Fas (human):Fc (human), (recombinant)

Citations: 14

View Online »

Ordering Information

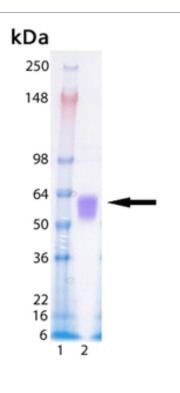
Order Online »

ALX-522-002-C050

50µg

Manuals, SDS & CofA

View Online »



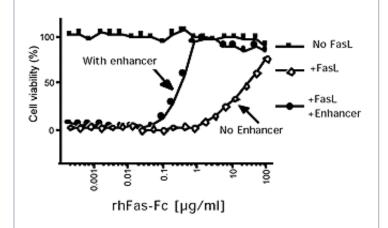


Figure: Inhibition of FasL, Soluble (human) (rec.) (Prod. No. <u>ALX-522-001</u>)-mediated lysis. Fas (human):Fc (human) (rec.) (Prod. No. ALX-522-002) exerts its inhibitory activity in a concentration range of 0.5-5μg/ml in the presence of the enhancer (1μg/ml).**Method:** Mouse A20 B lymphoma cells (50'000 cells in 100μl DMEM medium containing 5% fetal calf serum) were incubated with 0.2μg/ml FasL

Handling & Storage

Use/Stability Stable for at least 6 months after receipt when stored at -20°C.

Handling Avoid freeze/thaw cycles. After reconstitution, prepare aliquots and store at -20°C.

Long Term Storage -20°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name CD95, APO-1, TNFRSF 6

Application Notes ELISA: binds to FasL at 10-100 ng/ml.

Concentration 1mg/ml after reconstitution.

Endotoxin Content <0.1EU/µg purified protein (LAL test; Associates of Cape

Cod).

Formulation Lyophilized. Contains PBS.

MW ∼56 kDa

Purity ≥95% (SDS-PAGE)

Purity Detail Purified by multi-step chromatography.

Reconstitution Reconstitute with 50µl sterile water. Further dilutions can

be made with medium containing 5% fetal calf serum or a

carrier protein.

Source Produced in HEK 293 cells. The extracellular domain of

human Fas (CD95; APO-1) (aa 7-154) is fused to the Fc

portion of human IgG1.

Specificity Binds human, mouse and rat FasL (APO-1L; CD95L;

CD178).

Technical Info / Product Notes

Historical data has shown that Fas inhibits the activity of human and mouse soluble Fas (sFasL) and thereby sFasL-mediated lysis of Fas sensitive cells in a concentration range of 20-100µg/ml. When used with cross-linking enhancer (see Prod. No. ALX-203-001), the inhibitory activity of rhFas:Fc is increased approximately 20- to 50-fold. Concentrations of rhFas:Fc required to inhibit may vary depending on the cell viability and on the concentrations of sFasL used to kill the cells.

UniProt ID

P25445

Last modified: May 29, 2024

