

## HA-Tag antibody (mAb)

**Catalog Nos:** 39627, 39628

**RRID:** AB\_2793281

**Clone:** 3B9

**Isotype:** IgG1

**Application(s):** IP

**Reactivity:** Human, Not Species Specific

**Quantities:** 200 µg, 10 µg

**Purification:** Protein G Chromatography

**Host:** Mouse

**Concentration:** 1 µg/µl

**Molecular Weight:** Not Applicable

**Background:** 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 HA-Tag antibody was raised against a peptide containing multiple repeats of the HA-Tag epitope, YPYDVPDYA.

**Buffer:** Purified IgG in 70 mM Tris (pH 8), 105 mM NaCl, 31 mM glycine, 0.07 mM EDTA, 30% glycerol and 0.035% sodium azide. Sodium azide is highly toxic.

### Application Notes:

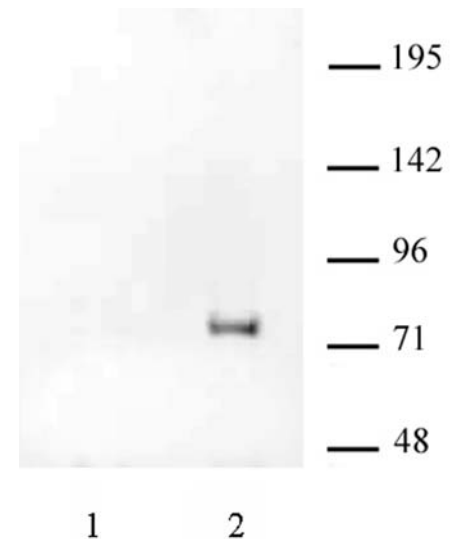
Applications Validated by Active Motif:

IP: 2 - 5 µg per IP

This antibody is also available as an AbFlex<sup>®</sup> engineered recombinant antibody. For details on the corresponding AbFlex Recombinant Antibody, see Catalog No. 91271.

**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.



### HA-Tag mAb tested by IP / Western blot.

Whole-cell lysate (400 µg) from U2OS cells expressing an HA-Tagged protein was immunoprecipitated with 2 µg of either control mouse IgG (lane 1) or HA-Tag mAb (lane 2). Immunoprecipitates were run on an SDS-PAGE gel and analyzed by Western blot using a second HA-Tag antibody.