

α -Synuclein polyclonal antibody

α -Synuclein accumulates in Lewy bodies, which are intraneuronal cytoplasmic inclusions present in the brains of sporadic Parkinson's disease patients, and is implicated in the pathogenesis of Parkinson's disease and related neurodegenerative disorders. α -Synuclein appears to associate with other proteins that aggregate and is found in β -amyloid plaques and neuritic tangles in Alzheimer's disease.

This antibody is covered by our [Worry-Free Guarantee](#).

Citations: 4

[View Online »](#)

Ordering Information

[Order Online »](#)

BML-SA3400-0025	25 μ l
BML-SA3400-0100	100 μ l

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Use/Stability	Vial contents, Storage and Use: Vial contains an immunoglobulin preparation, partially purified by chromatography from whole serum, suspended in phosphate-buffered saline (pH 7.4) containing 0.01M sodium azide. Store unopened vial at -20°C until required for use. AVOID REPEATED FREEZE-THAW CYCLES. Aliquot undiluted antibody into smaller volumes (not less than 10µL) prior to freezing if appropriate. The use of high quality 'antiserum-grade' plastic or glass vials is recommended. Store diluted antibody at 2-4°C (do not freeze) and use within 1 month. Dilute to working strength with 50mM PBS (pH 7.4) containing 1.5% sodium chloride and 1% normal goat serum (if a goat anti-rabbit IgG linker antibody is to be used).
Handling	Avoid freeze/thaw cycles. After opening, prepare aliquots and store at -20°C.
Long Term Storage	-20°C
Shipping	Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name	NACP, Non-A β component of senile plaques
Application	IHC, WB
Application Notes	<p>Western blot: The antibody has been characterized by single dimension SDS-PAGE using recombinant α- and β-synucleins as well as rat, mouse and human brain lysates with and without peptide blockade. Studies with the recombinant proteins demonstrated the absolute specificity of the antibody for α-synuclein. The antibody gives a strong band at a relative molecular weight of approximately 15kDa with human brain lysate which is wholly abolished by preincubation with the immunizing peptide (Prod. No. BML-SP9340).</p> <p>Immunohistochemistry: The antibody has been tested on de-waxed tissue sections of human brain fixed in neutral buffered formaldehyde. A standard protocol should be followed.</p>
Crossreactivity	Does not cross-react with mouse α -synuclein.
Formulation	Liquid. In PBS, pH 7.4, containing 10mM sodium azide.
GenBank ID	2136215
Host	Rabbit

Immunogen	Synthetic peptide corresponding to [Cys] + aa 117-131 of human α -synuclein conjugated to KLH. The sequence is conserved in rat α -synuclein except for two substitutions, namely Asp121Ser and Asn122Ser.
Purity Detail	Purified by caprylic acid and ammonium sulphate precipitation.
Recommendation Dilutions/Conditions	Western blot (1:1,000-1:20,000)Suggested dilutions/conditions may not be available for all applications.Optimal conditions must be determined individually for each application.
Source	Purified from rabbit serum.
Species Reactivity	Human, Rat
Specificity	This antibody has been shown to react strongly with human brain-derived α -synuclein. The antibody also reacts with rat brain-derived α -synuclein but reactivity appears substantially less than with human.
UniProt ID	P37840
Worry-free Guarantee	This antibody is covered by our Worry-Free Guarantee .



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
[info-
usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
[info-
eu@enzolifesciences.com](mailto:info-eu@enzolifesciences.com)

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
[info-
be@enzolifesciences.com](mailto:info-be@enzolifesciences.com)

France
Phone: +33 472 440 655
[info-
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany
Phone: +49 7621 5500 526
[info-
de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
[info-
uk@enzolifesciences.com](mailto:info-uk@enzolifesciences.com)