

APOL5 Antibody [#C14543]

Catalog Number:	C14543
Concentration:	1mg/ml
Swiss-Prot No.:	Q9BWW9
Other Names:	Apolipoprotein-L5; Apolipoprotein L-V; ApoL-V
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 APOL5.
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	APOL5 antibody detects endogenous levels of total APOL5 protein.
Reactivity:	Human (Identities = 100%, Positives = 100%)
Applications:	IHC: 1:50~1:100 IF: 1:500~1:1000 ELISA: 1:40000
References:	Page N.M., Genomics 74:71-78(2001). Dunham I., Nature 402:489-495(1999).