

MEKKK 4 Antibody [#C10242]

Catalog Number:	C10242
Concentration:	1mg/ml
Swiss-Prot No.:	O95819
Other Names:	EC 2.7.11.1, HPK/GCK-like kinase HGK, M4K4, MAP4K4, MAPK/ERK kinase kinase 4, MEK kinase kinase 4, MEKKK 4, Nck interacting kinase, hepatocyte progenitor kinase-like/germinal center kinase-like kinase, mitogen-activated protein kinase kinase kinase 4
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 MEKKK 4.
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	MEKKK 4 antibody detects endogenous levels of total MEKKK 4 protein.
Reactivity:	Human, Mouse
Applications:	WB: 1:500~1:1000 IF: 1:500~1:1000 ELISA: 1:5000
References:	Yao Z., J. Biol. Chem. 274:2118-2125(1999). Wright J.H., Mol. Cell. Biol. 23:2068-2082(2003). Ishikawa K., DNA Res. 5:169-176(1998).