

MEK1 (Ab-286) Antibody [#B0681]

Catalog Number:	B0681
Concentration:	1mg/ml
Swiss-Prot No.:	Q02750
Other Names:	Dual specificity mitogen-activated protein kinase kinase 1, EC 2.7.12.2, ERK activator kinase 1, MAP kinase kinase 1, MAP2K1, MAPK/ERK kinase 1, MAPKK 1, MAPKK1, MEK1, MP2K1, PRKMK1, kinase MEK1
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 MEK1 around the phosphorylation site of threonine 286 (A-E-T ^P -P-P).
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	MEK1 (Ab-286) antibody detects endogenous levels of total MEK1 protein.
Reactivity:	Human, Mouse, Rat
Applications:	WB: 1:500~1:1000 IHC: 1:50~1:100 ELISA: 1:10000
References:	Shunichi Murakami, Genes & Dev., Feb 2004; 18: 290. Ronald A. Rimerman, J. Biol. Chem., May 2000; 275: 14736. Anne C. Marinovic, J. Biol. Chem., May 2002; 277: 16673 - 16681. Rohit Lal, AACR Meeting Abstracts, Apr 2006; 2006: 687.