

# $\alpha$ -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.

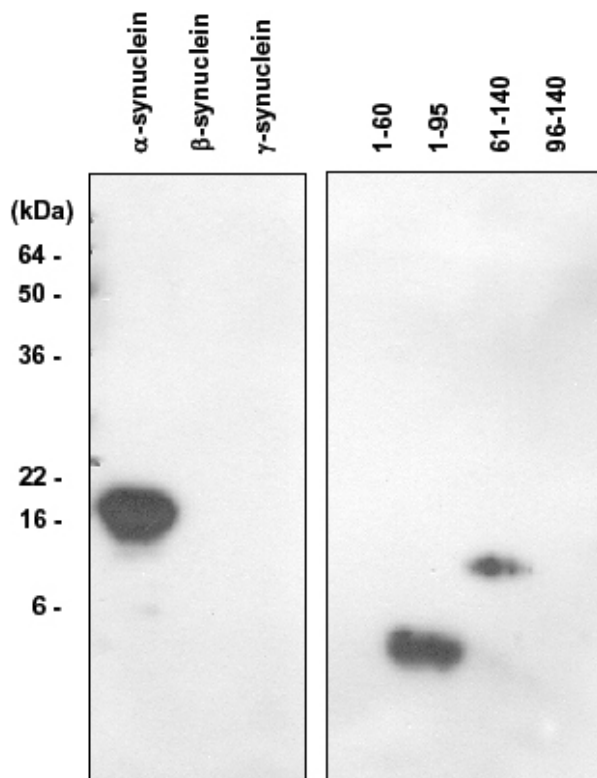
This antibody is covered by our [Worry-Free Guarantee](#).

Citations: 2

[View Online »](#)

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

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

**Regulatory Status** RUO - Research Use Only

## Product Details

<b>Application</b>	ELISA, WB
<b>Clone</b>	5C2
<b>Formulation</b>	Liquid. In PBS, pH 7.4, containing 0.02% sodium azide and 10% glycerol.
<b>Host</b>	Mouse
<b>Immunogen</b>	Recombinant human $\alpha$ -synuclein (aa 1-140).
<b>Isotype</b>	IgG1
<b>Purity Detail</b>	Protein G-affinity purified.
<b>Recommendation Dilutions/Conditions</b>	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.
<b>Species Reactivity</b>	Human
<b>Specificity</b>	Recognizes the NAC domain (aa 61-95) of human $\alpha$ -synuclein.
<b>UniProt ID</b>	P37840

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)