

## ZIC1/2/3 Antibody [#C10137]

<b>Catalog Number:</b>	C10137
<b>Concentration:</b>	1mg/ml
<b>Swiss-Prot No.:</b>	Q15915
<b>Other Names:</b>	Zinc finger protein ZIC 1, Zinc finger protein of the cerebellum 1, ZIC1
<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 ZIC1/2/3.
<b>Purification:</b>	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
<b>Specificity:</b>	ZIC1/2/3 antibody detects endogenous levels of total ZIC1/2/3 protein.
<b>Reactivity:</b>	Human, Mouse
<b>Applications:</b>	WB: 1:500~1:1000 IHC: 1:50~1:100 ELISA: 1:20000
<b>References:</b>	Yokota N., Cancer Res. 56:377-383(1996).