

TRRAP polyclonal antibody - Classic

Cat. No. C15410257

Type: Polyclonal

Source: Rabbit

Lot #: 41591

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

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

Results

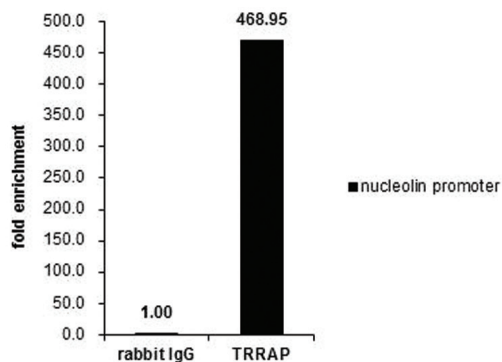


Figure 1. TRRAP ChIP results

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

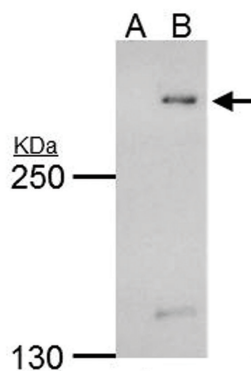


Figure 2. TRRAP IP results

TRRAP antibody immunoprecipitates TRRAP protein in IP experiments. IP samples: 293T whole cell extract
 A. Control with 4 µg of preimmune Rabbit IgG
 B. Immunoprecipitation of TRRAP protein by 4 µg TRRAP antibody (Cat. No. C15410257)
 5 % SDS-PAGE
 The immunoprecipitated TRRAP protein was detected by TRRAP antibody (C15410257) diluted 1:500.

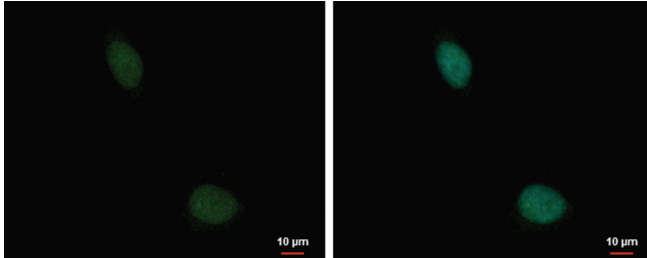


Figure 3. TRRAP IF results

TRRAP antibody detects TRRAP protein at nucleus by immunofluorescent analysis.

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

Green: TRRAP protein stained by TRRAP antibody [Cat. No. C15410257] diluted 1:1,000.

Blue: Hoechst 33342 staining.

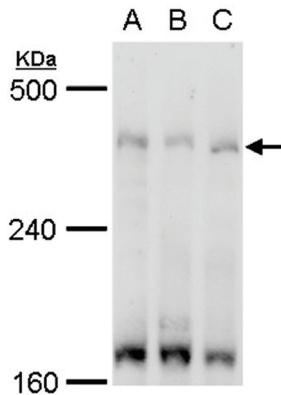


Figure 4. TRRAP Western blot results

TRRAP antibody detects TRRAP protein by western blot analysis.

A. 30 μ g Jurkat whole cell extract

B. 30 μ g Raji whole cell extract

C. 30 μ g NCI-H929 whole cell extract

5 % SDS-PAGE

TRRAP antibody [Cat. No. C15410257] dilution: 1:500

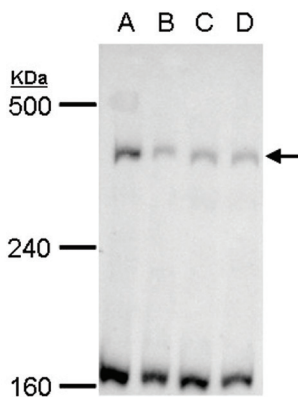


Figure 5. TRRAP western blot results

TRRAP antibody detects TRRAP protein by western blot analysis.

A. 30 μ g 293T whole cell extract

B. 30 μ g A431 whole cell extract

C. 30 μ g HeLa whole cell extract

D. 30 μ g HepG2 whole cell extract

5 % SDS-PAGE

TRRAP antibody [Cat. No. C15410257] dilution: 1:500