

CA181 Antibody [#C19639]

Catalog Number:	C19639
Concentration:	1mg/ml
Swiss-Prot No.:	Q9NWK9
Other Names:	HIT-type zinc finger-containing protein C1orf181, HIT-type zinc finger-containing protein C1orf181, C1orf181
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 CA181.
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	CA181 antibody detects endogenous levels of total CA181 protein.
Reactivity:	Human, Mouse
Applications:	WB: 1:500~1:1000 ELISA: 1:5000
References:	Ota T., Nat. Genet. 36:40-45(2004). Gregory S.G., Nature 441:315-321(2006). Olsen J.V., Cell 127:635-648(2006).