

# MEGACD40L<sup>®</sup> Protein (soluble) (human), (recombinant)

## High activity, high purity CD40L protein for co-stimulatory activation of an immune response

MEGACD40L<sup>®</sup> is a high activity protein in which two trimeric CD40 ligand molecules are artificially linked via the collagen domain of Adiponectin/ACRP30/AdipoQ. This protein very effectively simulates the natural membrane-assisted aggregation of CD40L *in vivo*. It provides a simple and equally potent alternative to [CD40L+enhancer] combinations (Prod. No. ALX-850-064).

Citations: 84

[View Online »](#)

## Ordering Information

[Order Online »](#)

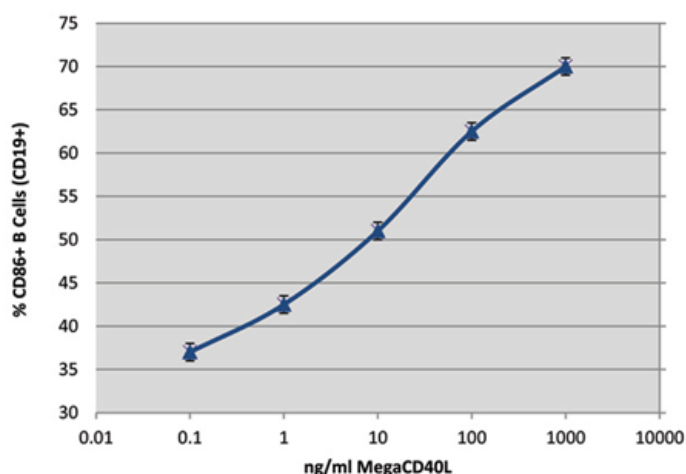
ALX-522-110-C010	10µg
------------------	------

## Manuals, SDS & CofA

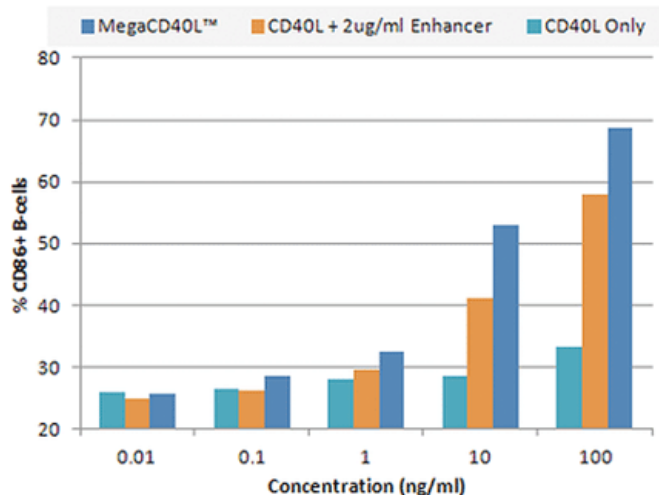
[View Online »](#)

- Highly active MEGACD40L<sup>®</sup> oligomer mimics *in vivo* membrane-assisted CD40L aggregation and stimulation without the need for enhancers
- High purity with low endotoxin levels
- Experienced manufacturing results in consistent and dependable lot-to-lot performance
- Optimized formulation for improved stability and enhanced activity

### Mimic Natural CD40L-mediated Immune Response

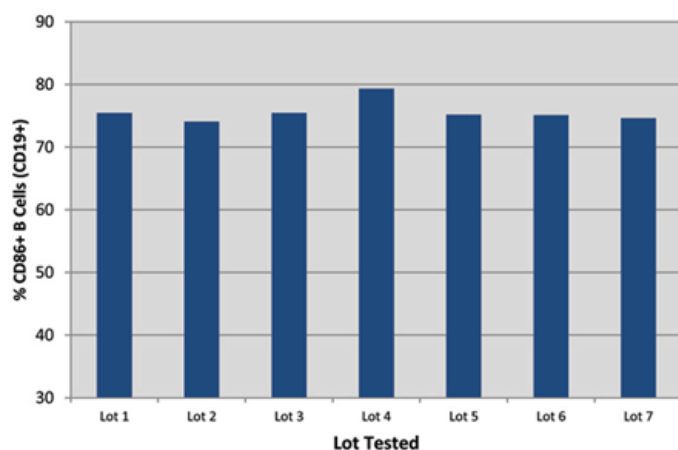


B cell lymphocyte activation by MEGACD40L<sup>®</sup> (Prod. No. ALX-522-110). Method: Peripheral Blood Mononuclear Cells (PBMCs) were incubated at 37°C, 5% CO<sub>2</sub> for 48 hours in 48 well plates (1 x 10<sup>6</sup> cells/well). Each well contained 200µl serum free test media with serially diluted MEGACD40L<sup>®</sup>. After treatment, cells were washed and dual-stained with mouse anti-human CD19–PE and mouse anti-human CD86–APC and analyzed by flow cytometry. The data are presented as the percent of CD19 positive B cells that co-stain as CD86 positive, at each concentration of MEGACD40L<sup>®</sup>.



