



## Recombinant Human $\alpha$ -Synuclein, Biotin Labeled

Revision Number: 2.0

Last updated: 18March2024

Catalog #	AS-55581
Size	200 $\mu$ g

Product Name:	Recombinant Human $\alpha$ -Synuclein, Biotin Labeled
Catalog Number:	AS-55581
Lot Number:	See label on vial
Amount:	200 $\mu$ g
Source:	The recombinant human $\alpha$ -synuclein (GenBank Accession # NP_000336) was expressed and purified from <i>E. coli</i> and conjugated with biotin.
Purity:	Greater than 90% as determined by SDS-PAGE and mass spectrometry.
DOS:	See label on the vial
Storage:	Biotinylated human $\alpha$ -synuclein is supplied frozen at 1 mg/ml in 10 mM sodium phosphate buffer (pH=7.4). Store at -80 °C, and avoid repeated freeze-thaw cycles.
Instructions:	Parkinson's disease is predominantly a movement disorder resulting from degeneration of dopaminergic neurons in the substantia nigra. The cause of the disease is unknown, but Substantial evidence suggests that the aggregation of $\alpha$ -synuclein is a critical step in the etiology of Parkinson's disease (PD). $\alpha$ -Synuclein is an abundant brain protein of 140 residues that is present in high concentration at presynaptic terminals and is found in both soluble and membrane-associated fractions of the brain. Several possible functions have been suggested, and it appears to be involved in vesicle release and trafficking. In vitro incubation in the presence of certain amount of salt (i.e. 0.1M NaCl) with agitation, $\alpha$ -synuclein can forms fibrilous structure.

### Related Products

Product Name	Cat. #
EndoClearPlus Recombinant human $\alpha$ -synuclein	AS-56081
<a href="#">Sensolyte® Anti-<math>\alpha</math>-Synuclein (Human) ELISA Kit</a>	AS-55550-H
<a href="#">Sensolyte® Anti-<math>\alpha</math>-Synuclein (Rat) ELISA Kit</a>	AS-55550-R
EndoClear Recombinant human $\alpha$ - synuclein	AS-55555
<a href="#">Recombinant human <math>\alpha</math> - synuclein, HiLyte Fluor™488 labeled</a>	AS-55457
EndoClear Recombinant mouse $\alpha$ - synuclein	AS-56082
EndoClear Recombinant rat $\alpha$ - synuclein	AS-56083

### References:

1. Trojanowski, J. Q. & Lee, V. M. (2003) *Ann. N. Y. Acad. Sci.* **991**, 107-110.
2. Masliah, E., et al. (2000) *Science* **287**, 1265-1269.
3. Van Der, P. H, et al. (2000) *J. Neurosci.* **20**, 6021-6029.
4. Feany, M. B. & Bender, W. W. (2000) *Nature* **404**, 394-398.
5. Weinreb, P. H., et al. (1996) *Biochemistry* **35**, 13709-13715.

**For in vitro research use only.**