results

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

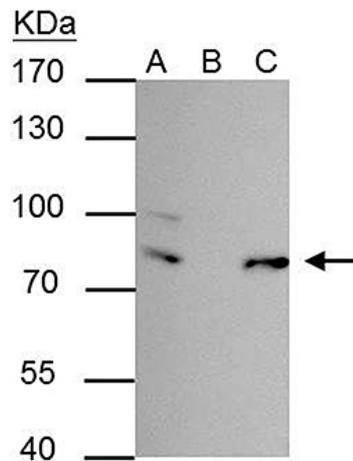


Figure 2. FUBP1 IP results

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

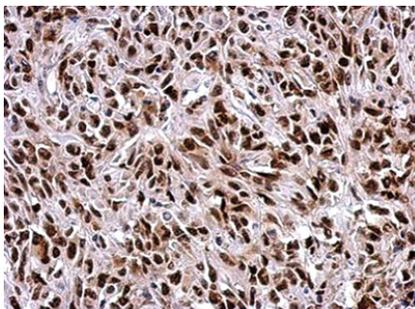


Figure 3. IHC

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

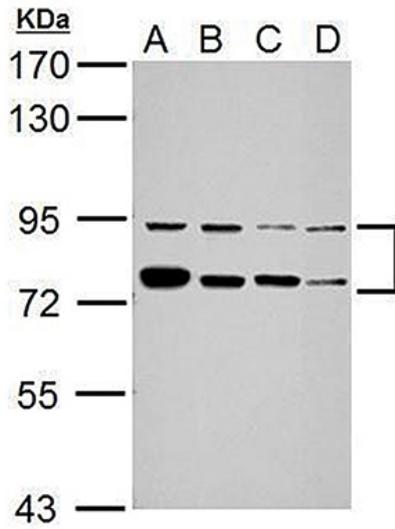


Figure 4. Western blot

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

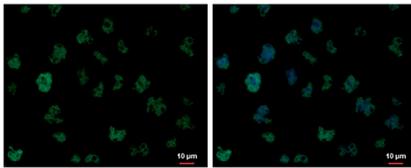


Figure 5. FUBP1 IF results

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.