

## AbFlex® HSP72 antibody (rAb)

**Catalog Nos:** 91303, 91304**Isotype:** IgG**Application(s):** WB**Reactivity:** Human**Quantities:** 100 µg, 10 µg**Purification:** Protein A Chromatography**Host:** Mouse**Concentration:** 1 µg/µl**Molecular Weight:** 72 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® HSP72 antibody was expressed in CHO cells and contains mouse immunoglobulin heavy and light chains.

HSP72 (Heat Shock Protein 72) is a member of the DnaK/HSP70 family of protein chaperones involved in protein homeostasis. HSP72 is expressed at low constitutive levels and is induced upon exposure to environmental stress. HSP72 released into the blood can serve as an alert for the immune system. Increased levels of HSP72 can repress expression of inflammatory response genes by inhibiting NF-κB.

**Immunogen:** This antibody was raised against full-length recombinant human HSPA1A heat shock protein family A (Hsp70) member 1A.

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

**Application Notes:**

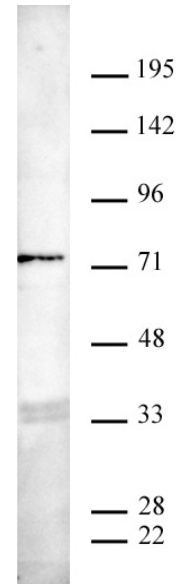
Applications Validated by Active Motif:

WB\*: 0.05 – 2 µg/ml

\*Note: Many chromatin-bound proteins are not soluble in a low salt nuclear extract and fractionate to the pellet. Therefore, we recommend a High Salt / Sonication Protocol when preparing nuclear extracts for Western blot.

**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® HSP72 antibody tested by Western blot. 20 µg of HEK293 nuclear extract\* was run on SDS-PAGE and probed with AbFlex HSP72 antibody at 2 µg/ml.