

Pol II S5p monoclonal antibody - Classic

Cat. No. C15200007-50

Type: Monoclonal Chip grade, ChIP-seq grade	Specificity: Human
Isotype: NA	Concentration: 1.0 µg/µl
Source: Mouse	Purity: Affinity purified monoclonal antibody in PBS containing 0.05% azide.
Lot No.: 001-12	Storage: Store at -20°C; for long storage, store at -80°C. Avoid multiple freeze-thaw cycles.
Size: 50 µg/50 µl	Precautions: This product is for research use only. Not for use in diagnostic or therapeutic procedures.

Last Data Sheet Update: January 17, 2017

Description

Alternative names: **POLR2A, RPB1, POLR2, RPOL2**

Monoclonal antibody raised in mouse against the YSPTSPS repeat in the B1 subunit of RNA polymerase II, phosphorylated at Ser5 of the repeat sequence.

Applications

Applications	Suggested dilution	References
ChIP *	1-2 µg/ChIP	Fig 1, 2
ELISA	1:3,000	Fig 3
Western Blotting	1:1,000	Fig 4, 5
Immunofluorescence	1:500	Fig 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

RNA polymerase II (pol II) is a key enzyme in the regulation and control of gene transcription. It is able to unwind the DNA double helix, synthesize RNA, and proofread the result. Pol II is a complex enzyme, consisting of 12 subunits, of which the B1 subunit (UniProt/Swiss-Prot entry P24928) is the largest. Together with the second largest subunit, B1 forms the catalytic core of the RNA polymerase II transcription machinery.

Validation Data

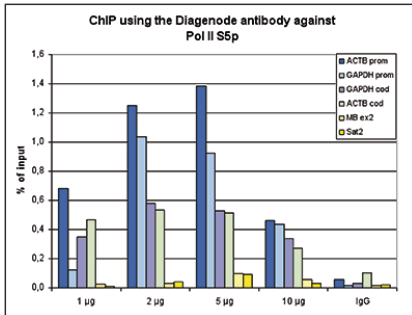


Figure 1. ChIP results obtained with the Diagenode monoclonal antibody directed against Pol II S5p

ChIP assays were performed using human HeLa cells, the Diagenode monoclonal antibody against Pol II S5p (Cat. No. C15200007) and optimized PCR primer pairs for qPCR. ChIP was performed with the “iDeal ChIP-seq” kit (Cat. No. C01010051), using sheared chromatin from 1 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 specific for the promoter and the coding region of the constitutively expressed GAPDH and ACTB genes, used as positive controls, and for exon 2 of the inactive myoglobin (MB) 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).

