

## AN32C Antibody [#C10101]

<b>Catalog Number:</b>	C10101
<b>Concentration:</b>	1mg/ml
<b>Swiss-Prot No.:</b>	O43423
<b>Other Names:</b>	Acidic leucine-rich nuclear phosphoprotein 32 family member C, Phosphoprotein 32-related protein 1, Tumorigenic protein pp32r1, ANP32C, PP32R1
<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 AN32C.
<b>Purification:</b>	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
<b>Specificity:</b>	AN32C antibody detects endogenous levels of total AN32C protein.
<b>Reactivity:</b>	Human
<b>Applications:</b>	WB: 1:500~1:1000 ELISA: 1:10000
<b>References:</b>	Kadkol S.S., Nat. Med. 5:275-279(1999). Kochevar G.J., Hum. Mutat. 23:546-551(2004). Matilla A., Cerebellum 4:7-18(2005).