

NRF1 monoclonal antibody

Other name: NRF-1, alpha PAL

Cat. No. C15200013

Type: Monoclonal ChIP-grade ChIP-seq grade

Isotype: IgG2a

Source: Mouse

Lot #: 001

Size: 50 µg/23 µl

Concentration: 2.2 µg/µl

Specificity: Human: positive. Other species: not tested.

Purity: Protein A purified monoclonal antibody in PBS containing 0.05% azide.

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: Monoclonal antibody raised in mouse against human NRF1 (Nuclear Respiratory Factor 1), using a recombinant protein.

Applications

	Suggested dilution	Results
ChIP*	2 µg per ChIP	Fig 1, 2
Western blotting	1:1,000	Fig 3

* Please note that the optimal antibody amount per IP should be determined by the end-user. We recommend testing 1-10 µg per IP.

Target description

NRF1 (UniProt/Swiss-Prot entry Q16656) is a transcription factor which activates the expression of some key metabolic genes regulating cellular growth and nuclear genes required for respiration, heme biosynthesis, and mitochondrial DNA transcription and replication. NRF1 has also been associated with the regulation of neurite outgrowth.

Results

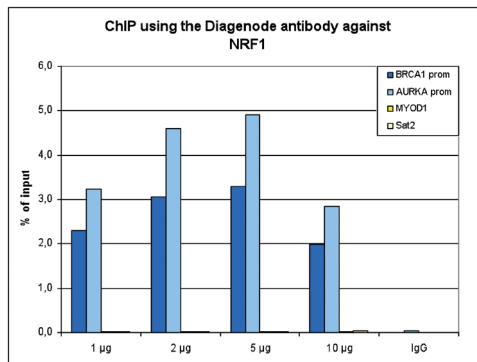


Figure 1. ChIP results obtained with the Diagenode monoclonal antibody directed against NRF1

ChIP was performed using K562 cells, the Diagenode monoclonal antibody against NRF1 (Cat. No. C15200013) and optimized PCR primer sets for qPCR. ChIP was performed with the “iDeal ChIP-seq” kit (cat. No. C01010055), using sheared chromatin from 4 million cells. A titration of the antibody consisting of 1, 2, 5 and 10 µg per ChIP experiment was analysed. IgG (2 µg/IP) was used as negative IP control. Quantitative PCR was performed with primers for the BRCA1 and AURKA promoters, used as positive controls, and for the coding region of the inactive MYOD1 gene and the Sat2 satellite repeat, used as negative controls. Figure 1 shows the recovery, expressed as a % of input (the relative amount of immunoprecipitated DNA compared to input DNA after qPCR analysis).

