

## Synuclein (Ab-129) Antibody [#B0582]

<b>Catalog Number:</b>	B0582
<b>Concentration:</b>	1mg/ml
<b>Swiss-Prot No.:</b>	P37840
<b>Other Names:</b>	NACP, Non-A beta component of AD amyloid, Non-A4 component of amyloid precursor, SYN, SYUA, alpha-synuclein, synuclein, alpha
<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 non-phosphopeptide derived from human Synuclein around the phosphorylation site of serine 129 (M-P-S <sup>P</sup> -E-E).
<b>Purification:</b>	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
<b>Specificity:</b>	Synuclein (Ab-129) antibody detects endogenous levels of total Synuclein protein.
<b>Reactivity:</b>	Human, Mouse
<b>Applications:</b>	WB: 1:500~1:1000 IHC: 1:50~1:100 ELISA: 1:5000
<b>References:</b>	Natalia Ninkina, Mol. Cell. Biol., Nov 2003; 23: 8233 - 8245. Julie Lotharius, J. Biol. Chem., Oct 2002; 277: 38884 - 38894. Deniz Kirik, PNAS, Mar 2003; 100: 2884. Zhong-Zong Pan, AACR Meeting Abstracts, Mar 2004; 2004: 1178.