

THOC1 polyclonal antibody - Classic

Cat. No. C15410243

Type: Polyclonal

Source: Rabbit

Lot #: 40163

Size: 25 µl/100 µl

Concentration: 1 µg/µl

Specificity: Human, mouse

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 3
IHC	1:100-1:1,000	Figure 4
Immunoprecipitation	1:100-1:500	Figure 2
Western blot	1:500-1:3,000	Figure 5

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

Results

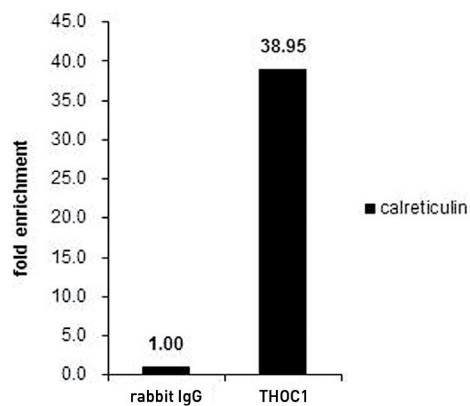


Figure 1. THOC1 ChIP results

ChIP was performed with HeLa chromatin extract and 5 µg of either control rabbit IgG or THOC1 antibody. The precipitated DNA was detected by PCR with primer set targeting to calreticulin.

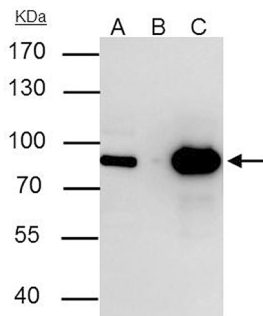


Figure 2. THOC1 IP results

THOC1 immunoprecipitates THOC1 (p84) protein in IP experiments. IP Sample: HepG2 whole cell extract A: 30 µg whole cell extract of THOC1 protein expressing HepG2 cells B: Control with 3 µg of pre-immune rabbit IgG C: Immunoprecipitation of THOC1 by 3 µg of THOC1 antibody. The immunoprecipitated THOC1 protein was detected by western blot with the THOC1 antibody diluted 1:1,000.

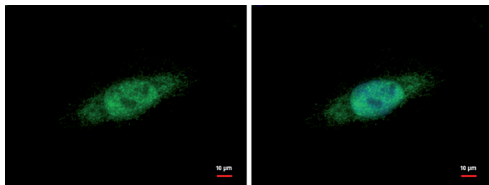


Figure 3. THOC1 IF results

THOC1 antibody detects THOC1 protein at cytoplasm and nucleus by immunofluorescent analysis. Sample: HeLa cells were fixed in 4% paraformaldehyde at RT for 15 min. Green: THOC1 protein stained by THOC1 antibody diluted 1:500. Blue: Hoechst 33342 staining.

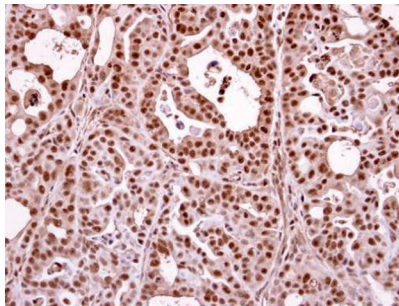


Figure 4. THOC1 IHC results

Immunohistochemical analysis of paraffin-embedded NCIN87 xenograft, using THOC1 antibody at a 1:500 dilution.

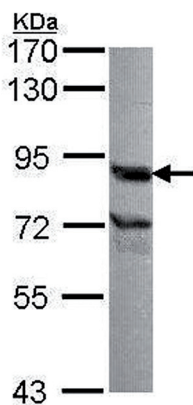


Figure 5. THOC1 Western blot results

Sample: 30 µg of A549 whole cell lysate
7.5% SDS PAGE
THOC1 antibody diluted 1:1,000