

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

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

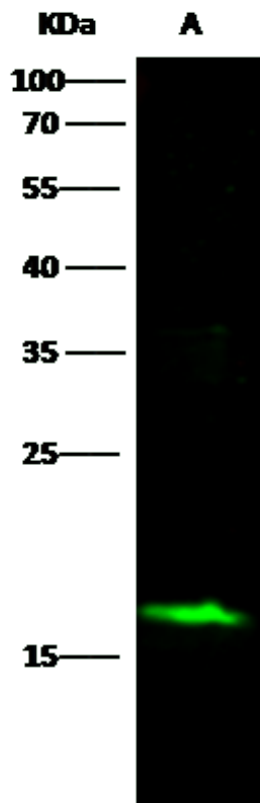
## Ordering Information

[Order Online »](#)

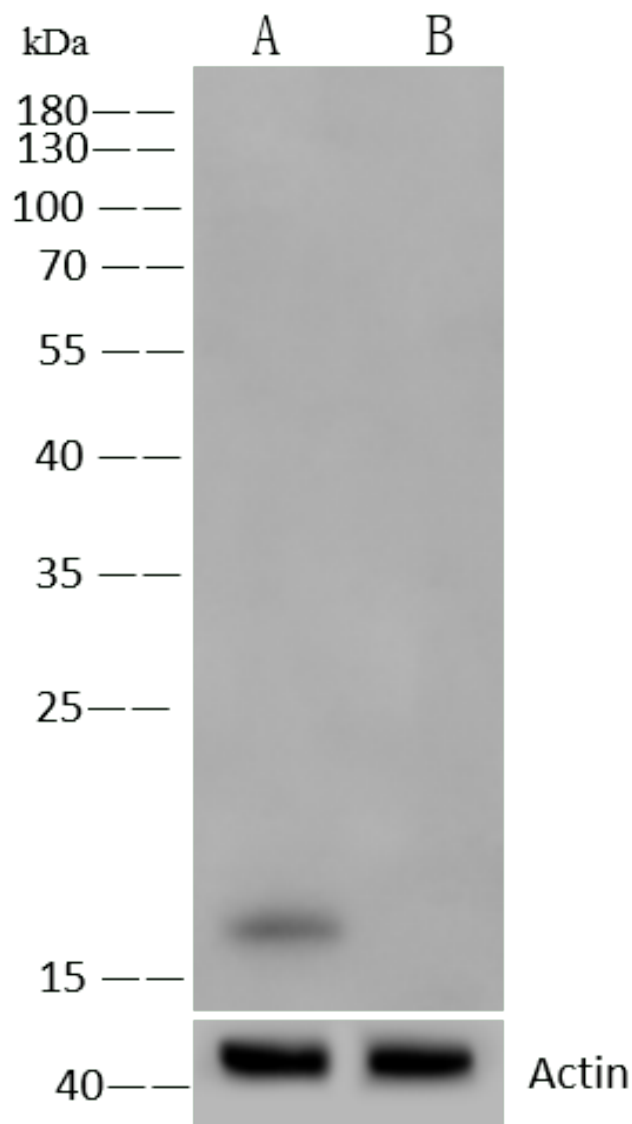
|                 |             |
|-----------------|-------------|
| ENZ-ABS989-0100 | 100 $\mu$ l |
|-----------------|-------------|

## 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 Dilutions/Conditions | Western Blot (1:500-1:1'000)ELISA (1:5'000-1:10'000)Suggested 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 <a href="#">Worry-Free Guarantee</a>  |

Last modified: September 17, 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)