

CKLF1 Antibody [#C15131]

Catalog Number:	C15131
Concentration:	1mg/ml
Swiss-Prot No.:	Q8IZ96
Other Names:	CKLF-like MARVEL transmembrane domain-containing protein 1; Chemokine-like factor superfamily member 1
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 CKLF1.
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	CKLF1 antibody detects endogenous levels of total CKLF1 protein.
Reactivity:	Human (Identities = 100%, Positives = 100%)
Applications:	IF: 1:500~1:1000 ELISA: 1:60000
References:	Han W., Genomics 81:609-617(2003). Wang L., Submitted (NOV-2002) to the EMBL/GenBank/DDBJ databases.