

## ATXN2 Antibody [#C14578]

<b>Catalog Number:</b>	C14578
<b>Concentration:</b>	1mg/ml
<b>Swiss-Prot No.:</b>	Q99700
<b>Other Names:</b>	ATX2; ATXN2; SCA2; Spinocerebellar ataxia type 2 protein
<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 ATXN2.
<b>Purification:</b>	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
<b>Specificity:</b>	ATXN2 antibody detects endogenous levels of total ATXN2 protein.
<b>Reactivity:</b>	Human, Mouse
<b>Applications:</b>	WB: 1:500~1:1000 ELISA: 1:20000
<b>References:</b>	Pulst S.-M., Nat. Genet. 14:269-276(1996). Sanpei K., Nat. Genet. 14:277-284(1996). Sahba S., Genomics 47:359-364(1998).