

BST1 Antibody [#C14339]

Catalog Number:	C14339
Concentration:	1mg/ml
Swiss-Prot No.:	Q10588
Other Names:	ADP-ribosyl cyclase 2; bone marrow stromal antigen 1; BST-1; cADPr hydrolase 2; CD157 antigen; cyclic ADP-ribose hydrolase 2; EC 3.2.2.5
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 BST1.
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	BST1 antibody detects endogenous levels of total BST1 protein.
Reactivity:	Human, Mouse, Rat
Applications:	WB: 1:500~1:1000 ELISA: 1:5000
References:	Kaisho T., Proc. Natl. Acad. Sci. U.S.A. 91:5325-5329(1994). Kalnine N., Submitted (OCT-2004) to the EMBL/GenBank/DDBJ databases. Hillier L.W., Nature 434:724-731(2005).