

ZAK (Ab-165) Antibody [#B8200]

Catalog Number:	B8200
Concentration:	1mg/ml
Swiss-Prot No.:	Q9NYL2
Other Names:	AZK; EC 2.7.11.25; mixed lineage kinase-related kinase; MLTK; MRK-beta; sterile alpha motif and leucine zipper containing kinase AZK
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 ZAK around the phosphorylation site of serine 165 (H-M-S ^P -L-V).
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	ZAK (Ab-165) antibody detects endogenous levels of total ZAK protein.
Reactivity:	Human (Identities = 100%, Positives = 100%); Mouse (Identities = 100%, Positives = 100%)
Applications:	WB: 1:500~1:1000 ELISA: 1:40000
References:	Liu T.-C., Biochem. Biophys. Res. Commun. 274:811-816(2000). Bloem L.J., J. Mol. Cell. Cardiol. 33:1739-1750(2001). The MGC Project Team; Genome Res. 14:2121-2127(2004).