

AbFlex[®] Histone H3, C-terminal antibody (rAb, Rabbit)

Catalog Nos: 91299, 91300

Isotype: IgG

Application(s): ChIP, WB

Reactivity: Human

Quantities: 100 µg, 10 µg

Purification: Protein A Chromatography

Host: Rabbit

Concentration: 1 µg/µl

Molecular Weight: 17 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, a Biotinylation Tag 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[®] Histone H3, C-term was expressed as full-length IgG with rabbit immunoglobulin heavy and light chains in CHO cells.

Histone H3 is one of the core components of the nucleosome. The nucleosome is the smallest subunit of chromatin and consists of 147 base pairs of DNA wrapped around an octamer of core histone proteins (two each of Histone H2A, Histone H2B, Histone H3 and Histone H4).

Immunogen: This antibody was raised against a peptide at the C-terminus of human Histone H3.

Buffer: Purified IgG in 140 mM Hepes, pH 7.5, 70 mM NaCl, 32 mM NaOAc, 0.035% sodium azide, 30% glycerol. Sodium azide is highly toxic.

Application Notes:

Applications Validated by Active Motif:

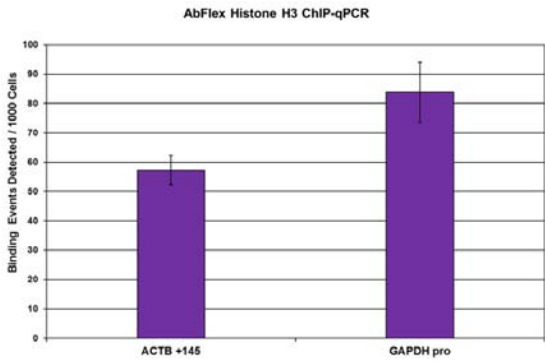
ChIP: 4 - 5 µg per ChIP

WB*: 0.01 – 0.5 µg/ml

*Note: Many chromatin-bound proteins are not soluble in a low salt nuclear extract and fractionate to the pellet. Therefore, we recommend a High Salt / Sonication Protocol when preparing nuclear extracts for Western Blot.

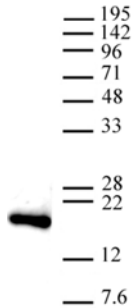
Storage and Guarantee: Some products may be shipped at room temperature. This will not affect their stability or performance. Avoid repeated freeze/thaw cycles by aliquoting items into single-use fractions for storage at -20°C for up to 2 years. Keep all reagents on ice when not in storage. This product is guaranteed for 12 months from date of receipt.

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



AbFlex® Histone H3 antibody (rAb) tested by ChIP

Chromatin immunoprecipitation (ChIP) was performed with 30 µg of chromatin from Farage human lymph-node derived B cells and 4.5 µg AbFlex Histone H3 antibody. ChIP DNA was used in qPCR with the gene-specific primers as indicated. Data are presented as Binding Events Detected per 1000 Cells using Active Motif's Epigenetic Services normalization scheme which accounts for primer efficiency and the amount of chromatin used in the ChIP reaction.



AbFlex® Histone H3, C-term antibody tested by Western blot.

20 µg of HeLa nuclear extract* was run on SDS-PAGE and probed with AbFlex® Histone H3, C-term antibody at 0.1 µg/ml.