

## LIMK2 (Ab-283) Antibody [#B8045]

<b>Catalog Number:</b>	B8045
<b>Concentration:</b>	1mg/ml
<b>Swiss-Prot No.:</b>	P53671
<b>Other Names:</b>	EC 2.7.11.1; kinase LIMK2; LIM domain kinase 2; LIMK-2
<b>Storage/Stability:</b>	Store at -20°C/1 year
<b>Form of Antibody:</b>	Rabbit IgG in phosphate buffered saline (without Mg <sup>2+</sup> and Ca <sup>2+</sup> ), pH 7.4, 150mM NaCl, 0.02% sodium azide and 50% glycerol.
<b>Immunogen:</b>	The antiserum was produced against synthesized non-phosphopeptide derived from human LIMK2 around the phosphorylation site of serine 283 (R-R-S <sup>P</sup> -L-R).
<b>Purification:</b>	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
<b>Specificity:</b>	LIMK2 (Ab-283) antibody detects endogenous levels of total LIMK2 protein.
<b>Reactivity:</b>	Human (Identities = 100%, Positives = 100%); Mouse (Identities = 100%, Positives = 100%); Rat (Identities = 100%, Positives = 100%)
<b>Applications:</b>	WB: 1:500~1:1000 ELISA: 1:5000
<b>References:</b>	Okano I., J. Biol. Chem. 270:31321-31330(1995). Osada H., Biochem. Biophys. Res. Commun. 229:582-589(1996). Collins J.E., Genome Biol. 5:RESEARCH84.1-RESEARCH84.11(2004).