

# TNF-R2 monoclonal antibody (80M2)

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

Citations: 1

[View Online »](#)

## Ordering Information

[Order Online »](#)

ALX-804-450-C100	100µg
------------------	-------

Manuals, SDS & CofA

[View Online »](#)

# Handling & Storage

Long Term Storage +4°C

Shipping Blue Ice

# Regulatory Status

RUO - Research Use Only

# Product Details

Alternative Name CD120b, p75, TNF-R 80kDa, TNFRSF 1B

Application ELISA, Flow Cytometry, FUNC, IHC (FS), IP, WB

Application Notes Not suitable for IHC (PS).  
Functional Application: non-agonistic receptor modulation – enhances in vitro TNF responses by increasing the affinity of soluble TNF-α to TNF-R2.

Clone 80M2

Formulation Liquid. 0.2µm-filtered solution in PBS containing 0.1% BSA.

Host Mouse

Immunogen Human TNF-R2.

Isotype IgG1

Positive Control Human lymph nodes can be used as positive control for immunohistochemistry, PHA activated T cells for flow cytometry.

Recommendation Dilutions/Conditions Typical starting dilution is 1:10Suggested dilutions/conditions may not be available for all applications.Optimal conditions must be determined individually for each application.

Species Reactivity Human, Rat

Specificity Recognizes the extracellular part of human and rat TNF-R2 and soluble human TNF-R2.

## Technical Info / Product Notes

The reactivity of the antibody with cell-bound TNF receptor is minimally inhibited by high concentrations of TNF- $\alpha$ .

## UniProt ID

P20333

## Worry-free Guarantee

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

Last modified: May 29, 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)