

## HSP72 antibody (mAb)

---

**Catalog Nos:** 61209, 61210

**RRID:** AB\_2793553

**Clone:** 3G7

**Isotype:** IgG2a

**Application(s):** ICC, IF, IP, WB

**Reactivity:** Human

**Quantities:** 100 µg, 10 µg

**Purification:** Protein G Chromatography

**Host:** Rat

**Concentration:** 1 µg/µl

**Molecular Weight:** 70 kDa

---

**Background:** 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 HSP72 antibody was raised against full-length recombinant human HSP72 protein.

**Buffer:** Purified IgG in PBS with 30% glycerol and 0.035% sodium azide. Sodium azide is highly toxic.

**Application Notes:**

For HSP72, we also offer AbFlex<sup>®</sup> HSP72 Recombinant Antibody (rAb). For details, see Catalog No. 91303.

**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.