

α -Synuclein (human) monoclonal antibody (5C2)

α -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 β component of amyloid plaques) domain (aa 61-95) of α -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 μ l
------------------	-------------

Manuals, SDS & CofA

[View Online »](#)

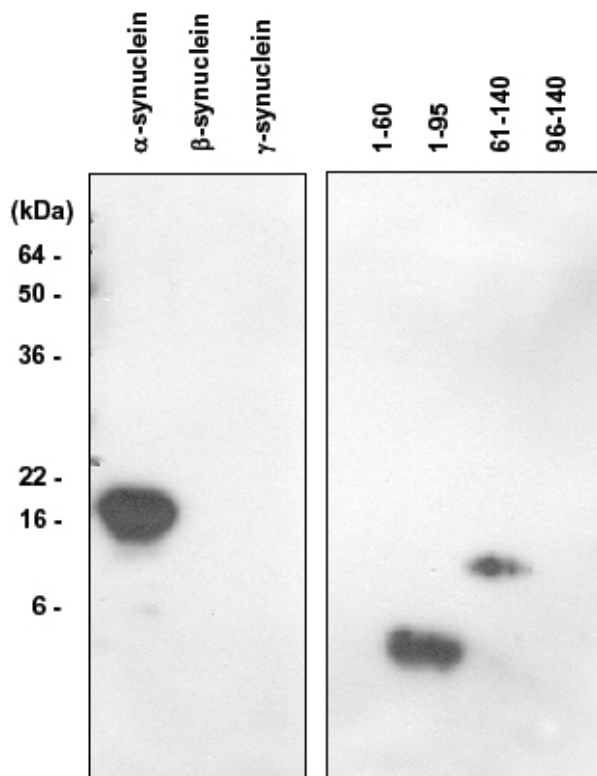


Figure: Western blot analysis. The recombinant human synuclein family (α -, β - and γ -) and α -synuclein domains (1-60, 1-95, 61-140 and 96-140) proteins were resolved by SDS-PAGE, transferred to PVDF membrane and probed with MAb to α -Synuclein (human) (5C2) (Prod. No. ALX-804-656) (1:1'000). Proteins 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 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



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
info-usa@enzolifesciences.com

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

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

France
Phone: +33 472 440 655
info-fr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info-de@enzolifesciences.com

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

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