FasL (soluble) (human), (recombinant)

Fas ligand

Citations: 44

View Online »

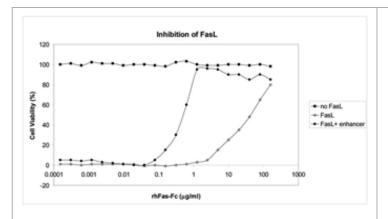
Ordering Information

Order Online »

ALX-522-001-C010	10µg
ALX-522-001-3010	SuperPack - 3x10µg

Manuals, SDS & CofA

View Online »



Activity Assay analysis: Inhibition of FasL, Soluble (human) (rec.) (Prod. No. ALX-522-001)-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).

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 APO-1L, CD95L, CD178, TNFSF 6

Application Notes ELISA: binds to Fas receptor at 1-100 ng/ml.

Biological Activity

Kills Jurkat cells in the absence of a cross-linker. Cross-

linking enhancer (see Set Prod. No. ALX-850-014) increases the activity of rhsFasL approx. 50-fold. **Attention:** Results using sFasL may differ from those

obtained with agonistic antibodies!

Concentration 0.1mg/ml after reconstitution.

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

Cod).

Formulation Lyophilized. Contains PBS.

MW ~32kDa (non-glycosylated), ~35kDa (glycosylated) (SDS-

PAGE).

Purity ≥95% (SDS-PAGE)

Purity Detail Purified by multi-step chromatography.

Reconstitution Reconstitute with 100µl sterile water. Further dilutions

should be made with medium containing 5% fetal calf

serum.

Source

Produced in HEK 293 cells. The extracellular domain of human FasL (APO-1L; CD95L; CD178) (aa 103-281) is fused at the N-terminus to a linker peptide (26 aa) and a FLAG[®]-tag. Glycosylation of rhsFasL is similar to that of natural human FasL.

Specificity

Binds to human, mouse and rat Fas (CD95; APO-1).

Technical Info / Product Notes

Historical lots have shown that FasL kills Jurkat cells at concentrations of >10ng/ml in the absence of a cross-linker. Historical lots have also shown an ED₅₀ of 50ng/ml (A20 cells).

UniProt ID

P48023

Last modified: May 29, 2024



Phone: 800.942.0430

usa@enzolifesciences.com

info-