

AbFlex® HA-Tag antibody (rAb)

Catalog Nos: 91271, 91272**RRID:** AB_2793828**Isotype:** IgG2a**Application(s):** IP**Reactivity:** Human, 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® HA-Tag antibody was expressed as full-length IgG with mouse immunoglobulin heavy and light chains (IgG2a isotype) in mammalian 293 cells.

HA-Tag has been derived from the hemagglutinin protein of the human influenza virus for use as an epitope for the detection and purification of tagged recombinant proteins expressed *in vivo*. The HA-Tag is a small epitope and generally does not interfere with the normal function of recombinant proteins.

Immunogen: This antibody was raised against a peptide containing multiple repeats of the HA tag epitope YPYDVPDYA.

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:

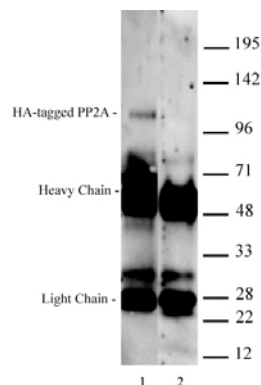
Applications Validated by Active Motif:

IP: 2 - 5 µg per IP

AbFlex® recombinant antibodies are genetically derived from DNA sequences of parental hybridoma clones. For details on the parental clone, see Catalog No. 39627.

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® HA-Tag antibody (rAb) tested by Immunoprecipitation.

Whole-cell lysate (400 µg) from U2OS cells expressing an HA-Tagged protein, was immunoprecipitated with 2 µg of AbFlex HA-Tag rAb (lane 1) or control IgG2a (lane 2). Immunoprecipitates were run on SDS-PAGE and analyzed by Western blot using HA-Tag mAb Cat. No. 39627 at 4 µg/ml.