

AbFlex[®] RNA pol II CTD phospho Ser5 antibody (rAb) (biotin)

Catalog Nos: 91117, 91118

RRID: AB_2793781

Isotype: IgG2a

Application(s): ChIP, ChIP-Seq, ELISA, ICC, IF, WB

Reactivity: Human

Quantities: 100 µg, 10 µg

Purification: Ni-NTA

Host: Mouse

Concentration: 1 µg/µl

Molecular Weight: 240 kDa

Background: AbFlex[®] antibodies are recombinant antibodies (rAbs) that have been generated using defined DNA sequences to produce highly specific, reproducible antibodies. Each AbFlex antibody contains a 6xHis-Tag, an avidin tag sequence for enzymatic biotin conjugation using the biotin ligase, BirA, and a sortase recognition motif (LPXTG) to attach a variety of labels directly to the antibody including fluorophores, enzymatic substrates (HRP, AP), peptides, drugs as well as solid supports.

AbFlex[®] RNA pol II CTD phospho Ser5 antibody (biotin) was expressed as full-length IgG with mouse immunoglobulin heavy and light chains (IgG2a isotype) in mammalian 293 cells. The antibody was directly labeled with biotin using the biotin ligase, BirA.

RNA pol II (RNA polymerase II) is responsible for synthesizing messenger RNA in eukaryotes. RNA pol II contains a carboxy terminal domain composed of heptapeptide repeats that are essential for polymerase activity. These repeats contain serine and threonine residues that are phosphorylated in actively transcribing RNA polymerase. In addition, RNA pol II, in combination with several other polymerase subunits, form the DNA binding domain of the polymerase, a groove in which the DNA template is transcribed into RNA.

During the transcription cycle, the CTD of the large subunit of RNA pol II is reversibly phosphorylated. RNA pol II containing unphosphorylated CTD is recruited to the promoter, whereas the hyperphosphorylated CTD form is involved in active transcription. Phosphorylation occurs at two sites within the heptapeptide repeat, at serine 2, serine 5 and serine 7. RNA pol II Serine 5 phosphorylation is confined to promoter regions and is necessary for the initiation of transcription.

Immunogen: This RNA pol II CTD phospho Ser5 antibody was raised against a peptide containing the RNA pol II CTD sequence phosphorylated at serine 5.

Buffer: Purified IgG in 50mM sodium phosphate pH 8.0, 150mM NaCl, and 100mM imidazole with 30% glycerol and 0.035% sodium azide.

Application Notes:

Validated Applications:

ChIP-Seq: 4 µg each

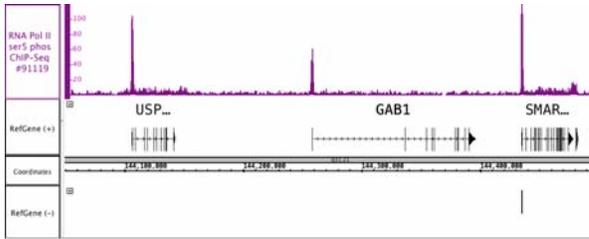
ICC/IF: 1 - 2 µg / mL

WB: 0.5 - 2 µg/mL

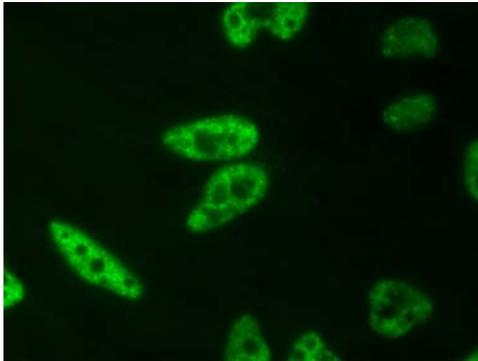
Bead-based ELISA: 0.1 - 5 µg/ml

Storage and Guarantee: Some products may be shipped at room temperature. This will not affect their stability or performance. Upon receipt, unconjugated antibodies may be stored at -20°C for up to 2 years. Fluorophore- & enzyme-conjugated antibodies should be stored at 4°C. Fluorophore-conjugated antibodies should be protected from light. Keep reagents on ice when not in storage; to avoid repeated freeze/thaw cycles, we recommend aliquoting items that will be stored frozen into single-use fractions prior to freezing. This product is guaranteed for 6 months from date of receipt.

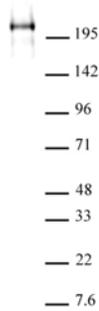
This product is for research use only and is not for use in diagnostic procedures.



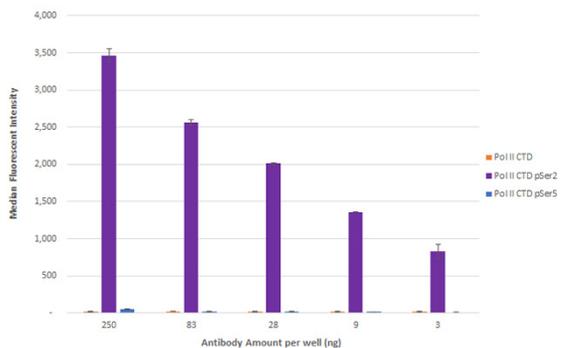
AbFlex[®] RNA pol II CTD phospho Ser5 antibody (rAb) tested by ChIP-Seq. ChIP was performed using the ChIP-IT[®] High Sensitivity Kit (Cat. No. 53040) with 20 µg of chromatin from HL60 cells and 4 µg of AbFlex[®] RNA pol II CTD phospho Ser5 antibody. ChIP DNA was sequenced on the Illumina HiSeq and 20 million sequence tags were mapped to identify RNA pol II phospho Ser5 binding. The image shows binding across a region of chromosome 4.



AbFlex[®] RNA pol II CTD phospho Ser5 antibody (rAb) tested by immunofluorescence. HeLa cell stained with 2µg /mL of AbFlex[®] RNA pol II CTD phospho Ser5 antibody (rAb) followed by anti-mouse-IgG-488.



AbFlex[®] RNA pol II CTD phospho Ser5 antibody (rAb) tested by Western blot. HeLa nuclear extract (20 µg per lane) probed with AbFlex[®] RNA pol II CTD phospho Ser5 antibody (0.5 µg/ml dilution).



AbFlex[®] RNA pol II CTD phospho Ser5 antibody (rAb) tested by Luminex bead-based specificity analysis.

Luminex bead-based specificity analysis was used to confirm the specificity of AbFlexTM RNA pol II CTD phospho Ser5 antibody (rAb) antibody for RNA pol II CTD peptides. RNA Pol II peptides were conjugated to MagPlex Luminex beads and incubated with various amounts of AbFlexTM RNA pol II CTD phospho Ser5 antibody (rAb). Peptide-bound antibody was detected with anti-mouse IgG-Phycoerythrin and read in a Luminex instrument.

Luminex[®] is a registered trademark of Luminex Corporation.