

## ACTR-1C Antibody [#C11368]

<b>Catalog Number:</b>	C11368
<b>Concentration:</b>	1mg/ml
<b>Swiss-Prot No.:</b>	Q8NER5
<b>Other Names:</b>	Activin receptor type-1C, EC 2.7.11.30, ACTR-1C, Activin receptor-like kinase 7, ALK-7, ACVR1C, ALK7
<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 ACTR-1C.
<b>Purification:</b>	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
<b>Specificity:</b>	ACTR-1C antibody detects endogenous levels of total ACTR-1C protein.
<b>Reactivity:</b>	Human (Identities = 100%, Positives = 100%); Mouse (Identities = 100%, Positives = 100%); Rat (Identities = 100%, Positives = 100%)
<b>Applications:</b>	IHC: 1:50~1:100 IF: 1:500~1:1000 ELISA: 1:20000
<b>References:</b>	Bondestam J., Cytogenet. Cell Genet. 95:157-162(2001). Roberts H.J., Biol. Reprod. 68:1719-1726(2003). Hillier L.W., Nature 434:724-731(2005).