

AbFlex[®] His-Tag antibody (rAb)

Catalog Nos: 91183, 91184

RRID: AB_2793793

Isotype: IgG2a

Application(s): WB

Reactivity: Not Species Specific

Quantities: 100 µg, 10 µg

Purification: Protein A Chromatography

Host: Mouse

Concentration: 1 µg/µl

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[®] His-Tag antibody was expressed as full-length IgG with mouse immunoglobulin heavy and light chains (IgG2a isotype) in mammalian 293 cells. The 6xHis-Tag that is normally present on the AbFlex antibody was removed to avoid self binding.

Immunogen: The His-Tag is comprised of a series of six or more histidine residues that are used to label the carboxy or amino terminus of a recombinant protein. This antibody is designed to recognize the His-Tag (HHHHHH).

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.

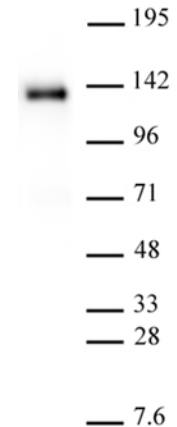
Application Notes:

Applications Validated by Active Motif:

WB: 0.5 - 2 µg/ml

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[®] His-Tag antibody (rAb) tested by Western blot. 250 ng of a His-Tagged protein (Lane 1) or Flag-tagged protein (Lane 2: Recombinant Cas9, Cat. No. 31457) was run under reducing conditions and probed with AbFlex[®] His-Tag antibody (0.5 µg/ml dilution). The AbFlex[®] His-Tag antibody specifically recognizes only the His-tagged protein.