

TNF- α (human) monoclonal antibody (TNF-D)

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

Citations: 1

[View Online »](#)

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Handling	After opening aliquot contents and freeze at -20°C. Avoid freeze/thaw cycles.
Long Term Storage	+4°C
Shipping	Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name	TNFSF 2, Tumor necrosis factor- α
Application	ELISA, Flow Cytometry, FUNC
Application Notes	Functional Application: selectively neutralizes human TNF- α bioactivity. TNF-D requires an incubation of antigen and antibody for 1 hour at 37 °C or 16 hours at 4 °C to completely neutralize TNF- α cytolytic activity. Not recommended for Western blot.
Clone	TNF-D
Crossreactivity	Does not cross-react with mouse TNF- α or human or mouse TNF- β .
Formulation	Liquid. In PBS.
Host	Mouse
Immunogen	Recombinant human TNF- α (tumor necrosis factor- α).
Isotype	IgG1
Recommendation Dilutions/Conditions	ELISA (use as capture or detection antibody in sandwich ELISA)Flow Cytometry (suitable as primary antibody for staining in FACS analysis; recommended concentration is 10 μ g/ml as determined on PBMC. Method: PBMC were incubated with PMA, ionomycin, and with monensin to inhibit secretion. Incubation for 20 minutes with TNF-D at 10 μ g/ml; stained with goat-anti-mouse IgG-FITC second step reagent for 15 minutes)Suggested dilutions/conditions may not be available for all applications.Optimal conditions must be determined individually for each application.

Species Reactivity

Human

UniProt ID

P01375

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)