

## CEP41 Antibody [#C15042]

<b>Catalog Number:</b>	C15042
<b>Concentration:</b>	1mg/ml
<b>Swiss-Prot No.:</b>	Q9BYV8
<b>Other Names:</b>	centrosomal protein of 41 kDa; CEP41; cep41 protein; testis-specific gene A14 protein
<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 internal of human CEP41.
<b>Purification:</b>	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
<b>Specificity:</b>	CEP41 antibody detects endogenous levels of total CEP41 protein.
<b>Reactivity:</b>	Human (Identities = 100%, Positives = 100%)
<b>Applications:</b>	IHC: 1:50~1:100   ELISA: 1:40000
<b>References:</b>	Yamada T., Gene 288:57-63(2002). Ota T., Nat. Genet. 36:40-45(2004). Andersen J.S., Nature 426:570-574(2003).