

CAGE1 Antibody [#C14912]

Catalog Number:	C14912
Concentration:	1mg/ml
Swiss-Prot No.:	Q8TC20
Other Names:	CAGE-1; cancer-associated gene 1; cancer/testis antigen 3; CTAG3
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 CAGE1.
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	CAGE1 antibody detects endogenous levels of total CAGE1 protein.
Reactivity:	Human (Identities = 100%, Positives = 100%)
Applications:	WB: 1:500~1:1000 IHC: 1:50~1:100 ELISA: 1:40000
References:	Park S., Biochim. Biophys. Acta 1625:173-182(2003). Ota T., Nat. Genet. 36:40-45(2004). Mungall A.J., Nature 425:805-811(2003).