

AbFlex® HP1β antibody (rAb)

Catalog Nos: 91407, 91408**RRID:** AB_3216363**Application(s):** ELISA, ICC, IF, IP, WB**Reactivity:** Human, Mouse**Quantities:** 100 µg, 10 µg**Purification:** Protein A Chromatography**Host:** Mouse**Isotype:** IgG2a**Molecular Weight:** 25 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®HP1 beta antibody was expressed as full-length IgG with mouse immunoglobulin heavy and light chains (IgG2a isotype) in mammalian 293 cells.

HP1 beta (Heterochromatin Protein 1 beta, Chromobox protein 1, CBX1) is one of the mammalian homologues of Drosophila HP1, first identified through its role in the maintenance of regions of the genome in a heterochromatic state. HP1 beta contains a chromo domain and a chromo shadow domain. Through the chromo domain, HP1 proteins are recruited to regions of chromatin that exhibit methylation of lysine 9 of histone H3. Interaction between HP1 beta and the lamin B receptor can facilitate the sequestration of heterochromatin to the nuclear envelope.

Immunogen: This HP1β antibody was raised against recombinant mouse HP1β.

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

Application Notes:

Applications Validated by Active Motif:

WB: 0.5-2 µg/ml dilution

AbFlex® recombinant antibodies are genetically derived from DNA sequences of parental hybridoma clones. For details on the parental antibody clone, see Catalog No. 39979.

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® HP1 beta antibody (rAb) tested by Western blot.

20 µg of HeLa whole cell extract was run on SDS-PAGE and probed with AbFlex HP1 beta antibody at 0.5 µg/ml.