

PGBD1 Antibody [#C11505]

Catalog Number:	C11505
Concentration:	1mg/ml
Swiss-Prot No.:	Q96JS3
Other Names:	PiggyBac transposable element-derived protein 1, Cerebral protein 4, PGBD1, hucep-4
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 PGBD1.
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	PGBD1 antibody detects endogenous levels of total PGBD1 protein.
Reactivity:	Human
Applications:	WB: 1:500~1:1000 ELISA: 1:20000
References:	Yoshimoto M., Submitted (OCT-1996) to the EMBL/GenBank/DDBJ databases. Totoki Y., Submitted (APR-2005) to the EMBL/GenBank/DDBJ databases. Mungall A.J., Nature 425:805-811(2003).