

GAB1 (Ab-659) Antibody [#B0788]

Catalog Number:	B0788
Concentration:	1mg/ml
Swiss-Prot No.:	Q13480
Other Names:	Growth factor receptor bound protein 2-associated protein 1
Storage/Stability:	Store at -20°C/1 year
Form of Antibody:	Rabbit IgG in phosphate buffered saline (without Mg^{2+} and Ca^{2+}), pH 7.4, 150mM NaCl, 0.02% sodium azide and 50% glycerol.
Immunogen:	The antiserum was produced against synthesized non-phosphopeptide derived from human GAB1 around the phosphorylation site of tyrosine 659 (V-D-Y ^P -V-V).
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	GAB1 (Ab-659) antibody detects endogenous levels of total GAB1 protein.
Reactivity:	Human, Mouse, Rat
Applications:	WB: 1:500~1:1000 IHC: 1:50~1:100 IF: 1:500~1:1000 ELISA: 1:10000
References:	Lisa S. Lock, Mol. Biol. Cell, Jun 2002; 13: 2132. Anders Kallin, J. Biol. Chem., Apr 2004; 279: 17897 - 17904. Mélanie Laramée, J. Biol. Chem., Mar 2007; 282: 7758 - 7769. Mariko Takahashi-Tezuka, Mol. Cell. Biol., Jul 1998; 18: 4109 - 4117.