TNF-α (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

100µg

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-

 $\boldsymbol{\alpha}$ 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-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).

lsotype lgG1

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 10μg/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 10μg/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



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

ENZO LIFE SCIENCES,

Phone: 800.942.0430

info-

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

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

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

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