

## ZHX2 polyclonal antibody - Classic

**Other name:** AFR1, RAF

**Cat. No.** C15410260

**Type:** Polyclonal ChIP-grade, ChIP-seq grade

**Source:** Rabbit

**Lot #:** 40107

**Size:** 25 µl/100 µl

**Concentration:** 0.86 µg/µl

**Specificity:** Human: positive

**Purity:** Affinity purified polyclonal antibody in 0.1 M Tris-HCl containing 0.1 M Glycine, 10% glycerol and 0.01% thimerosal.

**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 ZHX2 (Zinc Fingers And Homeoboxes 2), using a recombinant protein.

### Applications

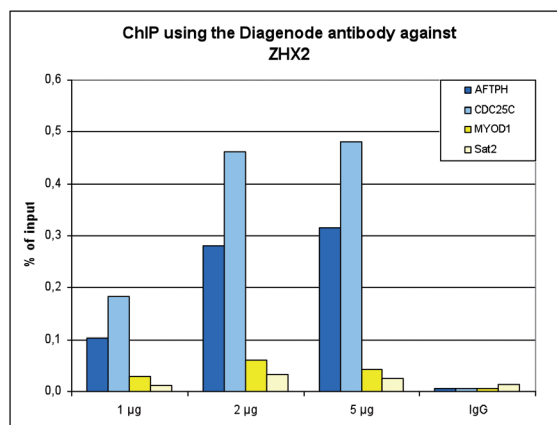
	Suggested dilution*	Results
ChIP*	2 - 5 µg per ChIP	Fig 1, 2
Western blotting	1:500 - 1:3,000	Fig 3
Immunofluorescence	1:100 - 1:1,000	Fig 4
Immunohistochemistry	1:100 - 1:1,000	Fig 5

\* 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

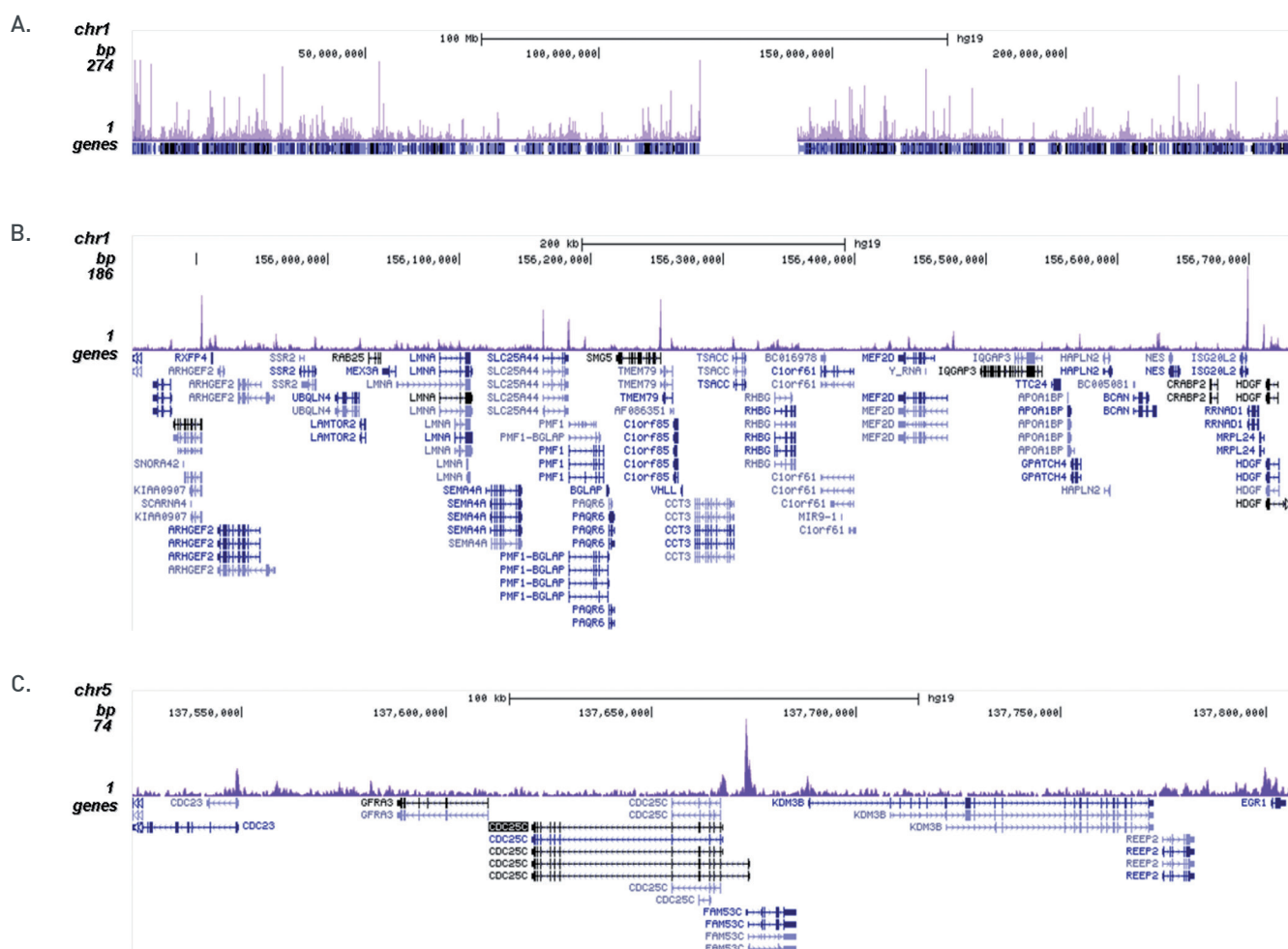
ZHX2 (UniProt/Swiss-Prot entry Q9Y6X8) is a transcriptional repressor that interacts with the A subunit of nuclear factor- $\kappa$ B (NF- $\kappa$ B). ZHX2 represses the promoter activity of the CDC25C gene.