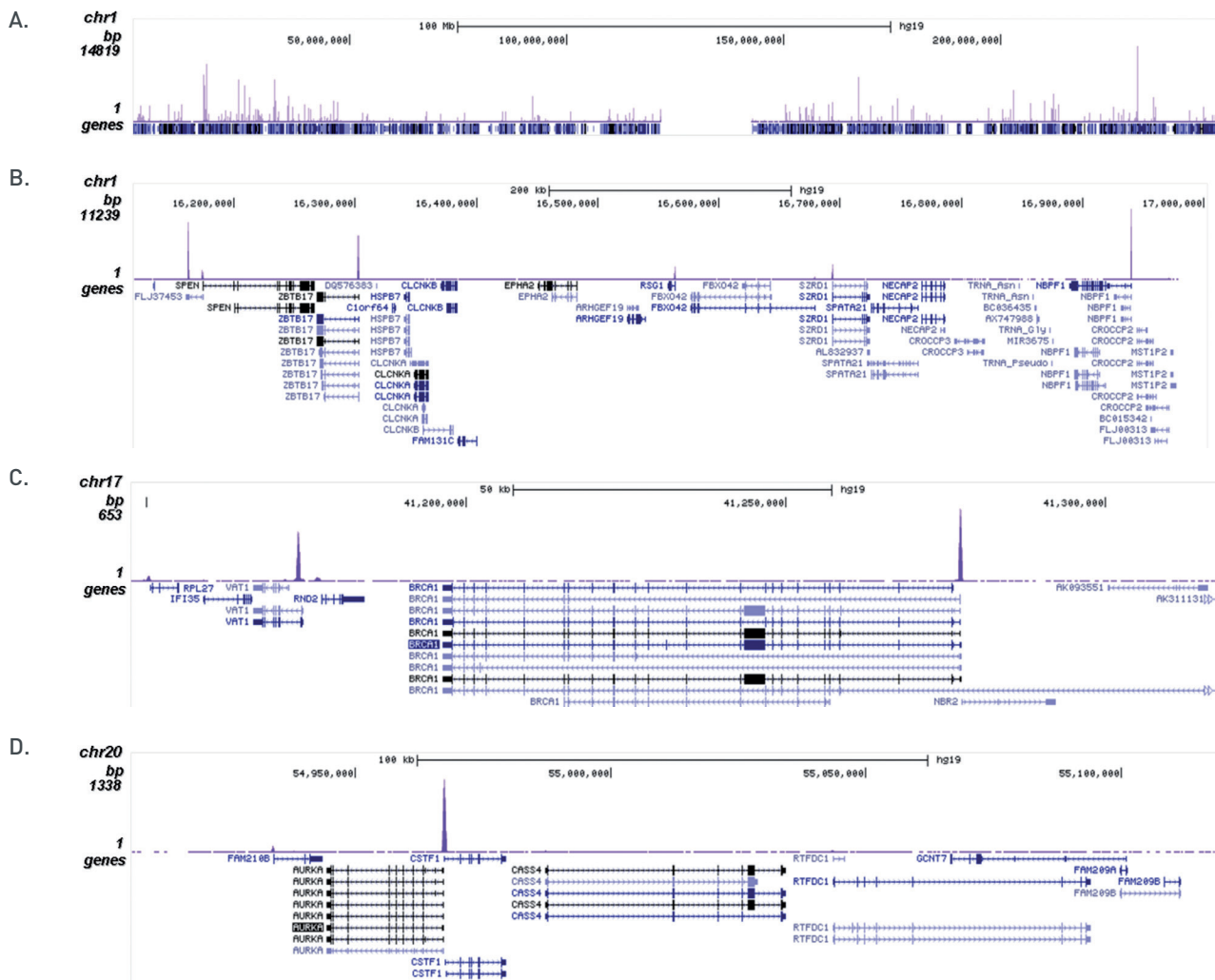


Figure 2. ChIP-seq results obtained with the Diagenode monoclonal antibody directed against NRF1
 ChIP was performed with 2 µg of the Diagenode antibody against NRF1 (Cat. No. C15200013) on sheared chromatin from 4 million K562 cells as described above. The IP'd DNA was subsequently analysed on an Illumina HiSeq 2000. Library preparation, cluster generation and sequencing were performed according to the manufacturer's instructions. The 50 bp tags were aligned to the human genome using the BWA algorithm. Figure 2 shows the peak distribution along the complete sequence and a 700 kb region of chromosome 1 (figure 2A and B) and in two regions surrounding the BRCA1 and AURKA positive control genes (figure 2C and D).

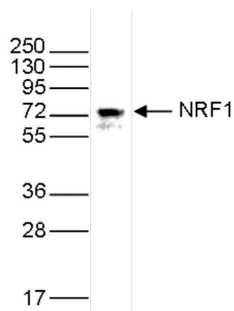


Figure 3. Western blot analysis using the Diagenode monoclonal antibody directed against NRF1
 Whole cell extracts from K562 cells were analysed by Western blot using the Diagenode antibody against NRF1 (Cat. No. C15200013) diluted 1:1,000 in TBST containing 5% milk powder. The position of the protein of interest is indicated on the right; the marker (in kDa) is shown on the left.