



## Rabbit (polyclonal) Anti- $\alpha$ -Synuclein

### PRODUCT ANALYSIS SHEET

<b>Catalog Number:</b>	AHB0401
<b>Lot Number:</b>	See product label
<b>Quantity:</b>	0.1 mL
<b>Form of Antibody:</b>	Purified immunoglobulin in phosphate buffered saline.
<b>Preservation:</b>	10 mM sodium azide (Caution: sodium azide is a poisonous and hazardous substance. Handle with care and dispose of properly.)
<b>Purification:</b>	Purified by caprylic acid and ammonium sulfate precipitation.
<b>Immunogen:</b>	A synthetic peptide, corresponding to amino acid residues 117-131 (PVDPDNEAVEMPSEE) of human $\alpha$ -synuclein, conjugated to keyhole limpet hemocyanin via an N-terminal cysteine residue.
<b>Specificity:</b>	<p><math>\alpha</math>-Synuclein (Genbank accession number 2136215; known alternatively as NACP, PARK1, and PD1 [Parkinson disease, familial 1]) is a member of the synuclein family of proteins that also includes <math>\beta</math>-synuclein and <math>\gamma</math>-synuclein. The gene for human <math>\alpha</math>-synuclein maps to chromosome 4 (cytogenetic band 4q21). Alternative splicing produces at least three variants. The major splice variant is a protein with <math>M_r=14.5</math> kDa, comprised of 140 amino acid residues. <math>\alpha</math>-Synuclein's N-terminus contains 11 positively charged amino acid repeat sequences that have the potential to form amphipathic alpha helices. <math>\alpha</math>-Synuclein's central region is hydrophobic, while its C-terminal region is highly acidic and contains several <math>\square</math>mygdal residues. In physiological solutions, <math>\alpha</math>-synuclein is found as three major forms: (1) soluble, natively unfolded monomers; (2) structured, beta pleated sheet-containing protofibrils; and (3) amyloid fibrils. Residues 71 and 81 of <math>\alpha</math>-synuclein are essential for filament assembly. <math>\alpha</math>-Synuclein undergoes several posttranslational modifications, including phosphorylation at serine 87, serine 129, and tyrosine 125, O-glycosylation, and nitrosylation.</p> <p><math>\alpha</math>-Synuclein is highly expressed in several areas of the brain, including the substantia nigra, thalamus, hippocampus, <math>\square</math>mygdale, corpus callosum, and caudate nucleus, appearing both in neurons where it concentrates at synapses, and in glial cells. Most other tissues express <math>\alpha</math>-synuclein at relatively low levels.</p> <p>Several neurodegenerative diseases are associated with the accumulation of <math>\alpha</math>-synuclein, including: (1) Parkinson's Disease, (2) Dementia with Lewy Bodies (DLB), (3) Lewy Body Variant of Alzheimer's disease (LBVAD), (4) Multiple Systems Atrophy (MSA), and (5) Neurodegeneration with Brain Iron Accumulation type-1 (NBIA-1). <math>\alpha</math>-Synuclein also appears as a minor protein within the neurofibrillary tangles of Alzheimer's disease.</p>
<b>Species Reactivity:</b>	Human. Slight reactivity observed with $\alpha$ -synuclein from rat brain. Does not react with mouse. Other species have not been tested.

**This product is for research use only. Not for use in diagnostic procedures.**

[www.invitrogen.com](http://www.invitrogen.com)

Invitrogen Corporation • 542 Flynn Rd • Camarillo • CA 93012 • Tel: 800.955.6288 • E-mail: [techsupport@invitrogen.com](mailto:techsupport@invitrogen.com)

PI AHB0401

(Rev 10/08) DCC-08-1089

**Important Licensing Information** - These products may be covered by one or more Limited Use Label Licenses (see the Invitrogen Catalog or our website, [www.invitrogen.com](http://www.invitrogen.com)). By use of these products you accept the terms and conditions of all applicable Limited Use Label Licenses. Unless otherwise indicated, these products are for research use only and are not intended for human or animal diagnostic, therapeutic or commercial use.