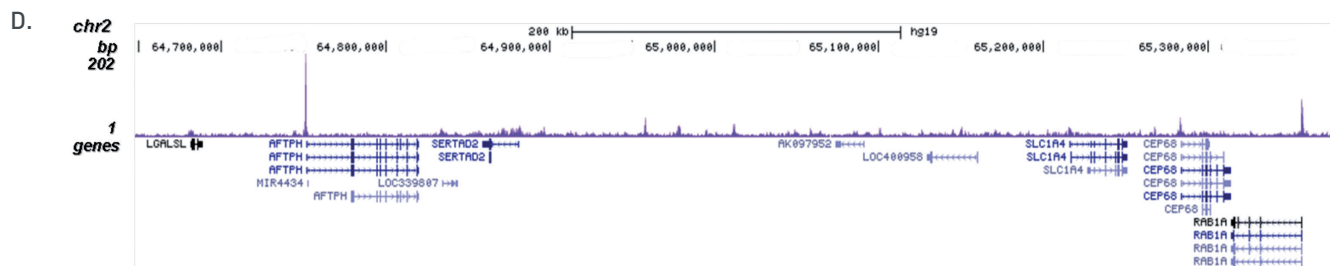
## TECHNICAL DATASHEET



**Figure 1. ChIP results obtained with the Diagenode antibody directed against ZHX2**

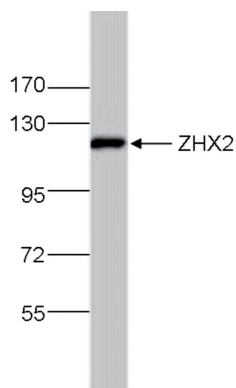
ChIP assays were performed using HepG2 cells, the Diagenode antibody against ZHX2 (Cat. No. C15410260) and optimized primer sets for qPCR. ChIP was performed with the "iDeal ChIP-seq" kit (Cat. No. C01010055), using sheared chromatin from 4 million cells. A titration of the antibody consisting of 1, 2 and 5 µg per ChIP experiment was analysed. IgG (1 µg/IP) was used as negative IP control. QPCR was performed with primers for the CDC25C and AFTPH promoters, used as positive controls, and for the 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).





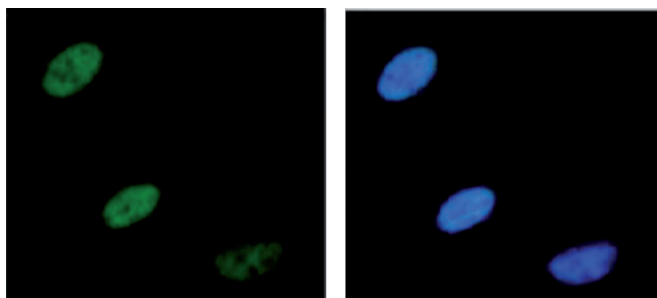
**Figure 2. ChIP-seq results obtained with the Diagenode antibody directed against ZHX2**

ChIP was performed on sheared chromatin from 5 million HepG2 cells using 5 µg of the Diagenode antibody against ZHX2 (Cat. No. C15410260). The IP'd DNA was subsequently analysed on an Illumina HiSeq. 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 enrichment along the complete sequence and a 1 Mb region of chromosomes 1 (fig 2A and B) and in two genomic regions surrounding the CDC25C and AFTPH genes on chromosome 5 and 2, respectively (fig 2C and D).



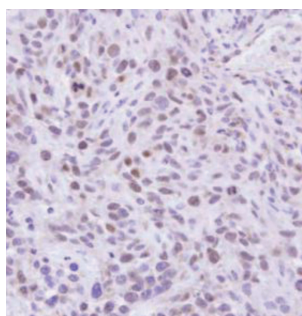
**Figure 3. Western blot analysis using the Diagenode antibody directed against ZHX2**

Whole cell extracts from HepG2 cells (30 µg) were analysed by Western blot using the Diagenode antibody against ZHX2 (Cat. No. C15410260) 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 4. Immunofluorescence with the Diagenode antibody directed against ZHX2**

HeLa cells were fixed with 4% formaldehyde for 15' at room temperature and stained with the Diagenode antibody against ZHX2 (Cat. C15410260) diluted 1:500 (left). The right picture shows costaining with Hoechst 33342 nucleic acid stain.



**Figure 5. Immunohistochemistry using the Diagenode antibody directed against ZHX2**

Formalin fixed paraffin embedded SAS xenograft tissue was stained with the Diagenode antibody against ZHX2 (Cat. No. C15410260) diluted 1:100 followed by a peroxidase labelled goat anti-rabbit secondary antibody.

**Diagenode sa. BELGIUM | EUROPE**

LIEGE SCIENCE PARK  
Rue Bois Saint-Jean, 3  
4102 Seraing (Ougrée) - Belgium  
Tel: +32 4 364 20 50  
Fax: +32 4 364 20 51  
orders@diagenode.com  
info@diagenode.com

**Diagenode Inc. USA | NORTH AMERICA**

400 Morris Avenue, Suite 101  
Denville, NJ 07834 - USA  
Tel: +1 862 209-4680  
Fax: +1 862 209-4681  
orders.na@diagenode.com  
info.na@diagenode.com

**Last update:** September 16, 2015