

## QSK Antibody [#C18604]

<b>Catalog Number:</b>	C18604
<b>Concentration:</b>	1mg/ml
<b>Swiss-Prot No.:</b>	Q9Y2K2
<b>Other Names:</b>	Serine/threonine-protein kinase QSK EC 2.7.11.1; Salt-inducible kinase 3; SIK-3; SIK3 QSK
<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 peptide derived from internal of human QSK.
<b>Purification:</b>	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
<b>Specificity:</b>	QSK antibody detects endogenous levels of total QSK protein.
<b>Reactivity:</b>	Human (Identities = 100%, Positives = 100%); Mouse (Identities = 100%, Positives = 100%)
<b>Applications:</b>	IHC: 1:50~1:100 ELISA: 1:40000
<b>References:</b>	Lizcano J.M., EMBO J. 23:833-843(2004). Nagase T., DNA Res. 6:63-70(1999). Nakajima D., DNA Res. 9:99-106(2002).