

## FGR (Ab-412) Antibody [#B8105]

<b>Catalog Number:</b>	B8105
<b>Concentration:</b>	1mg/ml
<b>Swiss-Prot No.:</b>	P09769
<b>Other Names:</b>	C-FGR; EC 2.7.10.2; FGR protein; kinase Fgr; P55-FGR; Proto-oncogene tyrosine-protein kinase FGR; SRC2
<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 non-phosphopeptide derived from human FGR around the phosphorylation site of tyrosine 412 (D-E-Y <sup>P</sup> -N-P).
<b>Purification:</b>	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
<b>Specificity:</b>	FGR (Ab-412) antibody detects endogenous levels of total FGR protein.
<b>Reactivity:</b>	Human (Identities = 100%, Positives = 100%)
<b>Applications:</b>	WB: 1:500~1:1000 ELISA: 1:40000
<b>References:</b>	Katamine S.,Mol. Cell. Biol. 8:259-266(1988). Gregory S.G.,Nature 441:315-321(2006). The MGC Project Team;Genome Res. 14:2121-2127(2004).