

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

Revision Number: 2.0

Last updated: 18March2024

<b>Catalog #</b>	<b>AS-55581</b>
<b>Size</b>	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 *E. coli* 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  $-80^{\circ}\text{C}$ , 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	<b>AS-56081</b>
<a href="#">Sensolyte® Anti-<math>\alpha</math>-Synuclein (Human) ELISA Kit</a>	<b>AS-55550-H</b>
<a href="#">Sensolyte® Anti-<math>\alpha</math>-Synuclein (Rat) ELISA Kit</a>	<b>AS-55550-R</b>
EndoClear Recombinant human $\alpha$ - synuclein	<b>AS-55555</b>
<a href="#">Recombinant human <math>\alpha</math> - synuclein, HiLyte Fluor™488 labeled</a>	<b>AS-55457</b>
EndoClear Recombinant mouse $\alpha$ - synuclein	<b>AS-56082</b>
EndoClear Recombinant rat $\alpha$ - synuclein	<b>AS-56083</b>

### 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.***