

## NRF2 polyclonal antibody - Classic

**Other names:** NFE2L2, HEBP1, TFIIIE-B, FE

**Cat. No.** C15410242

**Type:** Polyclonal **ChIP-grade**

**Source:** Rabbit

**Lot #:** 42004

**Size:** 25 µl/100 µl

**Concentration:** 0.51 µg/µl

**Specificity:** Human, rat: positive  
Other species: not tested

**Purity:** Affinity purified polyclonal antibody in PBS containing 20% glycerol and 0.025% ProClin 300

**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.

**Description:** Polyclonal antibody raised in rabbit against NRF2 (NF-E2-Related Factor 2), using a recombinant protein.

### Applications

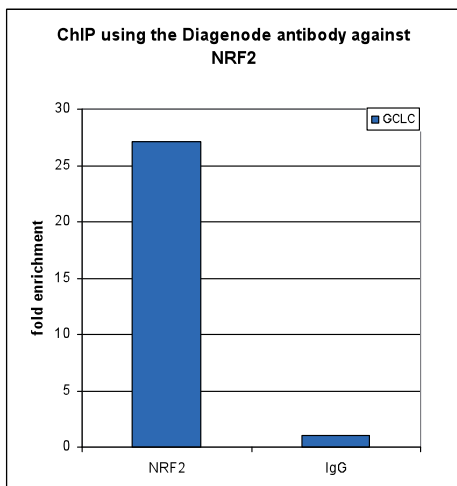
	Suggested dilution*	Results
ChIP*	5 µg per ChIP	Figure 1
IF	1:500	Figure 2
IHC	1:500	Figure 3
Western blot	1:500 - 1:1,000	Figure 4, 5
Immunoprecipitation	5 µg per IP	Figure 6

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

### Target description

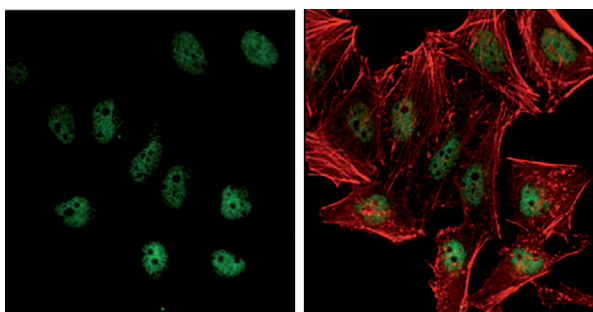
NRF2 (UniProt/Swiss-Prot entry Q16236) is a transcriptional factor which belongs to the family of basic leucine zipper proteins. It recognizes the antioxidant response (ARE) elements in the promoter regions of its target genes. As such, it regulates the response to injury, oxidative stress and inflammation,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

## Results



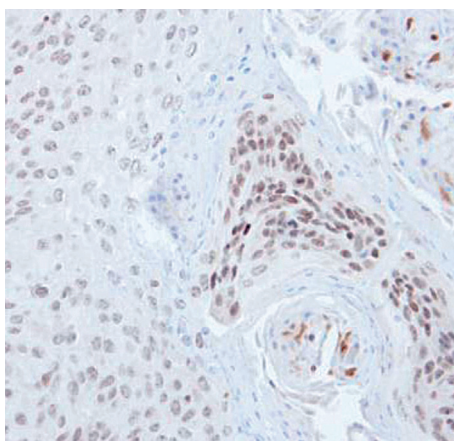
**Figure 1. ChIP results obtained with the Diagenode antibody directed against NRF2**

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.



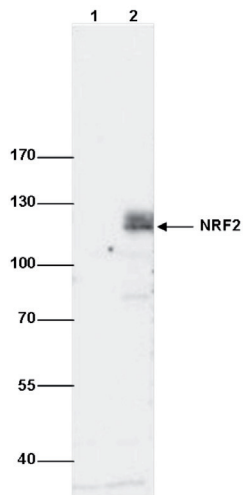
**Figure 2. Immunofluorescence using the Diagenode antibody directed against NRF2**

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



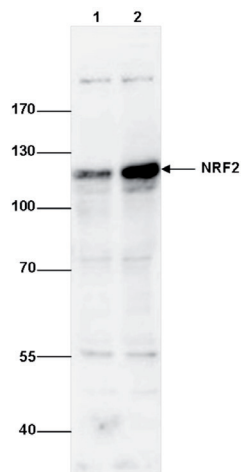
**Figure 3. Immunohistochemistry using the Diagenode antibody directed against NRF2**

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



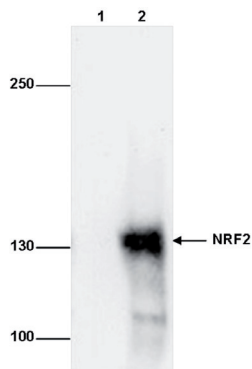
**Figure 4. Western blot analysis using the Diagenode antibody directed against NRF2**

Whole cell extracts from 293T cells (30 µg) 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 µg) treated with 30 µM 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 µg of the Diagenode antibody against NRF2 (Cat. No. C15410242) (lane 2) or with an equal amount of rabbit IgG (lane 1). The immunoprecipitated NRF2 protein was detected by western blot with the NRF2 antibody diluted 1:1,000.