

CFLAR Antibody [#C12187]

Catalog Number:	C12187
Concentration:	1mg/ml
Swiss-Prot No.:	O15519
Other Names:	CASP8 and FADD-like apoptosis regulator , Cellular FLICE-like inhibitory protein, c-FLIP, Caspase-eight-related protein, Casper, Caspase-like apoptosis regulatory protein, CLARP, MACH-related inducer of toxicity, MRIT, Caspase homolog, CASH, Inhibitor of FLICE, I-FLICE, FADD-like antiapoptotic molecule 1, FLAME-1, Usurpin, CASP8 and FADD-like apoptosis regulator subunit p43, CASP8 and FADD-like apoptosis regulator subunit p12, CASH, CASP8AP1, CLARP, MRIT
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 CFLAR.
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	CFLAR antibody detects endogenous levels of total CFLAR protein.
Reactivity:	Human (Identities = 100%, Positives = 100%)
Applications:	IHC: 1:50~1:100 ELISA: 1:20000
References:	Shu H.-B., Immunity 6:751-763(1997). Han D.K.M., Proc. Natl. Acad. Sci. U.S.A. 94:11333-11338(1997). Srinivasula S.M., J. Biol. Chem. 272:18542-18545(1997).