

EHF polyclonal antibody

Cat. No. C15410363

Type: Polyclonal	Specificity: Human: positive. Other species: not tested.
Size: 50 µg	Isotype: NA
Concentration: 0.8 µg/µl	Host: Rabbit
Lot No.: A2883P	Purity: Affinity purified polyclonal antibody.
Storage buffer: PBS containing 0.05% azide.	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.	

Last Data Sheet Update: June 20, 2019

Description

Other names: ESE3, ESE-3, ESE3B, ESEJ, HEHF

Polyclonal antibody raised in rabbit against human EHF (ETS Homologous Factor), using two synthetic peptides containing a sequence from the central part and the C-terminus of the protein, respectively.

Applications

Applications	Suggested dilution	References
ChIP/ChIP-seq*	1 – 2 µg per ChIP	Fig 1, 2
ELISA	1:1,000 - 1:50,000	Fig 3
Western Blotting	Not recommended	

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

Target Description

EHF (UniProtKB/Swiss-Prot entry Q9NZC4) belongs to the ETS transcription factor family. This transcriptional activator shows epithelial-specific expression and plays a role in regulating epithelial cell differentiation and proliferation. EHF may also act as a transcriptional repressor of a specific subset of ETS/AP-1-responsive genes and as a modulator of the nuclear response to mitogen-activated protein kinase signaling cascades. EHF is also a putative tumor suppressor gene and may be involved in carcinogenesis.

Validation data

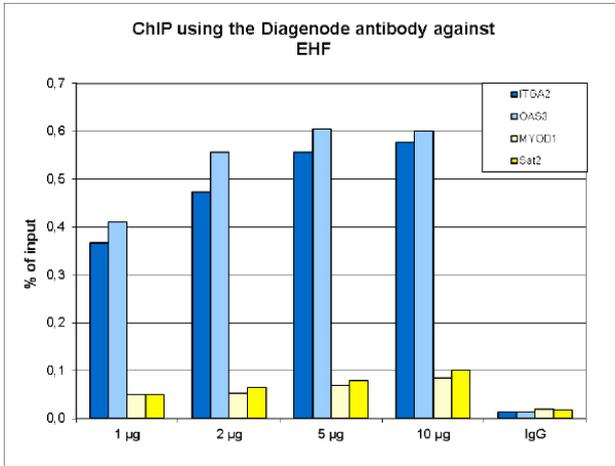
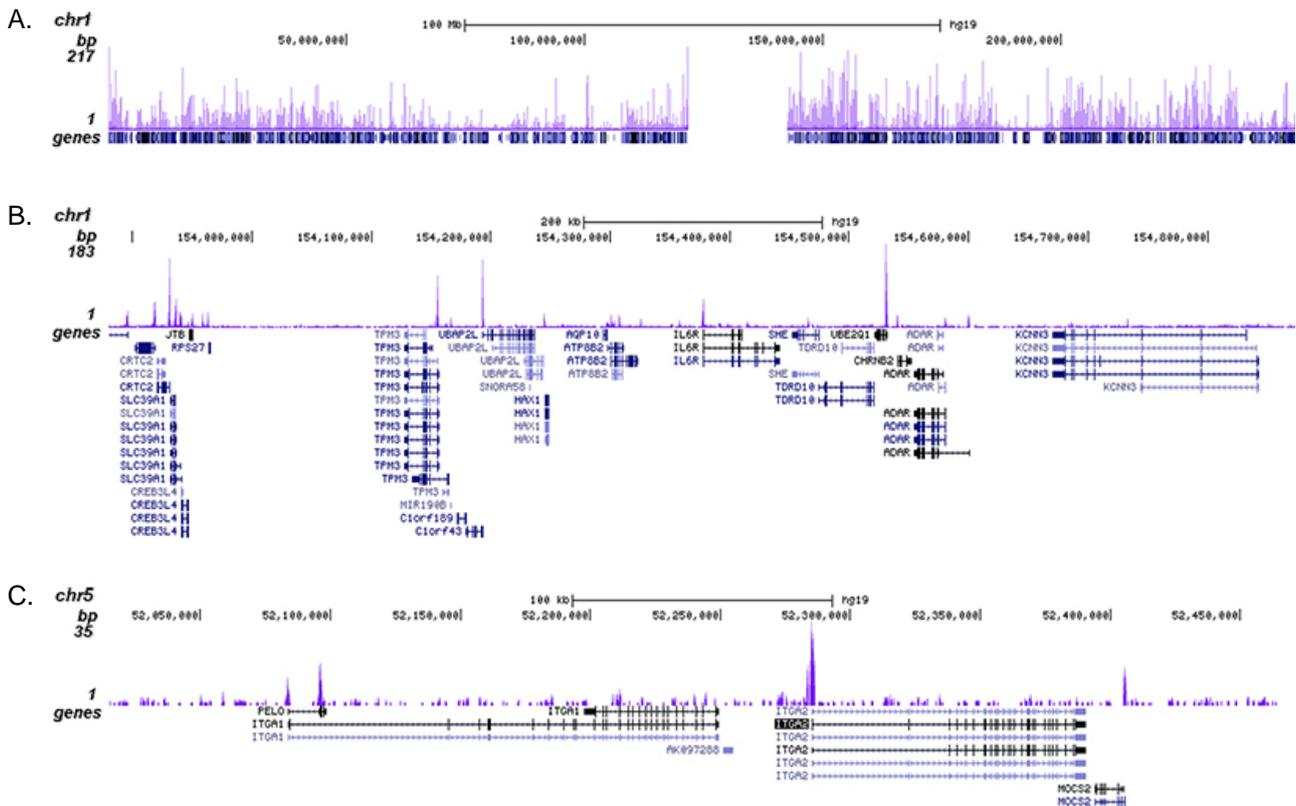


