

HSPB2 Antibody [#C12418]

Catalog Number:	C12418
Concentration:	1mg/ml
Swiss-Prot No.:	Q16082
Other Names:	DMPK-binding protein; heat-shock protein beta-2; HSP27; MKBP
Storage/Stability:	Store at -20°C/1 year
Form of Antibody:	Rabbit IgG in phosphate buffered saline (without Mg ²⁺ and Ca ²⁺), pH 7.4, 150mM NaCl, 0.02% sodium azide and 50% glycerol.
Immunogen:	The antiserum was produced against synthesized peptide derived from internal of human HSPB2.
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	HSPB2 antibody detects endogenous levels of total HSPB2 protein.
Reactivity:	Human, Mouse, Rat
Applications:	WB: 1:500~1:1000 IF: 1:500~1:1000 ELISA: 1:40000
References:	Iwaki A., Genomics 45:386-394(1997). Suzuki A., J. Cell Biol. 140:1113-1124(1998). Elisei R., J. Endocrinol. Invest. 16:533-540(1993).