

## EFNA3 Antibody [#C10709]

<b>Catalog Number:</b>	C10709
<b>Concentration:</b>	1mg/ml
<b>Swiss-Prot No.:</b>	P52797
<b>Other Names:</b>	Ephrin-A3 [Precursor], EPH-related receptor tyrosine kinase ligand 3, LERK-3, EHK1 ligand, EHK1-L, EFL-2, EFNA3, EFL2, EPLG3, LERK3
<b>Storage/Stability:</b>	Store at -20°C/1 year
<b>Form of Antibody:</b>	Rabbit IgG in phosphate buffered saline (without $Mg^{2+}$ and $Ca^{2+}$ ), pH 7.4, 150mM NaCl, 0.02% sodium azide and 50% glycerol.
<b>Immunogen:</b>	The antiserum was produced against synthesized peptide derived from C-terminal of human EFNA3.
<b>Purification:</b>	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
<b>Specificity:</b>	EFNA3 antibody detects endogenous levels of total EFNA3 protein.
<b>Reactivity:</b>	Human, Mouse, Rat
<b>Applications:</b>	WB: 1:500~1:1000   ELISA: 1:10000
<b>References:</b>	Kozlosky C.J., Oncogene 10:299-306(1995). Davis S., Science 266:816-819(1994). Gregory S.G., Nature 441:315-321(2006).