Figure 1. ChIP results obtained with the Diagenode antibody directed against EHF

ChIP assays were performed using T47D cells, the Diagenode antibody against EHF (Cat. No. C15410363) 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 consisting of 1, 2, 5 and 10 µg of antibody per ChIP experiment was analyzed. IgG (2 µg/IP) was used as a negative IP control. Quantitative PCR was performed with primers for the promoters of the ITGA2 and OAS3 genes, used as positive controls, and for the MYOD1 gene and the Sat2 satellite repeat, used as negative controls.



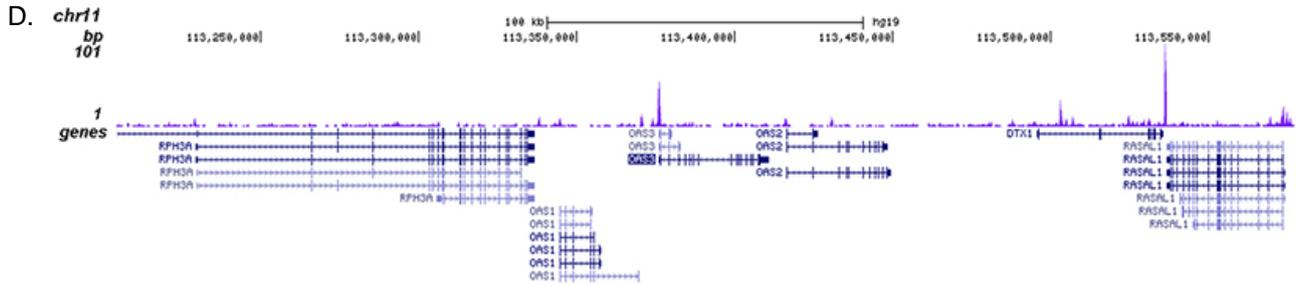


Figure 2. ChIP-seq results obtained with the Diagenode antibody directed against EHF

ChIP was performed on sheared chromatin from 4,000,000 T47D cells using 2 µg of the Diagenode antibody against EHF (Cat. No. C15410363) as described above. The IP'd DNA was subsequently analysed on an Illumina HiSeq 4000. 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 300 kb region of chromosome 1 (figure 2A and B) and in two genomic regions surrounding the ITGA2 and OAS3 positive control genes, respectively (figure 2C and D).

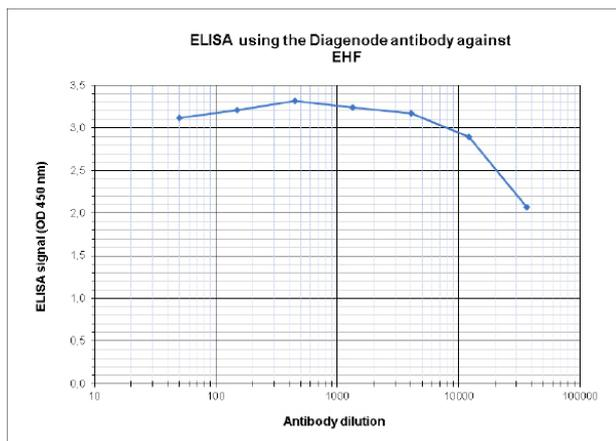


Figure 3. Determination of the antibody titer

To determine the titer of the antibody, an ELISA was performed using a serial dilution of Diagenode antibody directed against EHF (Cat. No. C15410363). The plates were coated with the peptides used for immunization of the rabbit. By plotting the absorbance against the antibody dilution (Figure 3), the titer of the antibody was estimated to be 1:98,000.