

SUPERFASLIGAND® Protein

(soluble) (human), (recombinant)

Fas ligand with improved stability providing significantly enhanced immune activation.

SUPERFASLIGAND® (soluble) (human), (recombinant) features an N-terminus linker shown to improve stability and enhance immune activation compared to recombinant ligands alone.

Citations: 72

[View Online »](#)

Ordering Information

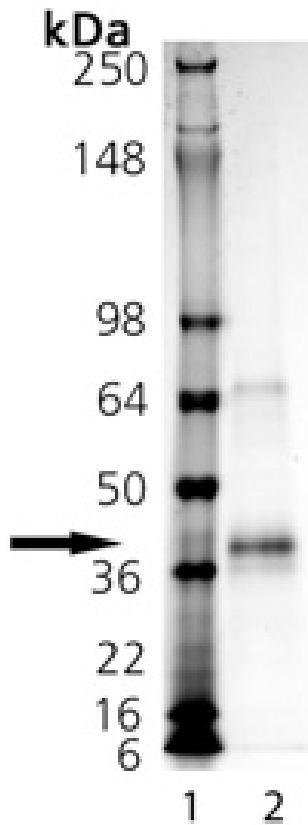
[Order Online »](#)

ALX-522-020-C005	5µg
ALX-522-020-3005	SuperPack - 3x5µg

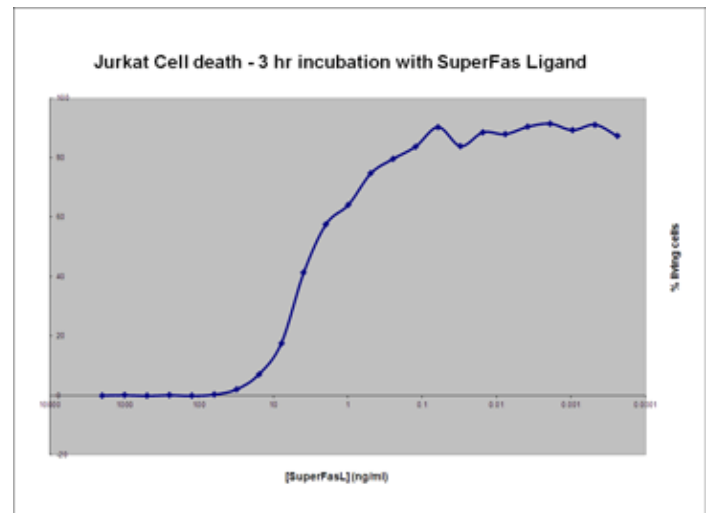
Manuals, SDS & CofA

[View Online »](#)

- Increased stability
- Enhanced immune activation compared to other recombinant ligands
- Mimics glycosylation of native human FasL



SDS-PAGE analysis: Lane 1: MW Marker, Lane 2: 1 μ g SuperFasL, stained with Imperial stain.



Jurkat Cell Death: 50,000 Jurkat cells per well are incubated with the indicated concentration of SuperFasL for 3 hours. Cell death is determined by Cell-titer AQueous one-solution cell proliferation reagent. Formazan product is allowed to develop for 6 hours before the plate is read at 490nm. % Viability is determined in comparison to control well with no SuperFasL.

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	FasL (oligomer), APO-1L (oligomer), CD95L (oligomer), CD178 (oligomer), TNFSF 6 (oligomer)
Application Notes	ELISA: binds to Fas receptor at 1-100 ng/ml.
Biological Activity	Kills Fas-sensitive cells. Note: Does not require enhancer.
Concentration	0.1mg/ml after reconstitution.
Formulation	Lyophilized. Contains PBS.
MW	~32kDa (nonglycosylated), ~35kDa (glycosylated).
Purity	≥95% (SDS-PAGE)
Reconstitution	Reconstitute with 50µl sterile water. Further dilutions should be made with cell culture 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 rhs SUPERFASLIGAND [®] is similar to natural human FasL.
Specificity	Binds to human, mouse and rat Fas (CD95; APO-1).

Technical Info / Product Notes

Historical data has shown that SUPERFASLIGAND[®] kills Fas-sensitive cells at concentrations of >1ng/ml without the use of enhancer. The ED₅₀ has been shown in previous data to be 1ng/ml (A20 cells).

Note: Results using rhsSUPERFASLIGAND[®] may differ from those obtained with agonistic antibodies!

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)