

## PKNOX1 polyclonal antibody - Classic

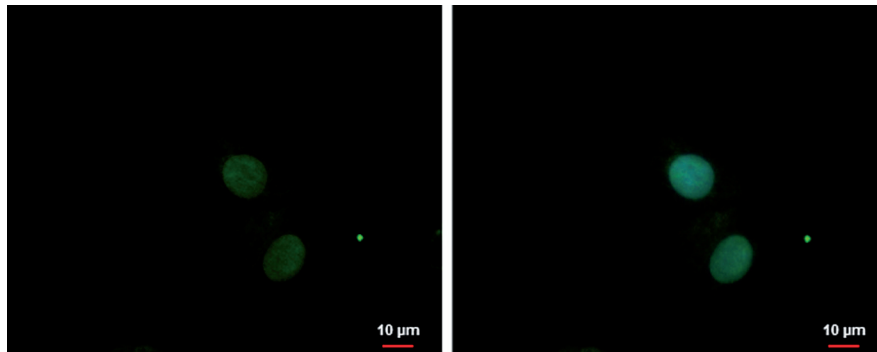
**Cat. No.** C15410267**Type:** Polyclonal**Source:** Rabbit**Lot #:** 40870**Size:** 25 µl/100 µl**Concentration:** 1 µg/µl**Specificity:** Human**Purity:** Purified by peptide affinity chromatography**Storage:** Store at -20°C; for long storage, store at -80°C.  
Avoid multiple freeze-thaw cycles.**Precautions:** This product is for research use only. Not for use in diagnostic or therapeutic procedures.

### Applications

	Suggested dilution*	Results
IF	1:100-1:1,000	Figure 1
Immunoprecipitation	1:100-1:1,000	Figure 2
Western blot	1:500-1:3,000	Figure 3

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

### Results



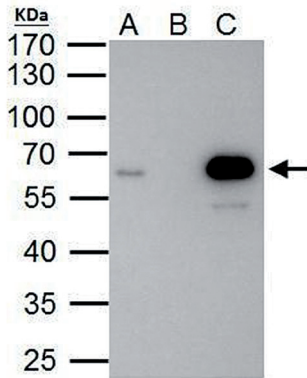
**Figure 1. PKNOX1 IF results**

PKNOX1 antibody detects PKNOX1 protein at nucleus by immunofluorescent analysis.

Sample: HepG2 cells were fixed in 4% paraformaldehyde at RT for 15 min.

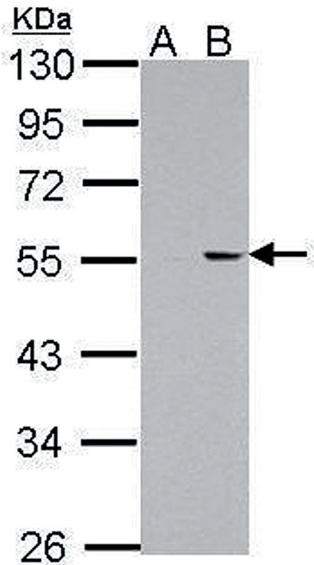
Green: PKNOX1 protein stained by PKNOX1 antibody diluted 1:1,000.

Blue: Hoechst 33342 staining.



**Figure 2. PKNOX1 Immunoprecipitation results**

PKNOX1 antibody immunoprecipitates PKNOX1 protein in IP experiments. IP Sample: HeLa whole cell extract A. 40 µg HeLa whole cell extract B. Control with 2 µg of preimmune rabbit IgG C. Immunoprecipitation of PKNOX1 protein by 2 µg of PKNOX1 antibody The immunoprecipitated PKNOX1 protein was detected by western blot with the PKNOX1 antibody diluted 1:1,000.



**Figure 3. PKNOX1 Western blot results**

Sample: 30 µg of whole cell lysate  
A: HepG2  
B: HepG2 nucleus  
10% SDS PAGE  
PKNOX1 antibody diluted 1:2,000