

CLIP3 Antibody [#C14917]

Catalog Number:	C14917
Concentration:	1mg/ml
Swiss-Prot No.:	Q96DZ5
Other Names:	CLIP-170-related 59 kDa protein; CLIP3; CLR59; cytoplasmic linker protein 170-related 59 kDa protein
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 CLIP3.
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	CLIP3 antibody detects endogenous levels of total CLIP3 protein.
Reactivity:	Human, Mouse
Applications:	WB: 1:500~1:1000 IF: 1:500~1:1000 ELISA: 1:40000
References:	Perez F., J. Cell Biol. 156:631-642(2002). The MGC Project Team; Genome Res. 14:2121-2127(2004). The German cDNA consortium; Submitted (SEP-1999) to the EMBL/GenBank/DDBJ databases.