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

This antibody is covered by our Worry-Free Guarantee.

Citations: 16

View Online »

**Ordering Information** 

Order Online »

ALX-804-200-C100

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**Not recommended for IHC (PS).

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%).

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.

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

