

S6K- α 2 Antibody [#C11074]

Catalog Number:	C11074
Concentration:	1mg/ml
Swiss-Prot No.:	Q15349
Other Names:	90 kDa ribosomal protein S6 kinase 2, EC 2.7.11.1, KS6A2, RPS6KA2, RSK-3, Ribosomal S6 kinase 3, Ribosomal protein S6 kinase alpha 2, S6K-alpha 2, kinase RSK3, p90-RSK 2, pp90RSK3
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 S6K- α 2.
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	S6K- α 2 antibody detects endogenous levels of total S6K- α 2 protein.
Reactivity:	Human (Identities = 100%, Positives = 100%); Mouse (Identities = 100%, Positives = 100%)
Applications:	IHC: 1:50~1:100 ELISA: 1:20000
References:	Zhao Y., Mol. Cell. Biol. 15:4353-4363(1995). Mungall A.J., Nature 425:805-811(2003). Denis N.J., Proteomics 7:868-874(2007).