

## EndoClearPlus Human Recombinant alpha-synuclein protein

Revision Number: 2.0 Last updated: 18March2024

Catalog #	AS-56081
Size	100 μg, 500 μg

Product Name: EndoClearPlus Human Recombinant alpha-synuclein protein (essentially endotoxin free)

Catalog Number: AS-56081

Lot Number: See label on the vial

Amount/size: 100 μg, 500 μg

Source: The sequence (Accession # NP\_000336) corresponding to the full-length human asynuclein protein plus an affinity tag was expressed in E. coli. The recombinant a-synuclein protein was purified from bacterial lysate using proprietary method, including tag cleavage. The resulting protein contains an additional Gly at the N-term of the a-synuclein sequence. The calculated molecular mass of the recombinant a-synuclein protein is ~14.5 kDa.

Activity: α-synuclein was assayed for the fibrillation properties using Thioflavin T assay.

Applications: Fibrillation, oligomerization, cell toxicity studies.

Purity: Greater than 90% as determined by SDS-PAGE.

Endotoxin: Less than 0.001 EU per μg/protein as tested by LAL method.

Storage: The purified  $\alpha$ -synuclein is supplied as lyophilized powder nearly salt free. Store at 2-4  $^{\circ}$ C for

immediate use within 1-2 weeks or at -80 °C for up to 12 months.

## Instructions:

A-synuclein is a major component of Lewy bodies in the affected neurons in Parkinson's disease.<sup>1-4</sup> This protein consists of a conserved degenerative amino-terminal domain and an acidic carboxyl-terminal with higher sequence divergence.<sup>1-4</sup> A-synuclein is predominantly expressed in brain: specifically in cerebellum, thalamus, neocortex, hippocampus, and striatum regions.<sup>1-2</sup> Other tissues express α-synuclein at very low levels.<sup>1-4</sup> The physiological role of α-synuclein is not yet well understood. However, the presence of imperfect KTKEGV lipid interacting repeats suggests that it may be involved in synaptic vesicle homeostasis.<sup>3</sup>

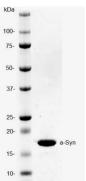


Figure 1. SDS-PAGE of human  $\alpha$ -synuclein. Purified protein was loaded onto 10-20% Tris-HCl gel at 3  $\mu$ g/well.



Figure 2. Western Immunoblot of human  $\alpha$ -synuclein. Purified  $\alpha$ -synuclein protein was probed with anti- $\alpha$ -synuclein specific antibody.

## **Related Products**

Product Name	Cat. #
Recombinant human a - synuclein, HiLyte Fluor™488 labeled	AS-55457
SensoLyte® Anti-a-Synuclein (Human) ELISA Kit	AS-55550-H
SensoLyte® Anti-a-Synuclein (Rat) ELISA Kit	AS-55550-R
EndoClear Recombinant human a - synuclein	AS-55555
Recombinant human a - synuclein, biotin labeled	AS-55581
EndoClear Recombinant mouse a - synuclein	AS-56082
EndoClear Recombinant rat a - synuclein	AS-56083

## **References:**

- 1. Rivers, R. et al. Protein Science 17, 887-898 (2008).
- 2. Bruening, W. et al. Cancer. 88, 9, 2154-2163 (2000).
- 3. George M. J. Genome Biology 3, 1, 1-6 (2001).
- 4. Latawiec, D. et al. PloS ONE 5, 2, 1-8 (2010).

For Research Use Only!