

# TNF-R1 (human) monoclonal antibody (H398)

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

Citations: 16

[View Online »](#)

Ordering Information

[Order Online »](#)

ALX-804-2000	100µg
--------------	-------

Manuals, SDS & CofA

[View Online »](#)

## Handling & Storage

Handling	Avoid freeze/thaw cycles. After opening, prepare aliquots and store at -20°C.
Long Term Storage	+4°C
Shipping	Blue Ice

**Regulatory Status** RUO - Research Use Only

## Product Details

Alternative Name	TNF-R 60kDa, CD120a, p55, TNFRSF 1A
Application	Flow Cytometry, FUNC, IHC (FS)
Application Notes	<p>Not recommended for IHC (PS).</p> <p>Functional Application: blocks TNF-R signalling. 10µg/ml H398 neutralizes TNF induced cytotoxicity on U937 (100%). Furthermore, 10µg/ml H398 neutralizes TNF induced MHC expression on K562, Colo 205 and IL2R on YT cells (always 100%).</p> <p>TNF induced cytotoxicity can be inhibited/neutralized by H398 on MCF-7, HeLa and L929. A working concentration could not be determined, since ascites antibody preparations have been used which exhibit an unspecific cytotoxic effect on the cells. Functional activity via the TNF-R1 is completely inhibited by H398 although on some cells the ligand still binds. Thus, the antibody H398 inhibits signalling via the receptor but not always the binding of TNF.</p>
Clone	H398
Crossreactivity	Does not cross-react with mouse TNF-R1.
Formulation	Liquid. In PBS.
Host	Mouse
Immunogen	Human TNF-R1.
Isotype	IgG2a
Purity Detail	Protein G affinity purified.

## Recommendation Dilutions/Conditions

Immunohistochemistry (frozen sections; stains acetone-fixed cryostat sections or cell smears. Recommended for H398 is the alkaline phosphatase-anti-alkaline-phosphatase (APAAP) or peroxidase anti-peroxidase (PAP) procedures or the three stage immunoperoxidase technique on acetone-fixed cryostat sections. H398 may be used at a concentration of 1-0.5mg/ml for cryostat sections) Suggested dilutions/conditions may not be available for all applications. Optimal conditions must be determined individually for each application.

## Species Reactivity

Human

## Specificity

Recognizes membrane-bound and soluble TNF-R1.

## UniProt ID

P19438

## 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)