SUPERFASLIGAND® Prof (soluble) (human), (recombinant)

Fas ligand with improved stability providing significantly enhanced immune activation.

SUPERFASLIGAND $^{\circledR}$ (soluble) (human), (recombinant) features an N-terminus linker shown to improve stability and enhance immune activation compared to recombinant ligands alone.

Citations: 71

View Online »

Ordering Information

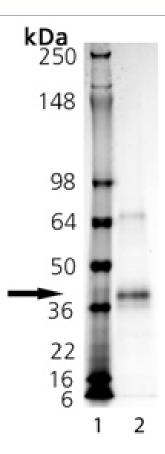
Order Online »

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

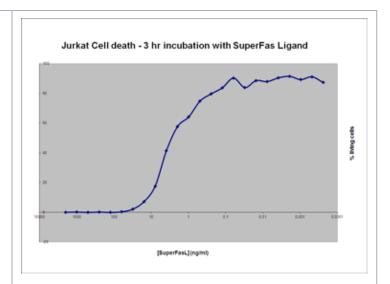
Manuals, SDS & CofA

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: 1µg 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 ActivityKills 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 ED $_{50}$ 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

