

EPHB4 Antibody [#C10732]

Catalog Number:	C10732
Concentration:	1mg/ml
Swiss-Prot No.:	P54760
Other Names:	EC 2.7.1.112, EC 2.7.10.1, EPB4, Ephrin type-B receptor 4 precursor, HTK, Tyrosine-protein kinase receptor HTK
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 EPHB4.
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	EPHB4 antibody detects endogenous levels of total EPHB4 protein.
Reactivity:	Human, Mouse
Applications:	WB: 1:500~1:1000 ELISA: 1:20000
References:	Bennett B.D., J. Biol. Chem. 269:14211-14218(1994). Wilson M.D., Nucleic Acids Res. 29:1352-1365(2001). Zhang Y., Mol. Cell. Proteomics 4:1240-1250(2005).