

# $\alpha$ -Synuclein (human) monoclonal antibody (5C2)

$\alpha$ -synuclein, an acidic neuronal protein of 140 aa, has been suggested to be implicated in the pathogenesis of Parkinson's disease and related neurodegenerative disorders, and to be an important regulatory component of vesicular transport in neuronal cells. The NAC (non-A $\beta$  component of amyloid plaques) domain (aa 61-95) of  $\alpha$ -synuclein is an highly amyloidogenic peptide which was first identified associated with senile plaques in the Alzheimer's disease brain.

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

Citations: 2

[View Online »](#)

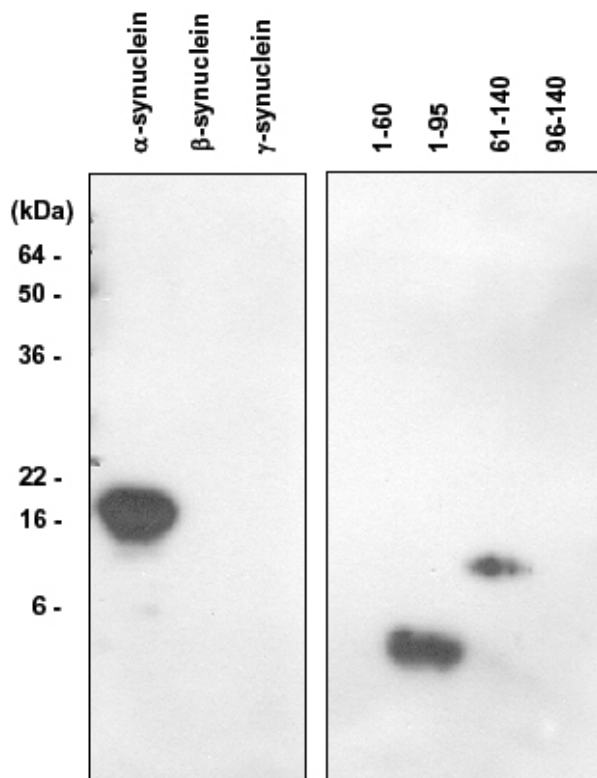
## Ordering Information

[Order Online »](#)

ALX-804-656-R100	100 $\mu$ l
------------------	-------------

## Manuals, SDS & CofA

[View Online »](#)



**Figure:** Western blot analysis. The recombinant human synuclein family ( $\alpha$ -,  $\beta$ - and  $\gamma$ -) and  $\alpha$ -synuclein domains (1-60, 1-95, 61-140 and 96-140) proteins were resolved by SDS-PAGE, transferred to PVDF membrane and probed with MAb to  $\alpha$ -Synuclein (human) (5C2) (Prod. No. ALX-804-656) (1:1'000). Proteins were visualized using a goat anti-mouse secondary antibody conjugated to HRP and an ECL detection system.

# Handling & Storage

Use/Stability	Stable for one month when stored at 4°C.
Handling	Avoid freeze/thaw cycles.
Short Term Storage	+4°C
Long Term Storage	-20°C
Shipping	Blue Ice

## Regulatory Status

RUO - Research Use Only

# Product Details

Application	ELISA, WB
Clone	5C2
Formulation	Liquid. In PBS, pH 7.4, containing 0.02% sodium azide and 10% glycerol.
Host	Mouse
Immunogen	Recombinant human α-synuclein (aa 1-140).
Isotype	IgG1
Purity Detail	Protein G-affinity purified.
Recommendation Dilutions/Conditions	Western Blot (1:500-1:2,000)Suggested dilutions/conditions may not be available for all applications.Optimal conditions must be determined individually for each application.
Species Reactivity	Human
Specificity	Recognizes the NAC domain (aa 61-95) of human α-synuclein.
UniProt ID	P37840

Worry-free Guarantee

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

Last modified: May 29, 2024



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)