

FasL (soluble) (human), (recombinant)

Fas ligand

Citations: 44

[View Online »](#)

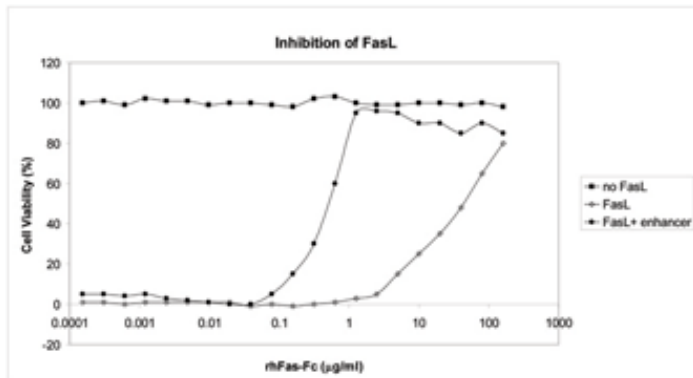
Ordering Information

[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	<p>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.</p> <p>Attention: Results using sFasL may differ from those obtained with agonistic antibodies!</p>
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	<p>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.</p>
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



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)