

Breast Tumor Kinase Antibody [#B0783]

Catalog Number:	B0783
Concentration:	1mg/ml
Swiss-Prot No.:	Q13882
Other Names:	Breast tumor kinase, EC 2.7.10.2, PTK6, Tyrosine-protein kinase 6, Tyrosine-protein kinase brk
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 Breast Tumor Kinase.
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	Breast Tumor Kinase antibody detects endogenous levels of total Breast Tumor Kinase protein.
Reactivity:	Human (Identities = 100%, Positives = 100%)
Applications:	IF: 1:500~1:1000 ELISA: 1:5000
References:	Mitchell P.J., Oncogene 9:2383-2390(1994). Lee H.-Y., Mol. Cells 8:401-407(1998). Ota T., Nat. Genet. 36:40-45(2004).