

# CD40L (soluble) (human), (recombinant)

Increases the stimulation of B cells

Citations: 37

[View Online »](#)

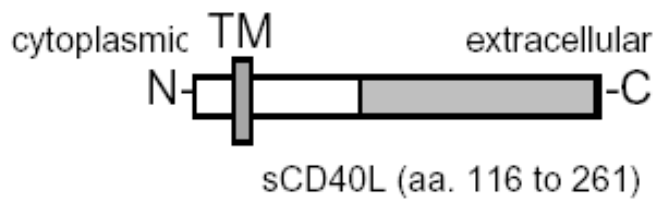
## Ordering Information

[Order Online »](#)

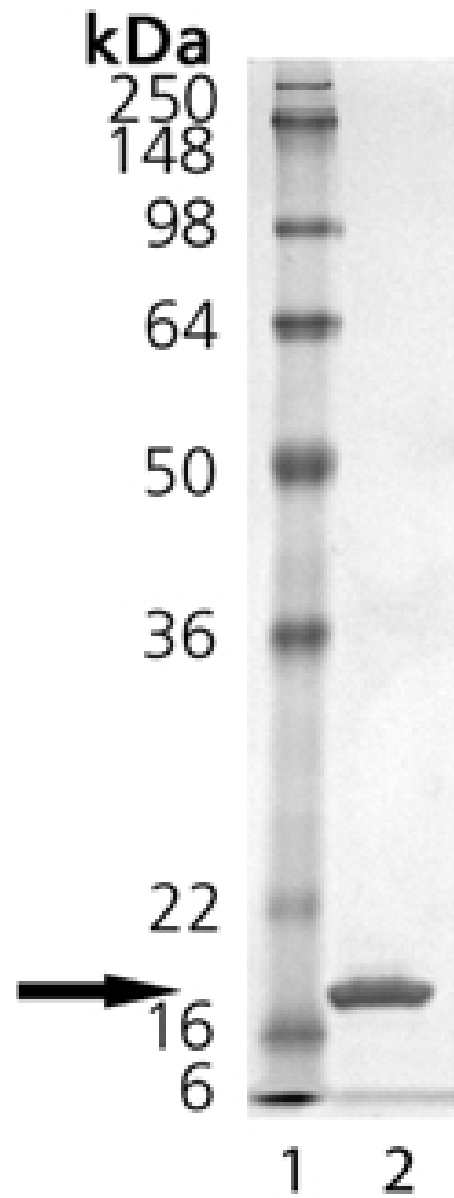
ALX-522-015-C010	10µg
ALX-522-015-2010	2x10µg
ALX-522-015-6010	SuperPack - 6x10µg

Manuals, SDS & CofA

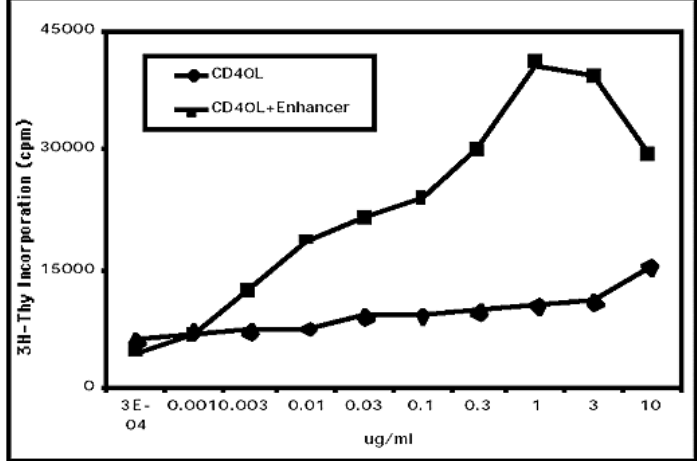
[View Online »](#)



*Schematic structure of human CD40L (peptide, aa. 116 to 261).*



SDS-PAGE Analysis of CD40L (soluble) (human), (recombinant). Lane 1: MW Marker, Lane 2: 1  $\mu$ g.



**Figure 2:** Stimulation of B cells by CD40L, Soluble (human) (rec.). Dose dependent costimulation of CD19+ cells by CD40L in the presence of 10ng/ml of IL-4 and of 2µg/ml anti-µ antibodies. Note that in the presence of enhancer (Prod. No. ALX-804-034)

## 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	CD154, TNFSF 5, gp39
Application Notes	ELISA: binds to CD40 receptor at 1-10 ng/ml.
Concentration	0.1mg/ml after reconstitution.
Endotoxin Content	<0.1EU/μg purified protein (LAL test; Associates of Cape Cod).
Formulation	Lyophilized. Contains PBS.
MW	~18kDa (SDS-PAGE).
Purity	≥95% (SDS-PAGE)
Purity Detail	Purified by Multi-Step Chromatography.
Reconstitution	Reconstitute each tube with 100μl sterile water. Further dilutions should be made with medium containing 5% fetal calf serum or a carrier protein.
Source	Produced in <i>E. coli</i> . The extracellular domain of human CD40L (CD154) (aa 116-261) is fused at the N-terminus to a linker peptide (6 aa) and a FLAG <sup>®</sup> -tag.
Specificity	Binds to human CD40.

## Technical Info / Product Notes

Historical data has shown that CD40L stimulates growth of B cells. The activity of rhsCD40L increases 1'000-fold (stimulation in the ng/ml range) in the presence of cross-linking enhancer (see Set Prod. No. ALX-850-064). For stimulation of mouse cells via CD40 use CD40L, Soluble (mouse) (recombinant) (Prod. No. ALX-522-070 or ALX-850-075).

FLAG is a registered trademark of Sigma-Aldrich Co.

UniProt ID

P29965

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)