

# LIGHT (soluble) (human), (recombinant)

Citations: 6

[View Online »](#)

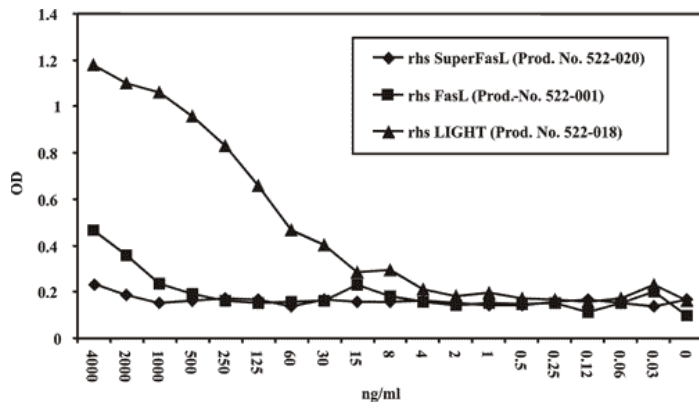
Ordering Information

[Order Online »](#)

ALX-522-0180	10µg
--------------	------

Manuals, SDS & CofA

[View Online »](#)



**Figure:** Binding of recombinant human DcR3:Fc to soluble human recombinant LIGHT.

## Handling & Storage

<b>Use/Stability</b>	Stable for at least 6 months after receipt when stored at -20°C.
<b>Handling</b>	Avoid freeze/thaw cycles. After reconstitution, prepare aliquots and store at -20°C.
<b>Long Term Storage</b>	-20°C
<b>Shipping</b>	Blue Ice

**Regulatory Status** RUO - Research Use Only

## Product Details

<b>Alternative Name</b>	HVEML, TNFSF 14, CD258
<b>Application Notes</b>	ELISA: binds to HVEM receptor at 1-10 ng/ml.
<b>Concentration</b>	0.1mg/ml after reconstitution.
<b>Endotoxin Content</b>	<0.1EU/μg purified protein (LAL test; Associates of Cape Cod).
<b>Formulation</b>	Lyophilized. Contains PBS.
<b>MW</b>	~25kDa (SDS-PAGE)
<b>Purity</b>	≥95% (SDS-PAGE)
<b>Reconstitution</b>	Reconstitute with 100μl sterile water. Further dilutions should be made with medium containing 5% fetal calf serum.
<b>Source</b>	Produced in HEK 293 cells. The extracellular domain of human LIGHT (aa 89-240) is fused at the N-terminus to a linker peptide (10 aa) and a FLAG <sup>®</sup> -tag.
<b>Specificity</b>	Binds to human and mouse LTβR and human HVEM and DcR3.
<b>Technical Info / Product Notes</b>	Historical data shows that the biological activity of LIGHT increases LNCaP (prostate cancer cell) growth.  FLAG is a registered trademark of Sigma-Aldrich Co.

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)