

Caspase 10 Antibody [#C0135]

Catalog Number:	C0135
Amount:	50µg/50µl, 100µg/100µl, 200µg/200µl
Swiss-Prot No. :	Q92851
All Names:	ALPS2; CASP10B; FLICE2, Caspase 10 isoform a preproprotein, Caspase 10 isoform b preproprotein, Caspase 10 isoform c preproprotein, Caspase 10 isoform d preproprotein, FADD-LIKE ICE 2, FADD-like ICE2, Fas-associated death domain protein, ICE-like apoptotic protease 4, MCH4 CASPASE 10, ISOFORM B, apoptotic protease MCH-4, caspase 10, apoptosis-related cysteine protease, interleukin-1B-converting enzyme 2
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.
Storage/Stability:	Store at -20°C/1 year
Immunogen:	The antiserum was produced against synthesized peptide derived from human Caspase 10.
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity/Sensitivity	Caspase 10 antibody detects endogenous levels of total Caspase 10 protein.
Reactivity:	Human
Applications:	WB: 1:300~1:1000 IHC: 1:50~1:100 IF: 1:500~1:1000 ELISA: 1:10000
References:	Frank C. Kischkel, J. Biol. Chem., Nov 2001; 276: 46639 - 46646. Ingo H. Engels, Mol. Cell. Biol., Apr 2005; 25: 2808 - 2818. Daniela Sciaky, J. Immunol., Apr 2000; 164: 3806. Jin Wang, PNAS, Nov 2001; 98: 13884.