

## Gab2 (Ab-623) Antibody [#B8350]

<b>Catalog Number:</b>	B8350
<b>Concentration:</b>	1mg/ml
<b>Swiss-Prot No.:</b>	Q9UQC2
<b>Other Names:</b>	GABJ; GRB2-associated binding protein 2; KIAA0571; pp100
<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 non-phosphopeptide derived from human Gab2 around the phosphorylation site of serine 623 (P-S-S <sup>P</sup> -P-S).
<b>Purification:</b>	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
<b>Specificity:</b>	Gab2 (Ab-623) antibody detects endogenous levels of total Gab2 protein.
<b>Reactivity:</b>	Human (Identities = 100%, Positives = 100%); Mouse (Identities = 91%, Positives = 91%); Rat (Identities = 91%, Positives = 91%)
<b>Applications:</b>	WB: 1:500~1:1000 ELISA: 1:40000
<b>References:</b>	Nishida K., Blood 93:1809-1816(1999). Nagase T., DNA Res. 5:31-39(1998). Salomon A.R., Proc. Natl. Acad. Sci. U.S.A. 100:443-448(2003).