

TNF- α polyclonal antibody

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

Citations: 1

[View Online »](#)

Ordering Information

[Order Online »](#)

ALX-210-3350	100 μ 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- α , TNFSF 2
Application	ELISA, FUNC, WB
Application Notes	Functional Application: Neutralizes cytotoxic activity of recombinant rat TNF- α . 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- α .
Formulation	Liquid. In PBS.
Host	Rabbit
Immunogen	Recombinant rat TNF- α (tumor necrosis factor- α).
Recommendation Dilutions/Conditions	Western Blot (10ng recombinant rat TNF- α by 40-60 μ g/ml PAb to TNF- α , 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 Worry-Free Guarantee



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430

info@enzolifesciences.com
www.enzolifesciences.com

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989

info@enzolifesciences.com
eu@enzolifesciences.com

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

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

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

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

Last modified: May 29, 2024