

CBX6 Antibody [#C10234]

Catalog Number:	C10234
Concentration:	1mg/ml
Swiss-Prot No.:	O95503
Other Names:	Chromobox protein homolog 6, CBX6
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 CBX6.
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	CBX6 antibody detects endogenous levels of total CBX6 protein.
Reactivity:	Human, Mouse, Rat
Applications:	WB: 1:500~1:1000 ELISA: 1:20000
References:	Dunham I., Nature 402:489-495(1999). The MGC Project Team, Genome Res. 14:2121-2127(2004).