

RFPL4A Antibody [#C18339]

Catalog Number:	C18339
Concentration:	1mg/ml
Swiss-Prot No.:	A6NLU0
Other Names:	Ret finger protein-like 4A; RING finger protein 210; RFPL4A; RFPL4, RNF210
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 RFPL4A.
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	RFPL4A antibody detects endogenous levels of total RFPL4A protein.
Reactivity:	Human
Applications:	WB: 1:500~1:1000 ELISA: 1:5000
References:	Grimwood J., Nature 428:529-535(2004). Gruetter C., FEBS Lett. 580:99-106(2006).