

BAGE3 Antibody [#C14676]

Catalog Number:	C14676
Concentration:	1mg/ml
Swiss-Prot No.:	Q86Y29
Other Names:	Cancer/testis antigen 2.3; CT2.3
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 BAGE3.
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	BAGE3 antibody detects endogenous levels of total BAGE3 protein.
Reactivity:	Human (Identities = 100%, Positives = 100%)
Applications:	IHC: 1:50~1:100 ELISA: 1:20000
References:	Ruault M., Eur. J. Hum. Genet. 10:833-840(2002). Ruault M., Genomics 81:391-399(2003).