

## EPHA7 Antibody [#C11078]

<b>Catalog Number:</b>	C11078
<b>Concentration:</b>	1mg/ml
<b>Swiss-Prot No.:</b>	Q15375
<b>Other Names:</b>	Ephrin type-A receptor 7 [Precursor], EC 2.7.10.1, Tyrosine-protein kinase receptor EHK-3, EPH homology kinase 3, Receptor protein-tyrosine kinase HEK11, EHK3, HEK11
<b>Storage/Stability:</b>	Store at -20°C/1 year
<b>Form of Antibody:</b>	Rabbit IgG in phosphate buffered saline (without Mg <sup>2+</sup> and Ca <sup>2+</sup> ), pH 7.4, 150mM NaCl, 0.02% sodium azide and 50% glycerol.
<b>Immunogen:</b>	The antiserum was produced against synthesized peptide derived from internal of human EPHA7.
<b>Purification:</b>	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
<b>Specificity:</b>	EPHA7 antibody detects endogenous levels of total EPHA7 protein.
<b>Reactivity:</b>	Human, Mouse, Rat
<b>Applications:</b>	WB: 1:500~1:1000 ELISA: 1:40000
<b>References:</b>	Fox G.M., Oncogene 10:897-905(1995). Mungall A.J., Nature 425:805-811(2003). Sjoebloom T., Science 314:268-274(2006).