

TNF-R2 monoclonal antibody (80M2)

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

Citations: 1

[View Online »](#)

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. 0.2 μ m-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 1:10. Suggested 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- α .

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

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

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

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

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

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