

AbFlex® HIF-1 alpha antibody (rAb)

Catalog Nos: 91139, 91140

RRID: AB_2793787

Application(s): ELISA

Reactivity: Bovine, Human, Mouse, Rat

Quantities: 100 µg, 10 µg

Purification: Ni-NTA

Host: Mouse

Isotype: IgG2a

Concentration: 1 µg/µl

Molecular Weight: 100 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® HIF-1 alpha antibody was expressed as full-length IgG with mouse immunoglobulin heavy and light chains (IgG2a isotype) in mammalian 293 cells.

HIF-1 (hypoxia inducible factor 1, alpha subunit, MOP1, PASD8, bHLHe78, PAS domain-containing protein 8) is a transcription factor found in mammalian cells cultured under reduced oxygen tension that plays an essential role in cellular and systemic homeostatic responses to hypoxia. HIF-1 is a heterodimer composed of an alpha subunit and a beta subunit. The beta subunit has been identified as the aryl hydrocarbon receptor nuclear translocator (ARNT). This gene encodes the alpha subunit of HIF-1. Overexpression of a natural antisense transcript (aHIF) of this gene has been shown to be associated with nonpapillary renal carcinomas. Two alternative transcripts encoding different isoforms have been identified.

Immunogen: This HIF-1 alpha antibody was raised against a partial protein containing amino acid residues 329-530 of human HIF-1 alpha.

Buffer: Purified IgG 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:

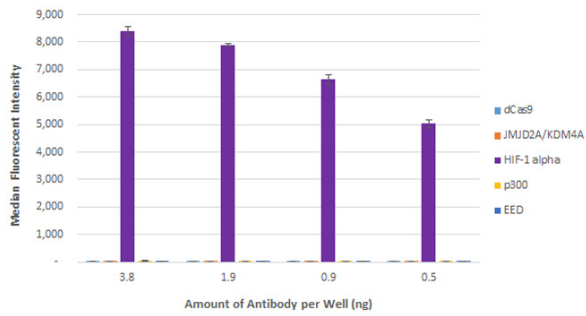
Applications Validated by Active Motif:

Bead-based ELISA: 10 - 76 ng/ml

AbFlex® recombinant antibodies are genetically derived from DNA sequences of parental hybridoma clones. For details on the parental clone, see Catalog No. 61275.

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[®] HIF-1 alpha antibody (rAb) tested by Luminex bead-based specificity analysis. Luminex bead-based specificity analysis was used to confirm the specificity of AbFlex[®] HIF-1 alpha antibody (rAb) antibody for HIF-1 alpha. Various proteins were conjugated to MagPlex Luminex beads and incubated with various amounts of AbFlex[®] HIF-1 alpha antibody (rAb). Protein-bound antibody was detected with anti-mouse IgG-Phycoerythrin and read in a Luminex instrument.

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