

NRF2 polyclonal antibody - Classic

Cat. No. C15410242-25

Type: Polyclonal	Specificity: Human, rat
Size: 25 µl	Isotype: NA
Concentration: 0.58 µg/µl	Source: Rabbit
Lot No.: 42347	Purity: Affinity purified polyclonal antibody.
Storage buffer: PBS containing 20% glycerol and 0.025% ProClin 300.	Storage conditions: 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.	

Description

Polyclonal antibody raised in rabbit against NRF2 (nuclear factor (erythroid-derived 2)-like 2), using a recombinant protein.

Applications

Applications	Suggested dilution	References
ChIP *	5 µg/ChIP	Fig 1
Immunofluorescence	1:500	Fig 2
Immunohistochemistry	1:500	Fig 3
Western Blotting	1:1,000	Fig 4, 5
Immunoprecipitation	5 µg/IP	Fig 6

* Please note that the optimal antibody amount per ChIP should be determined by the end-user. We recommend testing 1-5 µg per IP.

Target Description

NRF2 (UniProt/Swiss-Prot entry Q16236) is a transcriptional factor which belongs to the family of basic leucine zipper proteins. It recognizes the antioxidant response (ARE) elements in the promoter regions of its target genes. As such, it regulates the response to injury, oxidative stress and inflammation, including the production of free radicals. NRF2 may also be involved in the transcriptional activation of genes of the beta-globin cluster.

Validation Data

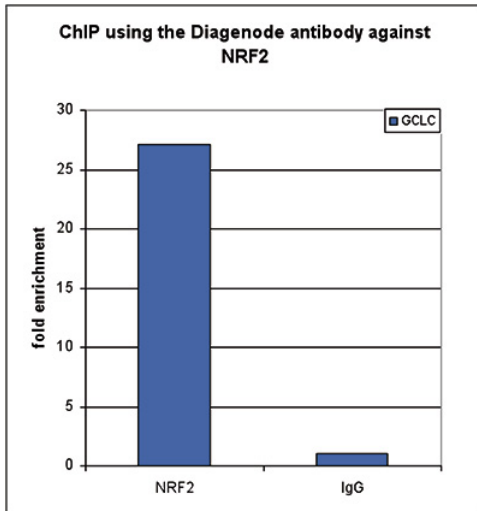


Figure 1. ChIP results obtained with the Diagenode antibody directed against NRF2

ChIP was performed on HeLa cells with 5 µg of the Diagenode antibody against NRF2 (Cat. No. C15410242). IgG (5 µg/IP) was used as negative IP control. QPCR was performed with primers for the GCLC gene. Figure 1 shows the fold enrichment over the IgG negative control.

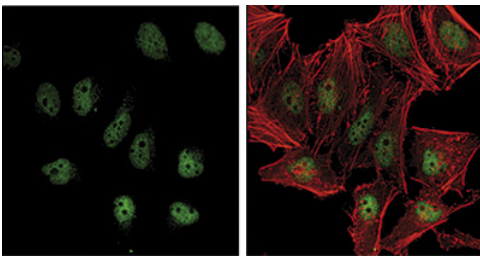


Figure 2. Immunofluorescence using the Diagenode antibody directed against NRF2

HeLa cells were stained with the Diagenode antibody against NRF2 (Cat. No. C15410242). Cells were fixed with 4% formaldehyde for 15 min at RT and blocked with PBS/TX-100 containing 5% normal goat serum and 1% BSA. The cells were immunofluorescently labeled with the NRF2 antibody (left) diluted 1:500 in blocking solution. The right panel shows costaining with Phalloidin, a cytoskeleton marker.

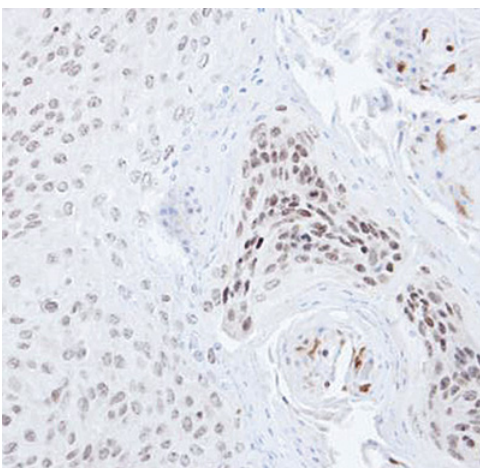


Figure 3. Immunohistochemistry using the Diagenode antibody directed against NRF2

Paraffin-embedded Cal27 xenograft was analysed by immunohistochemical analysis, using the Diagenode antibody against NRF2 (Cat. No. C15410242) diluted 1:500.

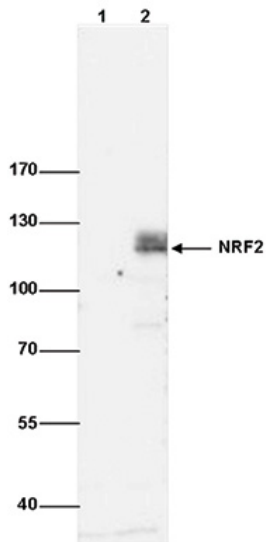


Figure 4. Western blot analysis using the Diagenode antibody directed against NRF2

Whole cell extracts from 293T cells (30 μ g) transfected with a 3xFlag-tagged NRF2 expression vector (lane 2) and untransfected control cells (lane 1) were analysed by western blot using the Diagenode antibody against NRF2 (Cat. No. C15410242) diluted 1:1,000. The position of the protein of interest is indicated on the right; the marker (in kDa) is shown on the left.

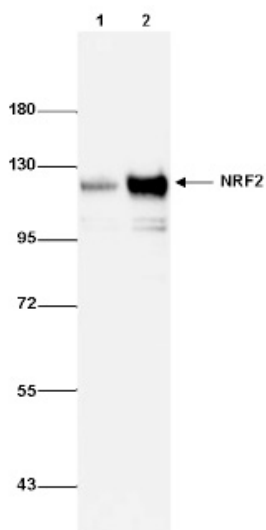


Figure 5. Western blot analysis using the Diagenode antibody directed against NRF2

Nuclear extracts from MDA-MB-231 cells (30 μ g) treated with 30 μ M tBHQ for 30' (lane 2) or untreated control cells (lane 1) were analysed by western blot using the Diagenode antibody against NRF2 (Cat. No. C15410242) diluted 1:500. The position of the protein of interest is indicated on the right; the marker (in kDa) is shown on the left.

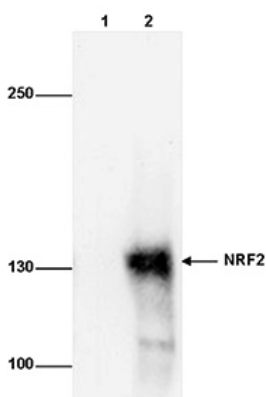


Figure 6. Immunoprecipitation using the Diagenode antibody directed against NRF2

Immunoprecipitation was performed on whole cell extracts from HepG2 cells using 5 μ g of the Diagenode antibody against NRF2 (Cat. No. C15410242) (lane 2) or with an equal amount of rabbit IgG (lane 1). The immunoprecipitated NRF22 protein was detected by western blot with the NRF2 antibody diluted 1:1,000.