

BCAS4 Antibody [#C14803]

Catalog Number:	C14803
Concentration:	1mg/ml
Swiss-Prot No.:	Q8TDM0
Other Names:	Breast carcinoma-amplified sequence 4; BCAS4
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 BCAS4.
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	BCAS4 antibody detects endogenous levels of total BCAS4 protein.
Reactivity:	Human
Applications:	WB: 1:500~1:1000 IHC: 1:50~1:100 ELISA: 1:20000
References:	Baerlund M., Genes Chromosomes Cancer 35:311-317(2002). Deloukas P., Nature 414:865-871(2001). Ota T., Nat. Genet. 36:40-45(2004).