

## SH3BGR Antibody [#C18718]

<b>Catalog Number:</b>	C18718
<b>Concentration:</b>	1mg/ml
<b>Swiss-Prot No.:</b>	P55822
<b>Other Names:</b>	SH3BGR protein; 21-glutamic acid-rich protein; 21-GARP
<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 SH3BGR.
<b>Purification:</b>	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
<b>Specificity:</b>	SH3BGR antibody detects endogenous levels of total SH3BGR protein.
<b>Reactivity:</b>	Human (Identities = 100%, Positives = 100%); Mouse (Identities = 91%, Positives = 100%); Rat (Identities = 91%, Positives = 100%)
<b>Applications:</b>	WB: 1:500~1:1000 ELISA: 1:10000
<b>References:</b>	Scartezzini P., Hum. Genet. 99:387-392(1997). Hattori M., Nature 405:311-319(2000). The MGC Project Team, Genome Res. 14:2121-2127(2004).