

## DBF4 Antibody [#C11821]

<b>Catalog Number:</b>	C11821
<b>Concentration:</b>	1mg/ml
<b>Swiss-Prot No.:</b>	Q9UBU7
<b>Other Names:</b>	Protein DBF4 homolog A, Activator of S phase kinase, DBF4-type zinc finger-containing protein 1, Chiffon homolog A, DBF4, ASK, DBF4A, ZDBF1
<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 peptide derived from N-terminal of human DBF4.
<b>Purification:</b>	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
<b>Specificity:</b>	DBF4 antibody detects endogenous levels of total DBF4 protein.
<b>Reactivity:</b>	Human, Mouse
<b>Applications:</b>	WB: 1:500~1:1000    ELISA: 1:10000
<b>References:</b>	Kumagai H., Mol. Cell. Biol. 19:5083-5095(1999). Jiang W., EMBO J. 18:5703-5713(1999). Ota T., Nat. Genet. 36:40-45(2004).