

TAK1 (Ab-184) Antibody [#B1123]

Catalog Number:	B1123
Concentration:	1mg/ml
Swiss-Prot No.:	O43318
Other Names:	EC 2.7.11.25, M3K7, MAP3K7, Mitogen-activated protein kinase kinase kinase 7, TGF-beta-activated kinase 1, Transforming growth factor-beta-activated kinase 1, kinase TAK1
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 TAK1 around the phosphorylation site of threonine 184 (I-Q-T ^P -H-M).
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	TAK1 (Ab-184) antibody detects endogenous levels of total TAK1 protein.
Reactivity:	Human, Mouse, Rat
Applications:	WB: 1:500~1:1000 ELISA: 1:10000
References:	Sakurai H., Biochem. Biophys. Res. Commun. 243:545-549(1998). Mungall A.J., Nature 425:805-811(2003). Shibuya H., Science 272:1179-1182(1996).