

# TNF- $\alpha$ polyclonal antibody

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

## Ordering Information

[Order Online »](#)

ALX-210-335-C100	100 $\mu$ g
------------------	-------------

## Manuals, SDS & CofA

[View Online »](#)

## Handling & Storage

Handling	Avoid freeze/thaw cycles. After opening, prepare aliquots and store at -20°C.
Long Term Storage	+4°C
Shipping	Blue Ice

## Regulatory Status

RUO - Research Use Only

## Product Details

Alternative Name	Tumor necrosis factor- $\alpha$ , TNFSF 2
Application	ELISA, FUNC, WB
Application Notes	Functional Application: Neutralizes cytotoxic activity of recombinant rat TNF- $\alpha$ . Neutralizing activity: $>2 \times 10^5$ U/mg. One unit is defined as the amount of antibody sufficient for neutralizing the cytotoxic activity of 1 unit recombinant rat TNF- $\alpha$ .
Formulation	Liquid. In PBS.
Host	Rabbit
Immunogen	Recombinant rat TNF- $\alpha$ (tumor necrosis factor- $\alpha$ ).
Recommendation Dilutions/Conditions	Western Blot (10ng recombinant rat TNF- $\alpha$ by 40-60 $\mu$ g/ml PAb to TNF- $\alpha$ , Prod. No. ALX-210-335)Suggested dilutions/conditions may not be available for all applications.Optimal conditions must be determined individually for each application.
Species Reactivity	Human, Mouse, Rat
UniProt ID	P16599
Worry-free Guarantee	This antibody is covered by our <a href="#">Worry-Free Guarantee</a>



Last modified: May 29, 2024

ENZO LIFE SCIENCES,  
INC.  
Phone: 800.942.0430  
[info@enzolifesciences.com](mailto:info@enzolifesciences.com)

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

Belgium, The Netherlands  
& Luxembourg  
Phone: +32 3 466 0420  
[info@enzolifesciences.com](mailto:info@enzolifesciences.com)  
[be@enzolifesciences.com](mailto:be@enzolifesciences.com)

France  
Phone: +33 472 440 655  
[info@enzolifesciences.com](mailto:info@enzolifesciences.com)  
[fr@enzolifesciences.com](mailto:fr@enzolifesciences.com)

Germany  
Phone: +49 7621 5500 526  
[info@enzolifesciences.com](mailto:info@enzolifesciences.com)  
[de@enzolifesciences.com](mailto:de@enzolifesciences.com)

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