

# $\alpha$ -Synuclein recombinant monoclonal antibody (017)

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

## Ordering Information

[Order Online »](#)

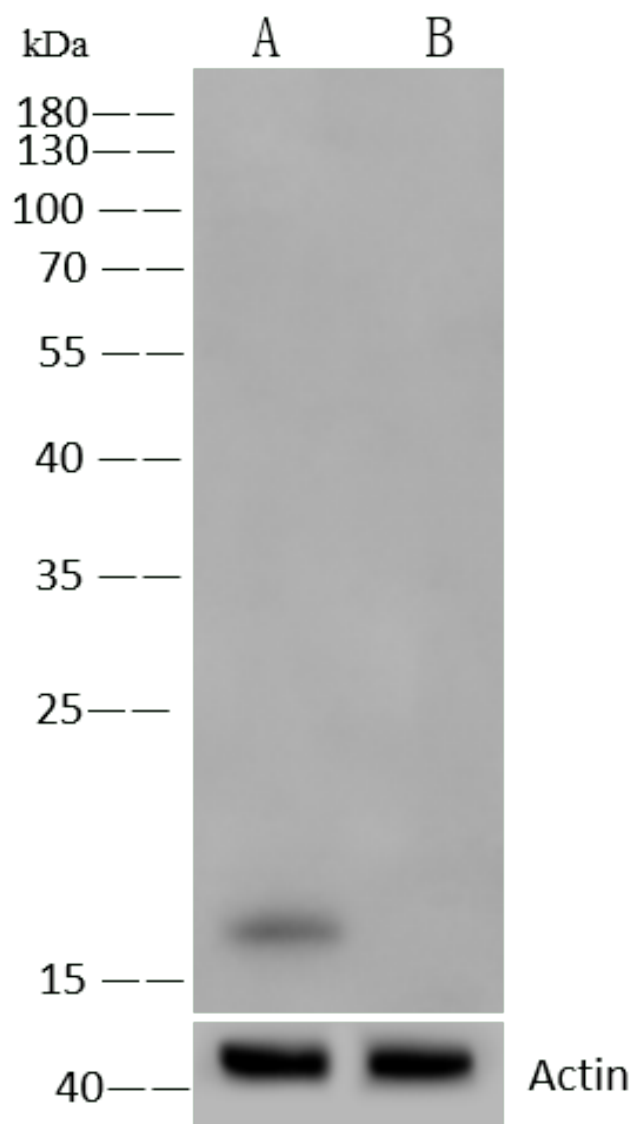
ENZ-ABS989-0100	100μl
-----------------	-------

## Manuals, SDS & CofA

[View Online »](#)



Anti-SNCA rabbit monoclonal antibody at 1:500 dilution  
 Lane A: 293T Whole Cell Lysate: Lysates/proteins at 30  $\mu$ g per lane. Secondary Goat Anti-Rabbit IgG H&L (Dylight800) at 1/10000 dilution. Developed using the Odyssey technique. Performed under reducing conditions. Observed band size:18 kDa.



Anti-alpha-Synuclein rabbit monoclonal antibody at 1:500 dilution. Lane A: Hela Whole Cell Lysate. Lane B: alpha-Synuclein konckout Hela Whole Cell Lysate. Lysates/proteins at 30  $\mu$ g per lane. Secondary Goat Anti-Rabbit IgG (H+L)/HRP at 1/10000 dilution. Developed using the ECL technique. Performed under reducing conditions. Observed band size:16 kDa

## Handling & Storage

**Handling** Protected from light. Avoid freeze-thaw cycles.

**Long Term Storage** -20°C

**Shipping** Blue Ice

**Regulatory Status** RUO - Research Use Only

## Product Details

**Alternative Name** SNCA

**Application** ELISA, WB

**Clone** 017

**Formulation** Liquid. 0.2 µm filtered solution in PBS. Preservative free.

**Host** Rabbit

**Immunogen** Recombinant human SNCA / alpha-Synuclein protein.

**Isotype** IgG

**Purity Detail** Protein A purified.

**Recommendation** Western Blot (1:500-1:1'000)ELISA (1:5'000-1:10'000)Suggested dilutions/conditions  
**Dilutions/Conditions** may not be available for all applications.Optimal conditions must be determined individually for each application.

**Species Reactivity** Human

**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-  
fr@enzolifesciences.com](mailto:info-fr@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)