

ACOT2 Antibody [#C14271]

Catalog Number:	C14271
Concentration:	1mg/ml
Swiss-Prot No.:	P49753
Other Names:	acyl-coenzyme A thioesterase 2; CTE-Ia; EC 3.1.2.2; peroxisomal acyl-coenzyme A thioester hydrolase 2a; peroxisomal long-chain acyl-coA thioesterase 2; ZAP128
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 ACOT2.
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	ACOT2 antibody detects endogenous levels of total ACOT2 protein.
Reactivity:	Human
Applications:	WB: 1:500~1:1000 ELISA: 1:40000
References:	Sherrington R., Nature 375:754-760(1995). Hunt M.C., FASEB J. 20:1855-1864(2006). Ota T., Nat. Genet. 36:40-45(2004).