

HSC70 / HSP7C antibody (mAb)

Catalog Nos: 61207, 61208

RRID: AB_2793552

Clone: 6E1

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: HSC70 / HSP7C (Heat Shock Cognate 70, HSP73, Heat Shock Protein 73) is a member of the DnaK/HSP70 family of protein chaperones involved in protein homeostasis. HSC70 is constitutively expressed and essential for cell viability and functions both in preventing or reversing protein aggregation and in disassembling protein complexes. HSC70 predominantly binds non-native polypeptides and folding intermediates that are either newly synthesized or stress-induced.

Immunogen: This HSC70 / HSP7C antibody was raised against full-length recombinant human HSC70 / HSP7C protein.

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

Application Notes:

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.