

## LKB1 (Ab-334) Antibody [#B8127]

<b>Catalog Number:</b>	B8127
<b>Concentration:</b>	1mg/ml
<b>Swiss-Prot No.:</b>	Q15831
<b>Other Names:</b>	AMPKK; EC 2.7.11.1; kinase LKB1; PJS; serine/threonine kinase 11 gene (LKB1); Serine/threonine kinase KPM; Serine/threonine protein kinase 11; STK11
<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 non-phosphopeptide derived from human LKB1 around the phosphorylation site of serine 334 (W-R-S <sup>P</sup> -M-T).
<b>Purification:</b>	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
<b>Specificity:</b>	LKB1 (Ab-334) antibody detects endogenous levels of total LKB1 protein.
<b>Reactivity:</b>	Human (Identities = 100%, Positives = 100%); Mouse (Identities = 100%, Positives = 100%)
<b>Applications:</b>	WB: 1:500~1:1000 ELISA: 1:20000
<b>References:</b>	Nezu J.; ubmitted (AUG-1996) to the EMBL/GenBank/DDBJ databases. Jenne D.E., at. Genet. 18:38-43(1998). Bignell G.R., ancer Res. 58:1384-1386(1998).