

EPHA6 Antibody [#C11829]

Catalog Number:	C11829
Concentration:	1mg/ml
Swiss-Prot No.:	Q9UF33
Other Names:	Ephrin type-A receptor 6 [Precursor], EC 2.7.10.1, Tyrosine-protein kinase receptor EHK-2, EPH homology kinase 2, EPHA6, EHK2
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 EPHA6.
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	EPHA6 antibody detects endogenous levels of total EPHA6 protein.
Reactivity:	Human (Identities = 100%, Positives = 100%); Mouse (Identities = 100%, Positives = 100%); Rat (Identities = 100%, Positives = 100%)
Applications:	IHC: 1:50~1:100 IF: 1:500~1:1000 ELISA: 1:20000
References:	The German cDNA consortium; Submitted (JAN-2005) to the EMBL/GenBank/DDBJ databases. Muzny D.M., Nature 440:1194-1198(2006). Hafner C., Clin. Chem. 50:490-499(2004).