

SMAD3 polyclonal antibody - Classic

Cat. No. C15410252

Type: Polyclonal

Source: Rabbit

Lot #: 40044

Size: 25 µl/100 µl

Concentration: 0.78 µg/µl

Specificity: Human.

Purity: Purified by antigen-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	1:500	Figure 5
Immunoprecipitation	1:500-1:1,000	Figure 2
Western blot	1:500-1:3,000	Figure 4

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

Results

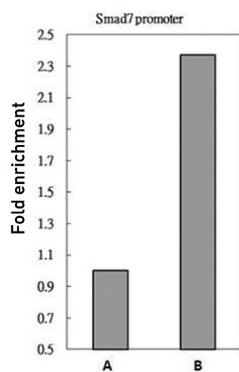


Figure 1. ChIP

SMAD3 antibody immunoprecipitates SMAD3 protein-DNA in ChIP experiments. ChIP Sample: HeLa chromatin extract A. 5 µg preimmune rabbit IgG B. 5 µg of SMAD3 antibody (C15410252). The precipitated DNA was detected by PCR with primer set targeting to SMAD7 promoter.

Figure 2. Immunoprecipitation

SMAD3 antibody immunoprecipitates SMAD3 protein in IP experiments. IP Sample: 1000 µg HeLa whole cell extract A. 50 µg HeLa whole cell extract B. Control with 2 µg of preimmune rabbit IgG C. Immunoprecipitation of SMAD3 protein by 2 µg of SMAD3 antibody (C15410252). The immunoprecipitated SMAD3 protein was detected by western blot with the SMAD3 antibody (C15410252) diluted 1:1,000.

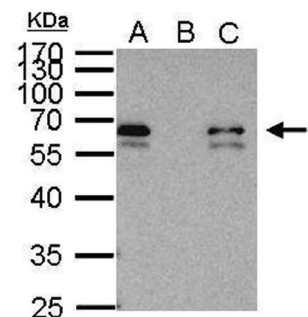




Figure 3. Immunofluorescence

Immunofluorescence analysis of paraformaldehyde-fixed HeLa cells using SMAD3 (C15410252) antibody at a 1:200 dilution.

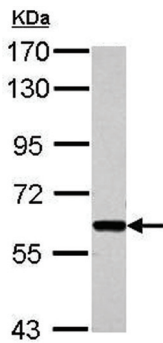


Figure 4. Western Blot

Sample: 30 µg of Jurkat whole cell lysate
7.5% SDS PAGE
C15410252 diluted 1:5,000

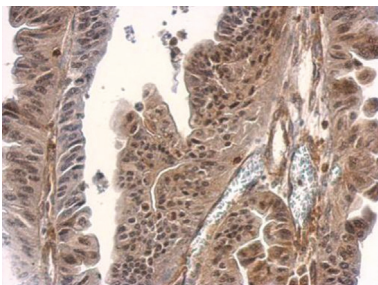


Figure 5. Immunohistochemistry

SMAD3 antibody detects SMAD3 protein at cytosol and nucleus on gastric carcinoma by immunohistochemical analysis.

Sample: Paraffin-embedded gastric carcinoma.
SMAD3 antibody (C15410252) dilution: 1:500.