

SGK (Ab-422) Antibody [#B0087]

Catalog Number:	B0087
Concentration:	1mg/ml
Swiss-Prot No.:	O00141
Other Names:	SGK1, SERUM/GLUCOCORTICOID-REGULATED KINASE
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 SGK around the phosphorylation site of serine 422 (G-F-S ^P -Y-A).
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	SGK (Ab-422) antibody detects endogenous levels of total SGK protein.
Reactivity:	Human, Mouse, Rat
Applications:	WB: 1:500~1:1000 IHC: 1:50~1:100 ELISA: 1:5000
References:	Nicola Perrotti, J. Biol. Chem., Mar 2001; 276: 9406 - 9412. Yun L. Ma, Learn. Mem., Mar 2006; 13: 114 - 118. Ge-Xin Wang, Am J Physiol Heart Circ Physiol, Aug 2004; 287: H533 - H544.