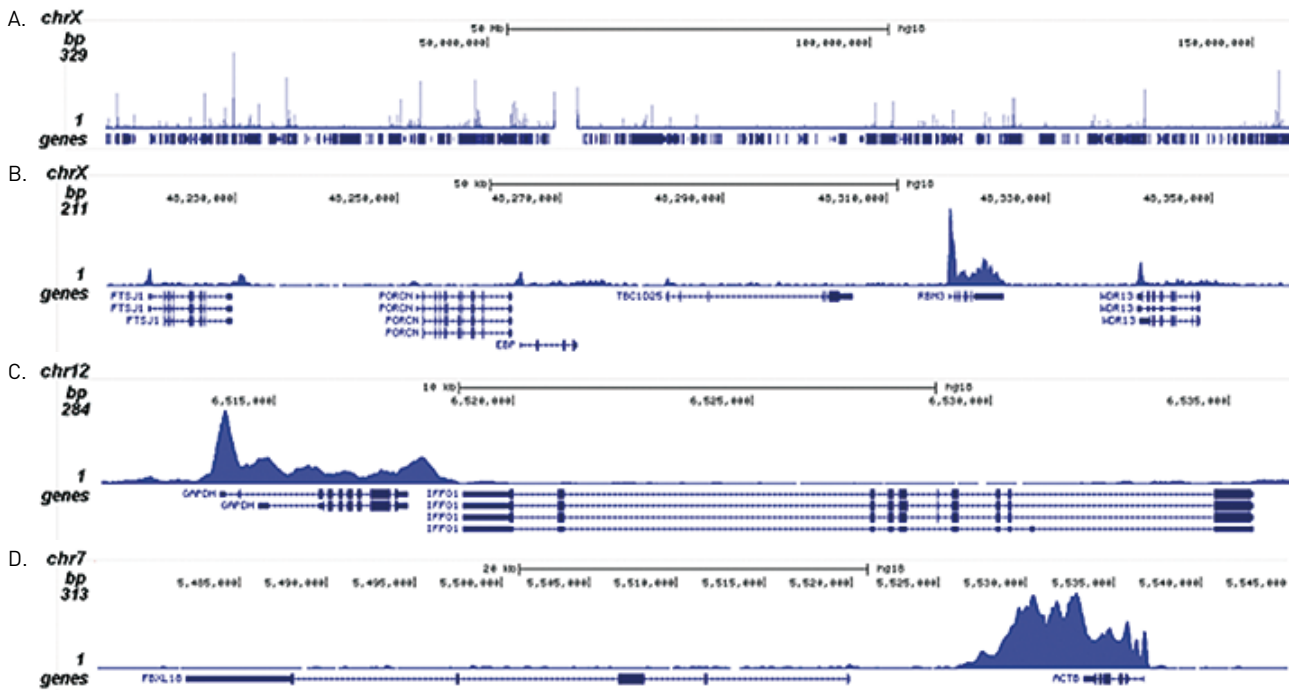


Figure 2. ChIP-seq results obtained with the Diagenode monoclonal antibody directed against Pol II S5p

ChIP was performed on sheared chromatin from 1 million HeLaS3 cells using 1 µg of the Diagenode antibody against Pol II S5p (Cat. No. C15200007) as described above. The IP'd DNA was subsequently analysed on an Illumina Genome Analyzer. Library preparation, cluster generation and sequencing were performed according to the manufacturer's instructions. The 36 bp tags were aligned to the human genome using the ELAND algorithm. Figure 2 shows the peak distribution along the complete sequence and a 150 kb region of the X-chromosome (figure 2A and B, respectively), and in a two genomic regions surrounding the GAPDH and ACTB positive control genes (figure 2C and D).

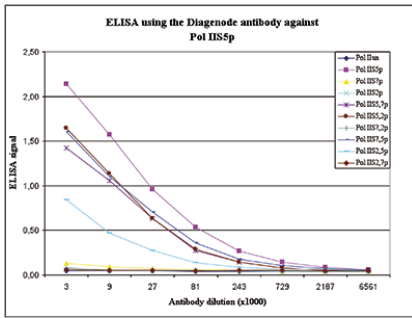


Figure 3. Cross reactivity of the Diagenode monoclonal antibody directed against Pol II S5p

To test the specificity an ELISA was performed using a serial dilution of the Diagenode monoclonal antibody against Pol II S5p (Cat. No. C15200007). The wells were coated with peptides containing the unmodified C-terminal repeat sequence as well as different phosphorylated peptides. Figure 3 shows the specificity of the antibody for the S5 phosphorylation.

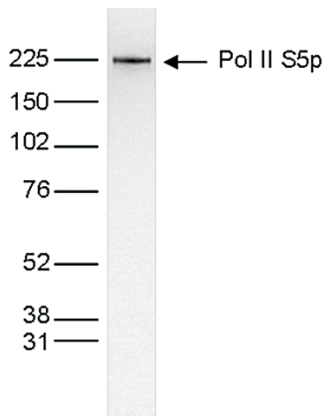


Figure 4. Western blot analysis using the Diagenode monoclonal antibody directed against Pol II S5p

Nuclear extracts (25 µg) from HeLa cells were analysed by Western blot using the Diagenode monoclonal antibody against Pol II S5p (Cat. No. C15200007) diluted 1:1,000 in TBS-Tween containing 5% skimmed milk. The position of the protein of interest is indicated on the right; the marker (in kDa) is shown on the left.

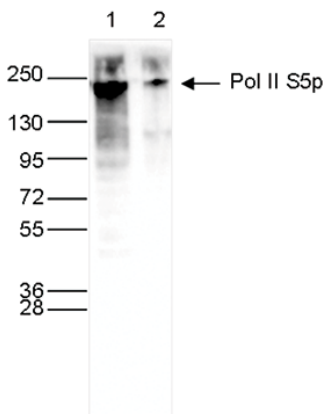


Figure 5. Western blot analysis using the Diagenode monoclonal antibody directed against Pol II S5p

Whole cell extracts (40 µg) from HeLa cells transfected with Pol II siRNA (lane 2) and from an untransfected control (lane 1) were analysed by Western blot using the Diagenode antibody against Pol II S5p (Cat. No. C15200007) diluted 1:1,000 in TBS-Tween containing 5% skimmed milk. The position of the protein of interest is indicated on the right; the marker (in kDa) is shown on the left.



Figure 6. Immunofluorescence using the Diagenode monoclonal antibody directed against Pol II S5p

HeLa cells were stained with the Diagenode antibody against Pol II S5p (Cat. No. C15200007) and with DAPI. Cells were fixed with methanol and blocked with PBS/TX-100 containing 5% normal goat serum and 1% BSA. The cells were immunofluorescently labelled with the Pol II S5p antibody (left) diluted 1:500 in blocking solution followed by an anti-mouse antibody conjugated to Alexa594. The middle panel shows staining of the nuclei with DAPI. A merge of the two stainings is shown on the right.