

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

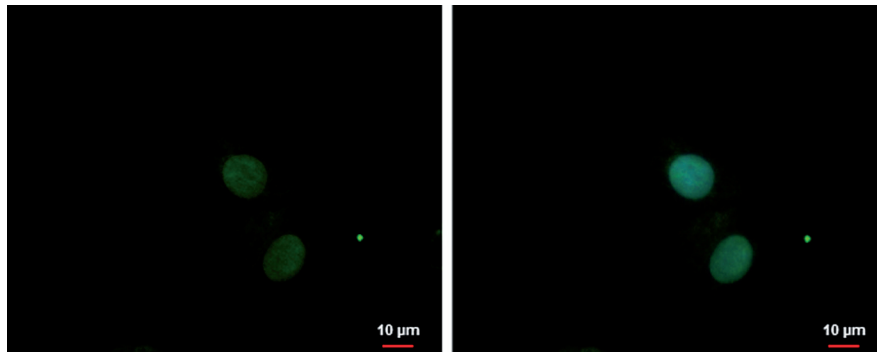


Figure 1. PKNOX1 IF results

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.

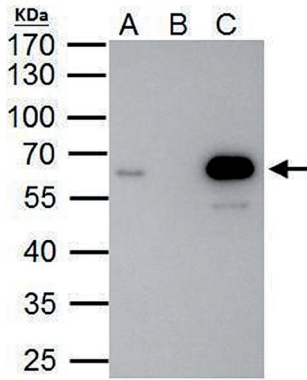


Figure 2. PKNOX1 Immunoprecipitation results

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.

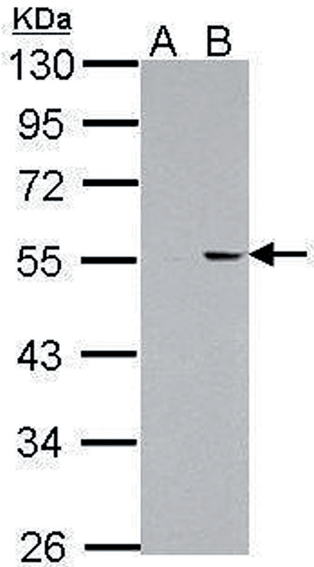


Figure 3. PKNOX1 Western blot results

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