

p38 MAPK (Ab-322) Antibody [#B0798]

Catalog Number:	B0798
Concentration:	1mg/ml
Swiss-Prot No.:	Q16539
Other Names:	CRK1, CSAID binding protein, CSBP, CSBP1, CSBP2, Cytokine suppressive anti- inflammatory drug binding protein, EC 2.7.11.24, MAP kinase MXI2, MAP kinase p38, MAPK14, MAX-interacting protein 2, MK14, MXI2, Mitogen-activated protein kinase 14, Mitogen-activated protein kinase p38, kinase p38-alpha
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 p38 MAPK around the phosphorylation site of tyreonine 322 (D-P-Y ^P -D-Q).
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	p38 MAPK (Ab-322) antibody detects endogenous levels of total p38 MAPK protein.
Reactivity:	Human, Mouse, Rat
Applications:	WB: 1:500~1:1000 IHC: 1:50~1:100 IF: 1:500~1:1000 ELISA: 1:5000
References:	Nasreen Khalil, J. Biol. Chem., Dec 2005; 280: 43000 - 43009. Monica Ambrose, AACR Meeting Abstracts, Apr 2005; 2005: 241. Hiroe Shiratsuchi, Am J Physiol Renal Physiol, Oct 2002; 283: 678.