

DcR3 (human):Fc (human), (recombinant)

Citations: 2

[View Online »](#)

Ordering Information

[Order Online »](#)

ALX-522-055-C050	50µg
------------------	------

Manuals, SDS & CofA

[View Online »](#)

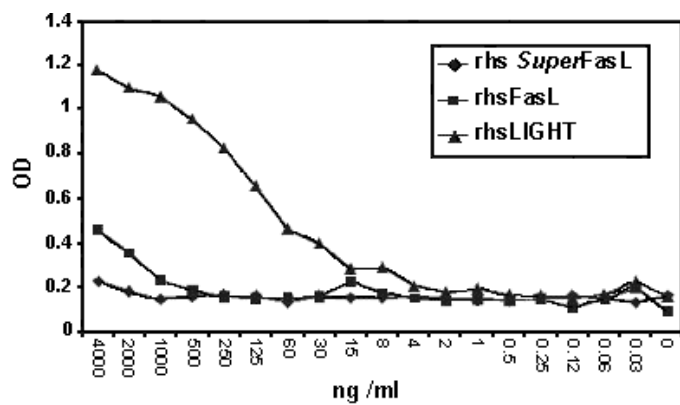


Figure:

Method: A 96-well ELISA plate was coated O/N with 50 ng rh DcR3:Fc (Prod. No. [ALX-522-055](#)) per well. After blocking, the indicated concentrations of rhsFasL (Prod. No. **ALX-522-001**)

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	TNFRSF 6B, Decoy receptor 3
Application Notes	ELISA: binds to LIGHT. Detection of human LIGHT and TL1A (VEGI) in combination with PAb to human IgG1.
Concentration	1mg/ml after reconstitution.
Endotoxin Content	<0.1EU/µg purified protein (LAL test; Bio Whittaker).
Formulation	Lyophilized. Contains PBS.
MW	~58kDa (SDS-PAGE).
Purity	≥90% (SDS-PAGE)
Reconstitution	Reconstitute with 50µl sterile water.
Source	Produced in HEK 293 cells. The extracellular domain of human DcR3 (decoy receptor 3) (aa 1-246) is fused to the Fc portion of human IgG1.
Specificity	Binds human LIGHT. Does not bind human FasL.
UniProt ID	O95407



Last modified: May 29, 2024

ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
usa@enzolifesciences.com

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
info-be@enzolifesciences.com

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
info-fr@enzolifesciences.com

France
Phone: +33 472 440 655
info-fr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info-de@enzolifesciences.com

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
info-uk@enzolifesciences.com