

## Synuclein-pan Antibody [#C0336]

- Catalog Number:** C0336-1, C0336-2, C0336  
**Amount:** 50µg/50µl, 100µg/100µl, 200µg/200µl  
**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 peptide derived from human synuclein.
- Purification:** The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
- Specificity/Sensitivity** Synuclein-pan antibody detects endogenous levels of total Synuclein protein.
- Reactivity:** Human, Mouse, Rat
- Applications:** WB: 1:500~1:1000    IHC: 1:50~1:100    ELISA: 1:5000
- References:** Eirene Kontopoulou, Hum. Mol. Genet., Oct 2006; 15: 3012 - 3023.  
Heather Snyder, J. Biol. Chem., Mar 2005; 280: 7562 - 7569.  
Wanli W. Smith, J. Neurosci., Jun 2005; 25: 5544 - 5552.  
Yuxin Fan, Hum. Mol. Genet., Oct 2006; 15: 3002 - 3011.