

AF9 polyclonal antibody

Other names: MLLT3, YEATS3

Cat. No. C15310266

Type: Polyclonal ChIP-grade / ChIP-seq grade

Source: Rabbit

Lot #: A1474-001

Size: 100 µl

Concentration: Not determined

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

Purity: Whole antiserum from rabbit 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: Polyclonal antibody raised in rabbit against human AF9 (Super Elongation Complex Subunit) using a KLH-conjugated synthetic peptide containing a sequence from the central region of the protein.

Applications

	Suggested dilution	Results
ChIP	2 µl/ChIP	Fig 1, 2
ELISA	1:100 - 1:500	Fig 3

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

Target description

AF9 (UniProtKB/Swiss-Prot entry P42568) is a component of the super elongation complex which is required to increase the catalytic rate of RNA polymerase II transcription by suppressing transient pausing by the polymerase at multiple sites along the DNA.

Results

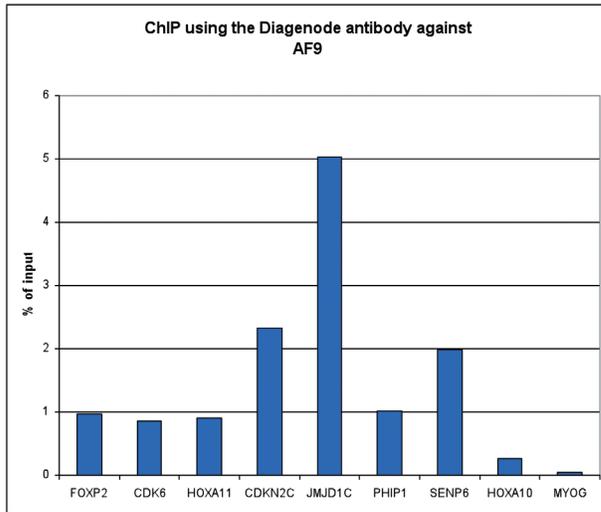


Figure 1. ChIP results obtained with the Diagenode antibody directed against AF9

ChIP was performed on THP-1 cells using the Diagenode antibody against AF9 (Cat. No. C15310266). Sheared chromatin from 1 million cells and 2 μ l of antibody were used per ChIP experiment. QPCR was performed using primers specific for the indicated genes. Figure 1 shows the recovery (the relative amount of immunoprecipitated DNA compared to input DNA).

