

p70 S6 Kinase (Ab-371) Antibody [#B0532]

Catalog Number:	B0532
Concentration:	1mg/ml
Swiss-Prot No.:	P23443
Other Names:	EC 2.7.11.1, KS6B1, P70-S6K, RPS6KB1, Ribosomal protein S6 kinase, Ribosomal protein S6 kinase, 70kDa, polypeptide 1, S6K, kinase p70S6K, p70-S6K
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 p70 S6 Kinase around the phosphorylation site of serine 371 (V-D-S ^P -P-D).
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	p70 S6 Kinase (Ab-371) antibody detects endogenous levels of total p70 S6 Kinase protein.
Reactivity:	Human, Mouse, Rat
Applications:	WB: 1:500~1:1000 IHC: 1:50~1:100 IF: 1:500~1:1000 ELISA: 1:20000
References:	O. Jameel Shah, J. Biol. Chem., Jan 2002; 277: 2525 - 2533. Chuanshu Huang, Cancer Res., Oct 2002; 62: 5689 - 5697. O. Jameel Shah, Am J Physiol Endocrinol Metab, Jul 2000; 279: E74 - E82. Isabelle Briaud, Diabetes, Apr 2003; 52: 974.