



Product Data Sheet

Product Name:	Recombinant Human α-Synuclein (1-140), Hilyte Fluor™ 488 Labeled
Catalog Number:	AS-55457
Lot Number:	See label on vial
Amount:	200 μ g
Source:	The recombinant human α -synuclein (1-140) (GenBank Accession # NP_000336) was expressed and purified from <i>E. coli</i> and conjugated with the fluorescence dye Hilyte Fluor™ 488.
Purity:	Greater than 90% as determined by SDS-PAGE.
Fluorescence:	Green fluorescence. Excitation/Emission wavelengths= 490nm/525nm
DOS:	See label on the vial
Storage:	Hilyte Fluor™ 488 labeled human α -synuclein is supplied frozen at 1 mg/ml in 10 mM sodium phosphate buffer (pH=7.0). Store at 2-4 °C for immediate use within 1 week or at -80 °C for long-term storage. Keep in dark and avoid repeated freeze-thaw cycles.

Description: 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 α -synuclein is a critical step in the etiology of Parkinson's disease (PD). α -Synuclein is an abundant brain protein of 140 residues that 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.

Related Products

Product Name	Cat. #
EndoClearPlus Recombinant human α -synuclein (1-140)	AS-56081
Recombinant Human β - Synuclein (1 - 134)	AS-55458
Sensolyte® Anti-α-Synuclein (Human) ELISA Kit	AS-55550
EndoClear Recombinant human α - synuclein (1 - 140)	AS-55555
Recombinant human α - synuclein (1 - 140), biotin labeled	AS-55581

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.