

## $\alpha$ -Synuclein (Ab-133) Antibody [#B7235]

- Catalog Number:** B7235-1, B7235-2, B7235  
**Amount:** 50 $\mu$ g/50 $\mu$ l, 100 $\mu$ g/100 $\mu$ l, 200 $\mu$ g/200 $\mu$ l  
**Swiss-Prot No. :** P37840  
**All Names:** NACP, Non-A beta component of AD amyloid, Non-A4 component of amyloid precursor, SYN, SYUA, alpha-synuclein, synuclein alpha  
**Form of Antibody:** 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.  
**Storage/Stability:** Store at -20°C/1 year  
**Immunogen:** The antiserum was produced against synthesized non-phosphopeptide derived from human  $\alpha$ -Synuclein around the phosphorylation site of tyrosine 133 (E-G-Y-Q-D).  
**Purification:** The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.  
**Specificity/Sensitivity**  $\alpha$ -Synuclein (Ab-133) antibody detects endogenous levels of total  $\alpha$ -Synuclein protein.  
**Reactivity:** Human, Mouse, Rat  
**Applications:** WB: 1:500~1:1000 ELISA: 1:20000  
**References:** Yangfu Jiang, Cancer Res., Jul 2004; 64: 4539 - 4546.  
Alessandro Negro, FASEB J, Dec 2001; 10.1096.  
ALESSANDRA RECCHIA, FASEB J, Apr 2004; 18: 617 - 626.  
Christopher E. Ellis, J. Biol. Chem., Feb 2001; 276: 3879 - 3884.