

AWAT1 Antibody [#C14257]

Catalog Number:	C14257
Concentration:	1mg/ml
Swiss-Prot No.:	Q58HT5
Other Names:	acyl-CoA wax alcohol acyltransferase 1; AWAT1; DGA2; diacyl-glycerol acyltransferase 2; diacylglycerol O-acyltransferase 2-like 3; EC 2.3.1.75; long-chain-alcohol O-fatty-acyltransferase 1
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 AWAT1.
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	AWAT1 antibody detects endogenous levels of total AWAT1 protein.
Reactivity:	Human
Applications:	WB: 1:500~1:1000 ELISA: 1:5000
References:	Turkish A.R., J. Biol. Chem. 280:14755-14764(2005). Ross M.T., Nature 434:325-337(2005). The MGC Project Team; Genome Res. 14:2121-2127(2004).