

HIF1 alpha polyclonal antibody - Classic

Cat. No. C15410234

Type: Polyclonal

Source: Rabbit

Lot #: 41185

Size: 25 µg/100 µg

Concentration: 1 µg/µl

Specificity: Human, mouse

Purity: Purified by and 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 7
IHC	1:100-1:1,000	Figure 3
Immunoprecipitation	1:100-1:500	Figure 2
WB	1:1,000-1:10,000	Figure 4, 5, 6

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

Results

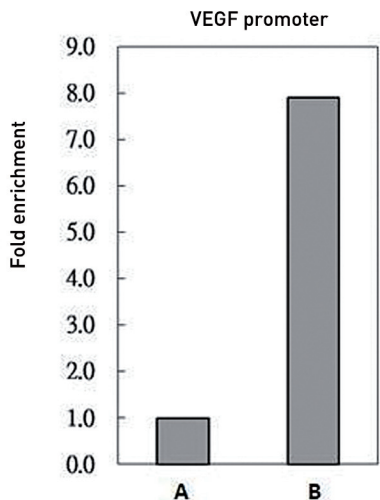


Figure 1. HIF1 alpha ChIP results

HIF1 alpha antibody immunoprecipitates HIF1 alpha protein-DNA complex in ChIP experiments. ChIP Sample: HepG2 whole chromatin extract under hypoxia (1%O₂) treatment 24hr A: 5 µg preimmune rabbit IgG. B: 5 µg of HIF1 alpha antibody The precipitated DNA was detected by PCR with primer set targeting to VEGF promoter.

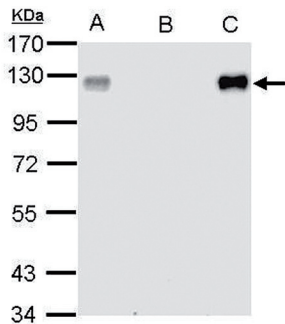


Figure 2. HIF1 alpha IP results

HIF1 alpha antibody immunoprecipitates HIF1 alpha protein in IP experiments. IP Sample: 950 µg whole cell extract of HIF1 alpha protein expressing HeLa cells. A. 20 µg whole cell extract of transfected HeLa cells B. Control with 2.5 µg of preimmune rabbit IgG C. Immunoprecipitation of HIF1 alpha protein by 2.5 µg of HIF1 alpha antibody. The immunoprecipitated HIF1 alpha protein was detected by western blot with the HIF1 alpha antibody diluted 1:10,000.

Figure 3. HIF1 alpha IHC results

HIF1 alpha antibody detects HIF1 alpha protein at nucleus on human kidney cancer by immunohistochemical analysis. Sample: Paraffin-embedded human kidney cancer. HIF1 alpha antibody dilution: 1:500.

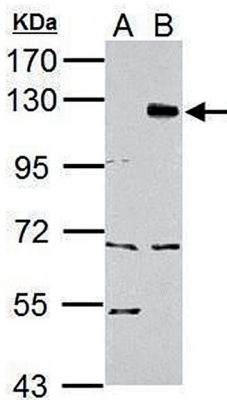
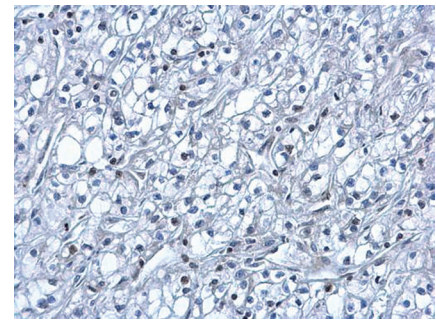


Figure 4. HIF1 alpha Western blot results

HIF1 alpha antibody detects HIF1 alpha protein in CoCl₂-treated samples by Western blot analysis.

A: 30 µg of HeLa cell lysate (normoxia)

B: 30 µg of HeLa cell lysate (500 µM CoCl₂ treatment for 24 hrs)

7.5% SDS-PAGE

HIF1 alpha antibody dilution: 1:1,000

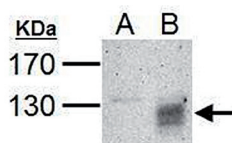


Figure 5. HIF1 alpha Western blot results

HIF1 alpha antibody detects HIF1 alpha protein in CoCl₂-treated samples by Western blot analysis.

A. 30 µg NIH-3T3 whole cell extract (untreated)

B. 30 µg NIH-3T3 whole cell extract (200 µM CoCl₂ treatment for 24hr)

HIF1 alpha antibody dilution: 1:1,000

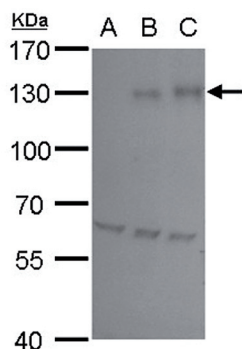


Figure 6. HIF1 alpha Western blot results

HIF1 alpha antibody detects HIF1 alpha protein by western blot analysis.

A. 30 ug HepG2 whole cell extract (untreated)

B. 30 μg HepG2 whole cell extract (200 μM CoCl₂ treatment for 24 hr)

C. 30 μg HepG2 whole cell extract (500 μM CoCl₂ treatment for 24 hr)

7.5 % SDS-PAGE

HIF1 alpha antibody dilution: 1:5000

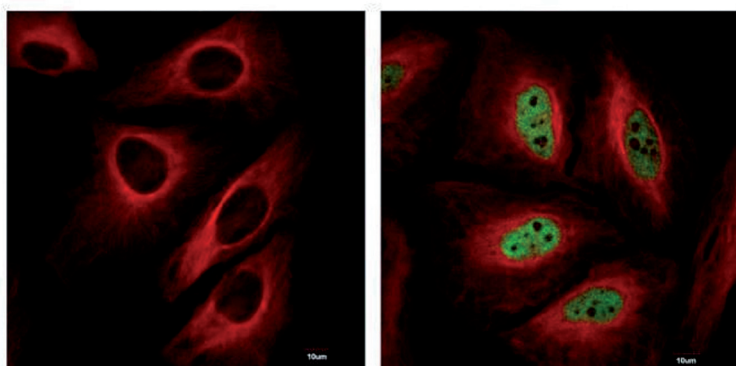


Figure 7. HIF1 alpha IF results

HIF1 alpha antibody detects HIF1 alpha protein at nucleus in CoCl₂-treated samples by immunofluorescent analysis.

Samples: Untreated HeLa cells (left panel) and HeLa cells treated with 500 μM CoCl₂ for 24 hrs (right panel) were fixed in 4% paraformaldehyde for 15 min.

Green: HIF1alpha protein stained by HIF1 alpha antibody diluted 1:1,000

Red: Alpha-tubulin, a cytoskeleton marker.