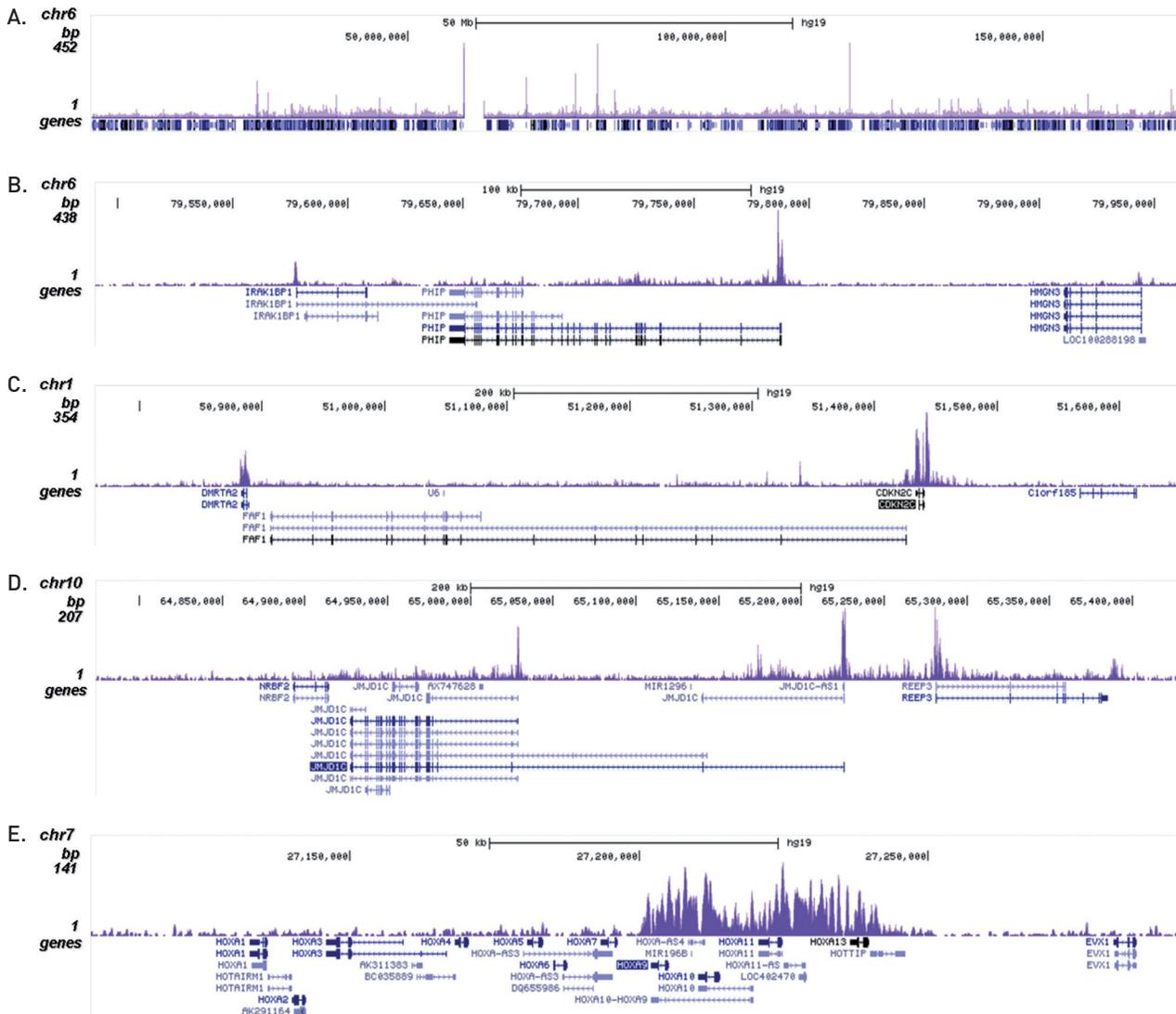


Figure 2. ChIP-seq results obtained with the Diagenode antibody directed against AF9

ChIP was performed as described above. The IP'd DNA of 5 ChIP's was pooled and 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 400 kb region of chromosome 6 containing the PHIP positive control gene (fig 2A and B), and in 3 genomic regions surrounding the CDKN2C gene, the JMJD1C gene and HOX cluster on chromosome 1, 10 and 7, respectively (fig 2C, D and E).

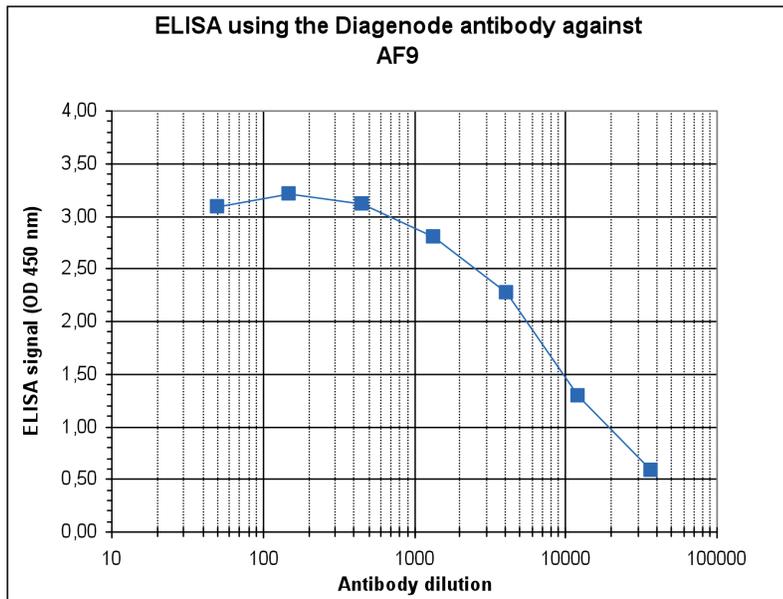


Figure 3. Determination of the antibody titer

To determine the titer of the antibody, an ELISA was performed using a serial dilution of the Diagenode antibody directed against AF9 (Cat. No. C15310266). The plates were coated with the peptide 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 1:8,000.

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: May 4, 2017