

## FRK Antibody [#C10609]

<b>Catalog Number:</b>	C10609
<b>Concentration:</b>	1mg/ml
<b>Swiss-Prot No.:</b>	P42685
<b>Other Names:</b>	EC 2.7.1.112, EC 2.7.10.2, Nuclear tyrosine protein kinase RAK, Tyrosine-protein kinase FRK, fyn-related kinase
<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 internal of human FRK.
<b>Purification:</b>	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
<b>Specificity:</b>	FRK antibody detects endogenous levels of total FRK protein.
<b>Reactivity:</b>	Human, Mouse, Rat
<b>Applications:</b>	WB: 1:500~1:1000 IHC: 1:50~1:100 IF: 1:500~1:1000 ELISA: 1:5000
<b>References:</b>	Lee J., Gene 138:247-251(1994). Cance W.G., Cell Growth Differ. 5:1347-1355(1994). Zhang Y., Mol. Cell. Proteomics 4:1240-1250(2005).