

MRP7 Antibody [#C16727]

Catalog Number:	C16727
Concentration:	1mg/ml
Swiss-Prot No.:	Q5T3U5
Other Names:	ATP-binding cassette sub- family C member 10; MRP7; multidrug resistance-associated 7; SIMRP7
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 MRP7.
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	MRP7 antibody detects endogenous levels of total MRP7 protein.
Reactivity:	Human
Applications:	WB: 1:500~1:1000 ELISA: 1:10000
References:	Kao H.-H., J. Biomed. Sci. 10:98-110(2003). Ota T., Nat. Genet. 36:40-45(2004). Mungall A.J., Nature 425:805-811(2003).