

Fas (human):Fc (human), (recombinant)

Citations: 14

[View Online »](#)

Ordering Information

[Order Online »](#)

ALX-522-002-C050	50µg
------------------	------

Manuals, SDS & CofA

[View Online »](#)

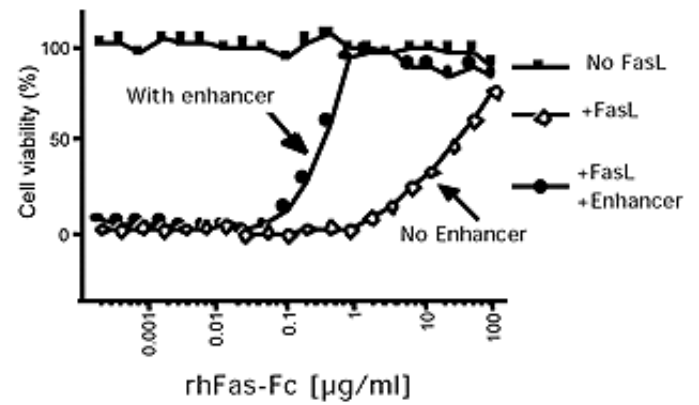
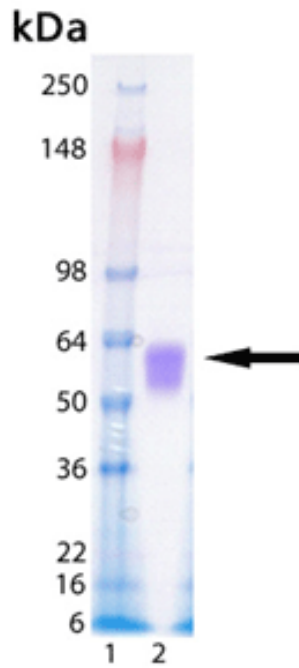


Figure: Inhibition of FasL, Soluble (human) (rec.) (Prod. No. [ALX-522-001](#))-mediated lysis. Fas (human):Fc (human) (rec.) (Prod. No. ALX-522-002) exerts its inhibitory activity in a concentration range of 0.5-5μg/ml in the presence of the enhancer (1μg/ml). **Method:** Mouse A20 B lymphoma cells (50'000 cells in 100μl DMEM medium containing 5% fetal calf serum) were incubated with 0.2μg/ml FasL

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	CD95, APO-1, TNFRSF 6
Application Notes	ELISA: binds to FasL at 10-100 ng/ml.
Concentration	1mg/ml after reconstitution.
Endotoxin Content	<0.1EU/μg purified protein (LAL test; Associates of Cape Cod).
Formulation	Lyophilized. Contains PBS.
MW	~56 kDa
Purity	≥95% (SDS-PAGE)
Purity Detail	Purified by multi-step chromatography.
Reconstitution	Reconstitute with 50μl sterile water. Further dilutions can be made with medium containing 5% fetal calf serum or a carrier protein.
Source	Produced in HEK 293 cells. The extracellular domain of human Fas (CD95; APO-1) (aa 7-154) is fused to the Fc portion of human IgG1.
Specificity	Binds human, mouse and rat FasL (APO-1L; CD95L; CD178).

Technical Info / Product Notes

Historical data has shown that Fas inhibits the activity of human and mouse soluble Fas (sFasL) and thereby sFasL-mediated lysis of Fas sensitive cells in a concentration range of 20-100µg/ml. When used with cross-linking enhancer (see Prod. No. ALX-203-001), the inhibitory activity of rhFas:Fc is increased approximately 20- to 50-fold. Concentrations of rhFas:Fc required to inhibit may vary depending on the cell viability and on the concentrations of sFasL used to kill the cells.

UniProt ID

P25445

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)