<b>Applications:</b>	This antibody is suitable for use in Western blotting and immunohistochemistry with formalin-fixed, paraffin-embedded tissue sections.
<b>Suggested Working Dilutions:</b>	In Western blotting, a 1:1,000 to 1:20,000 dilution is recommended. The optimal concentration should be determined for each specific application.
<b>Recommended Positive Control:</b>	Human brain temporal cortex exhibiting Dementia with Lewy Body pathology.
<b>Storage:</b>	Store at -20°C. Upon initial thawing, apportion into working aliquots and store at -20°C. Avoid repeated freeze-thaw cycles to prevent denaturing the antibody.
<b>Expiration Date:</b>	Expires one year from date of receipt when stored as instructed.
<b>References:</b>	<p>Ancolio, K. et al. (2000) <math>\alpha</math>-Synuclein and the Parkinson's disease-related mutant Ala53Thr-alpha-synuclein do not undergo proteasomal degradation in HEK293 and neuronal cell. <i>Neurosci. Lett.</i> 285:79-82.</p> <p>Bruening, W., et al. (2000) Synucleins are expressed in the majority of breast and ovarian carcinomas and in preneoplastic lesions of the ovary. <i>Cancer</i> 88(9):2154-2163.</p> <p>Dawson, T.M. and V.L. Dawson (2003) Molecular pathways of neurodegeneration in Parkinson's disease. <i>Science</i> 302:819-822.</p> <p>Dev K.K., et al. (2003) Part II: alpha-synuclein and its molecular pathophysiological role in neurodegenerative disease. <i>Neuropharmacology</i> 45(1):14-44.</p> <p>Glasson, B.I. and V. M.-Y. Lee (2003) Are ubiquitination pathways central to Parkinson's disease? <i>Cell</i> 114:1-8.</p> <p>Gorner, K., et al. (2003) Differential effects of Parkinson's disease-associated mutations on stability and folding of DJ-1. <i>J. Biol. Chem.</i> 279:6943-6951.</p> <p>Kaplan, B., et al. (2003) Alpha-synuclein: its biological function and role in neurodegenerative diseases. <i>J. Mol. Neurosci.</i> 20(2):83-92.</p> <p>Lim K.L., et al. (2003) The cast of molecular characters in Parkinson's disease: felons, conspirators, and suspects. <i>Ann. N.Y. Acad. Sci.</i> 991:80-92.</p> <p>McLean, P.J., et al. (2000) Membrane association and protein conformation of <math>\alpha</math>-synuclein in intact neurons. <i>J. Biol. Chem.</i> 275 (12):8812-8816.</p> <p>Okochi, M., et al. (2000) Constitutive phosphorylation of the Parkinson's disease associated <math>\alpha</math>-synuclein. <i>J. Biol. Chem.</i> 275(1):390-397.</p> <p>Richfield, E.K., et al. (2002) Behavioral and neurochemical effects of wild-type and mutated human <math>\alpha</math>-synuclein in transgenic mice. <i>Exp. Neurol.</i> 175:35-48.</p> <p>Singleton, A.B., et al. (2003) <math>\alpha</math>-Synuclein locus triplication causes Parkinson's disease. <i>Science</i> 308:841.</p> <p>Trojanowski, J.Q. and V.M. Lee (2003) Parkinson's disease and related alpha-synucleinopathies are brain amyloidoses. <i>Ann. N. Y. Acad. Sci.</i> 991:107-110.</p>

**This product is for research use only. Not for use in diagnostic procedures.**

[www.invitrogen.com](http://www.invitrogen.com)

Invitrogen Corporation • 542 Flynn Rd • Camarillo • CA 93012 • Tel: 800.955.6288 • E-mail: [techsupport@invitrogen.com](mailto:techsupport@invitrogen.com)

PI AHB0401

(Rev 10/08) DCC-08-1089

**Important Licensing Information** - These products may be covered by one or more Limited Use Label Licenses (see the Invitrogen Catalog or our website, [www.invitrogen.com](http://www.invitrogen.com)). By use of these products you accept the terms and conditions of all applicable Limited Use Label Licenses. Unless otherwise indicated, these products are for research use only and are not intended for human or animal diagnostic, therapeutic or commercial use.