

Anti-HSP27

Mouse Monoclonal Antibody

Catalog # H30-60M

Lot # B3216-49

Cited Applications

ELISA, WB

Ideal working dilutions for each application should be empirically determined by the investigator.

Specificity

Recognizes the HSP27 protein

Cross Reactivity

Human

Host/Isotype/Clone#

Mouse, IgG1

Immunogen

The antibody was produced against prokaryotic recombinant protein corresponding to the full length human HSP27 protein

Formulation

0.02 M Potassium Phosphate, 0.15 M Sodium Chloride, pH 7.2 + 0.01% (w/v) Sodium Azide

Stability

1yr at -20°C from date of shipment

Scientific Background

HSP27 is a member of heat shock proteins (HSPs) which are synthesized in cells in response to heat shock and other metabolic stresses and provide a transient state of thermotolerance. HSP27 plays a major role in the increased thermal resistance acquired by cells after exposure to HSP inducers (1). The level of HSP27 phosphorylation is significantly elevated after exposure of cells to heat shock, sodium arsenite, IL-1 and TNF- α . MAPKAPK2 and MAPKAPK3 are both activated by these conditions and can phosphorylate HSP27 on serine residues (2).

References:

1. Landry, J. et al: Heat shock resistance conferred by expression of the human HSP27 gene in rodent cells. J Cell Biol. 1989 Jul; 109(1): 7-15.
2. Clifton, A D. et al: A comparison of the substrate specificity of MAPKAP kinase-2 and MAPKAP kinase-3 and their activation by cytokines and cellular stress. FEBS Lett. 1996; 392(3): 209-14.

Sample Data



Western Blot using Anti-HSP27 (1:1000) antibody to detect human HSP27. 10 µg of HeLa whole cell lysate loaded per lane

Anti-HSP27

Mouse Monoclonal Antibody

Catalog # **H30-60M**

Lot # **B3216-49**

Purification Protein A Chromatography

Stability 1yr at -20°C from date of shipment

Storage & Shipping Store product at -20°C. For optimal storage, aliquot antibody into smaller quantities after centrifugation and store at recommended temperature. For optimal performance, avoid repeated handling and multiple freeze/thaw cycles. Product shipped on ice packs.