

MAP3K4 Antibody [#C11995]

Catalog Number:	C11995
Concentration:	1mg/ml
Swiss-Prot No.:	Q9Y6R4
Other Names:	EC 2.7.11.25, KIAA0213, M3K4, MAP three kinase 1, MAP3K4, MAPK/ERK kinase kinase 4, MAPKKK4, MEK kinase 4, MEKK 4, MTK1, Mitogen-activated protein kinase kinase kinase 4, kinase MEKK4
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 peptide derived from internal of human MAP3K4.
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	MAP3K4 antibody detects endogenous levels of total MAP3K4 protein.
Reactivity:	Human (Identities = 100%, Positives = 100%); Mouse (Identities = 100%, Positives = 100%)
Applications:	IHC: 1:50~1:100 ELISA: 1:10000
References:	Wong C.K., J. Biol. Chem. 279:39366-39373(2004). Mita H., Mol. Cell. Biol. 22:4544-4555(2002). Takekawa M., EMBO J. 16:4973-4982(1997).