

APOF Antibody [#C14533]

Catalog Number:	C14533
Concentration:	1mg/ml
Swiss-Prot No.:	Q13790
Other Names:	Apolipoprotein F; Apo-F; Lipid transfer inhibitor protein; LTIP
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 APOF.
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	APOF antibody detects endogenous levels of total APOF protein.
Reactivity:	Human (Identities = 100%, Positives = 100%)
Applications:	WB: 1:500~1:1000 ELISA: 1:40000
References:	Day J.R., Biochem. Biophys. Res. Commun. 203:1146-1151(1994). Wang X., J. Biol. Chem. 274:1814-1820(1999). The MGC Project Team; Genome Res. 14:2121-2127(2004).