

## SGK(Phospho-Ser422)Antibody [#8A0087]

<b>Catalog Number:</b>	8A0087
<b>Concentration:</b>	1mg/ml
<b>Origin of Antibody:</b>	Rabbit
<b>Swiss-Prot No.:</b>	O00141
<b>Other Names:</b>	SGK1, SERUM/GLUCOCORTICOID-REGULATED KINASE
<b>All Sites:</b>	Human: Ser422; Mouse: Ser422; Rat: Ser421
<b>Storage/Stability:</b>	Store at -20°C/1 year
<b>Form of Antibody:</b>	Rabbit IgG in phosphate buffered saline (without $Mg^{2+}$ and $Ca^{2+}$ ), pH 7.4, 150mM NaCl, 0.02% sodium azide and 50% glycerol.
<b>Immunogen:</b>	The antiserum was produced against synthesized phosphopeptide derived from human SGK around the phosphorylation site of serine 422 (G-F-S <sup>P</sup> -Y-A).
<b>Purification:</b>	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.
<b>Specificity:</b>	SGK (phospho-Ser422) antibody detects endogenous levels of SGK only when phosphorylated at serine 422.
<b>Reactivity:</b>	Human, Mouse, Rat.
<b>Applications:</b>	WB: 1:500~1:1000 IHC: 1:50~1:100 ELISA: 1:40000
<b>References:</b>	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. Bao-Hong Zhang, J. Biol. Chem., Aug 2001; 276: 31620 - 31626.