

ADA2L Antibody [#C10181]

Catalog Number:	C10181
Concentration:	1mg/ml
Swiss-Prot No.:	O75478
Other Names:	Transcriptional adapter 2-like, ADA2-like protein, TADA2L, KL04P
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 ADA2L.
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	ADA2L antibody detects endogenous levels of total ADA2L protein.
Reactivity:	Human, Mouse, Rat
Applications:	WB: 1:500~1:1000 ELISA: 1:10000
References:	Candau R., Mol. Cell. Biol. 16:593-602(1996). Ogryzko V.V., Cell 94:35-44(1998). Henriksson A., Mol. Cell. Biol. 17:3065-3073(1997).