

FasL, soluble (human) (recombinant) set

Citations: 25

[View Online »](#)

Ordering Information

[Order Online »](#)

ALX-850-014-KI02	1Set
------------------	------

Manuals, SDS & CofA

[View Online »](#)

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	FUNC
Application Notes	Kills Fas-sensitive cells at concentrations of >1ng/ml in the presence of cross-linking enhancer.
Concentration	FasL: 0.1mg/ml after reconstitutionEnhancer: 1mg/ml
Contents	10µg of FasL, Soluble (human) (recombinant) (Prod. No. ALX-522-001) and 4×50µg of Enhancer for Ligands (Prod. No. ALX-804-034) which increases the biological activity (induction of apoptosis) by ~50-fold to a concentration range of 1-5ng/ml.
Formulation	FasL: Lyophilized. Contains PBS. Enhancer: Liquid. In PBS, pH7.4, containing 50% glycerol. Contains no preservatives.
Purity	≥95% (SDS-PAGE)
Reconstitution	Reconstitute FasL with 100µl sterile water. Further dilutions should be made with medium containing 5% fetal calf serum or carrier protein.
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 FasL, Soluble (human) (recombinant) is similar to natural human FasL.

Specific ActivityED₅₀: 1ng/ml (A20 cells)**Specificity**

FasL binds to human, mouse and rat Fas (CD95).

Technical Info / Product Notes

FLAG is a registered trademark of Sigma-Aldrich Co.

UniProt ID

P48023

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)