

ACOT8 Antibody [#C14273]

Catalog Number:	C14273
Concentration:	1mg/ml
Swiss-Prot No.:	O14734
Other Names:	ACTE-III; ACTEIII; acyl-CoA thioesterase 8; acyl-coenzyme A thioesterase 8; choloyl-coenzyme A thioesterase; EC 3.1.2.27; HIV-Nef-associated acyl coA thioesterase; peroxisomal acyl-coenzyme A thioester hydrolase 1; peroxisomal long-chain acyl-coA thioesterase 1; PTE-1; PTE-2; TE; thioesterase II
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 ACOT8.
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	ACOT8 antibody detects endogenous levels of total ACOT8 protein.
Reactivity:	Human, Mouse, Rat
Applications:	WB: 1:500~1:1000 IF: 1:500~1:1000 ELISA: 1:40000
References:	Watanabe H., Biochem. Biophys. Res. Commun. 238:234-239(1997). Liu L.X., J. Biol. Chem. 272:13779-13785(1997). Jones J.M., J. Biol. Chem. 274:9216-9223(1999).