

CKLF2 Antibody [#C15132]

Catalog Number:	C15132
Concentration:	1mg/ml
Swiss-Prot No.:	Q8TAZ6
Other Names:	chemokine-like factor 2; CKLF-like MARVEL transmembrane domain containing 2; CKLF2; CKLFSF2
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 CKLF2.
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	CKLF2 antibody detects endogenous levels of total CKLF2 protein.
Reactivity:	Human
Applications:	WB: 1:500~1:1000 ELISA: 1:40000
References:	Han W., Genomics 81:609-617(2003). Ota T., Nat. Genet. 36:40-45(2004). The MGC Project Team; Genome Res. 14:2121-2127(2004).