

BRP44L Antibody [#C14781]

Catalog Number:	C14781
Concentration:	1mg/ml
Swiss-Prot No.:	Q9Y5U8
Other Names:	BR44L; brain protein 44-like
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 BRP44L.
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	BRP44L antibody detects endogenous levels of total BRP44L protein.
Reactivity:	Human, Mouse, Rat
Applications:	WB: 1:500~1:1000 IHC: 1:50~1:100 IF: 1:500~1:1000 ELISA: 1:20000
References:	Zhang Q.-H., Genome Res. 10:1546-1560(2000). Lai C.-H., Genome Res. 10:703-713(2000). Mungall A.J., Nature 425:805-811(2003).