

p70 S6 Kinase(Phospho-Thr389)Antibody [#8A0533]

Catalog Number:	8A0533
Concentration:	1mg/ml
Origin of Antibody:	Rabbit
Swiss-Prot No.:	P23443
Other Names:	KS6B1, P70-S6K, RPS6KB1, Ribosomal protein S6 kinase, Ribosomal protein S6 kinase, 70kDa, polypeptide 1, S6K, kinase p70S6K, p70-S6K
All Sites:	Human: Thr389/412; Mouse: Thr412; Rat:Thr412
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 phosphopeptide derived from human p70 S6 Kinase around the phosphorylation site of threonine 389/412 (G-F-T _P -Y-V).
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific phosphopeptide. The antibody against non-phosphopeptide was removed by chromatography using non-phosphopeptide corresponding to the phosphorylation site.
Specificity:	p70 S6 Kinase (Phospho-Thr389/412) antibody detects endogenous levels of p70 S6 Kinase only when phosphorylated at threonine 389/412.
Reactivity:	Human, Mouse, Rat.
Applications:	WB: 1:500~1:1000 IHC: 1:50~1:100 IF: 1:100~1:500 ELISA: 1:1000
References:	Grove J., Mol. Cell. Biol. 11:5541-5550(1991). Gout I., J. Biol. Chem. 273:30061-30064(1998). Olsen J.V., Cell 127:635-648(2006).