

HSP27 (pSer78) Polyclonal Antibody

(Catalog No. A-7427)

Background

HSP27 is a member of the small heat shock proteins (sHsp) family, along with alphaB-crystallin. It is involved in stress resistance and actin organization and is associated with CRYAB, alpha- and beta-tubulin, and microtubules. HSP27 can also interact with other heat shock proteins, such as HSPB8 and HSPBAP1. It is usually expressed in response to certain environmental stresses like heat shock, or as reaction to estrogen stimulation in the MCF-7 cells. Expression is seen throughout most tissues with the highest amounts being in the heart and in the striated and smooth muscle tissue. Overexpression of HSP27 has been linked to certain cancers including breast, prostate, gastric, ovarian, and urinary bladder cancers.

Description

Rabbit polyclonal antibody to HSP27 (pSer78)

Formulation

Liquid. In PBS containing 50% glycerol and 0.09% sodium azide.

Immunogen

Synthetic peptide corresponding to a portion of human HSP27 phosphorylated at Ser78.

Specificity

Human, Rat, Bovine, Guinea pig, Monkey

Purification

Protein A- affinity purified

Storage

Store at -20°C. Avoid multiple freeze/thaw cycles.

Application

IHC, WB (1:1000, colorimetric)

Ordering Information

Products

HSP27 (pSer78) Polyclonal Antibody

Size

50 µg
200 µg

Cat. No.

A-7427-050
A-7427-200



▲ **Fig. 1.** Western blot analysis of HSP27 (pSer78) Polyclonal Antibody: Lane 1 MW marker, Lane 2 phospho HSP27, Lane 3 non-phospho HSP27 and Lane 4 MDBK (weak reactivity).

This product is for research purposes only. Not intended for use in diagnostic procedures.