

# TNF- $\alpha$ (human) monoclonal antibody (TNF-D)

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

Citations: 1

[View Online »](#)

## Ordering Information

[Order Online »](#)

|                  |             |
|------------------|-------------|
| ALX-804-199-C100 | 100 $\mu$ g |
|------------------|-------------|

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- $\alpha$   |
| Application                            | ELISA, Flow Cytometry, FUNC  |
| Application Notes                      | Functional Application: selectively neutralizes human TNF- $\alpha$ 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- $\alpha$ cytolytic activity.<br>Not recommended for Western blot.  |
| Clone                                  | TNF-D  |
| Crossreactivity                        | Does not cross-react with mouse TNF- $\alpha$ or human or mouse TNF- $\beta$ .   |
| Formulation                            | Liquid. In PBS.  |
| Host                                   | Mouse  |
| Immunogen                              | Recombinant human TNF- $\alpha$ (tumor necrosis factor- $\alpha$ ).  |
| Isotype                                | IgG1   |
| Recommendation<br>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 $\mu$ 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 $\mu$ 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](#).



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)