

HSP20 Antibody [#B0488]

Catalog Number:	B0488
Concentration:	1mg/ml
Swiss-Prot No.:	O14558
Other Names:	Heat-shock 20 kDa like-protein p20; HSPB6
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 HSP20.
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	HSP20 antibody detects endogenous levels of total HSP20 protein.
Reactivity:	Human (Identities = 100%, Positives = 100%); Mouse (Identities = 100%, Positives = 100%); Rat (Identities = 100%, Positives = 100%)
Applications:	IHC: 1:50~1:100 ELISA: 1:5000
References:	Ota T., Nat. Genet. 36:40-45(2004). Grimwood J., Nature 428:529-535(2004). Kato K., J. Biol. Chem. 269:15302-15309(1994).