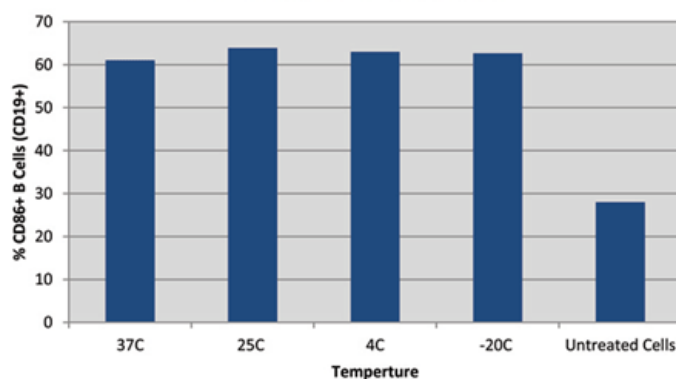
PBMCs (peripheral blood mononuclear cells) were incubated at 37°C, 5% CO<sub>2</sub> for 48 hours in 48 well plates (1 x 10<sup>6</sup> cells per well) containing 200µl serum free test media with serially diluted MEGACD40L<sup>®</sup> (Prod. No. ALX-522-110), CD40L (Prod. No. ALX-522-015), or CD40L + 2µg/ml Enhancer (Prod. No. ALX-804-034). After treatment the cells were washed 3X with PBS + 1% BSA. The cells were dual stained for 2 hours at 4°C with mouse anti-human CD19 (PE conjugate) and mouse anti-human CD86 (APC conjugate). The cells were washed 2X with PBS + 1% BSA, then re-suspended in PBS. Samples were analyzed on a BD FACS Calibur flow cytometer. The data is presented as the percent of CD86 positive cells per CD19 positive B cells at each concentration.

**Reliable Manufacturing Yields Consistent Results**

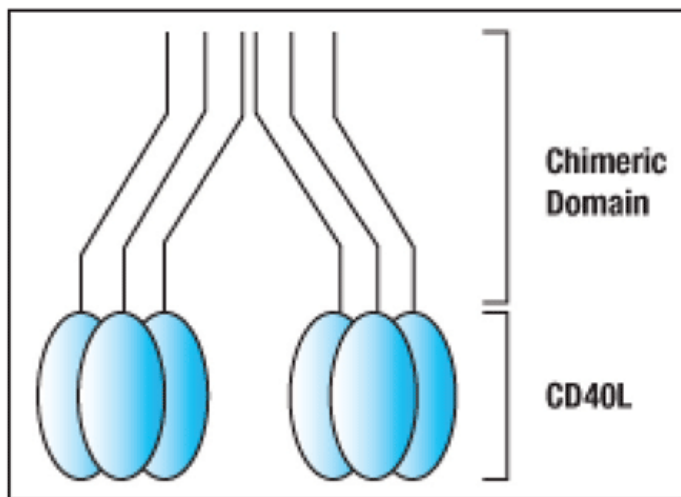


MEGACD40L<sup>®</sup> from 7 consecutive manufacturing lots was tested using a B-cell activation assay at 1000ng/mL.

**High-Stability Protein Construct**



MEGACD40L<sup>®</sup> was subjected to stability testing by storing the protein at the indicated temperature for 2 months. The protein was then tested for activity using a B cell activation assay. Results indicate MEGACD40L<sup>®</sup> retains activity under even harsh storage conditions.



Graphical representation of MEGACD40L<sup>®</sup>, Soluble (human) (recombinant) (Prod. No. ALX-522-110).

# 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	CD40L:ACRP30headless, CD154:ACRP30headless, TNFSF 5:ACRP30headless, gp39:ACRP30headless
Application Notes	Can be used to mediate immune and inflammatory responses, as a protein tool for pre-clinical studies in immunotherapy, to monitor patients undergoing chemotherapy and to mimic autoimmune phenotypes <i>in vivo</i> .
Biological Activity	Binds to human CD40. Induces B cells activation (as demonstrated by dose-dependent upregulation of CD86).
Concentration	0.1mg/ml after reconstitution.
Endotoxin Content	≤0.01EU/μg purified protein (LAL test; Cape Cod Associates).
Formulation	Lyophilized. Contains PBS.
MW	~38kDa
Purity	≥90% (SDS-PAGE)
Purity Detail	Purified by multi-step chromatography.
Reconstitution	Reconstitute with 100μl sterile water. Further dilutions should be made with medium containing 5% fetal calf serum.

<b>Source</b>	Produced in CHO cells. The extracellular domain of human CD40L (CD154) (aa 116-261) is fused at the N-terminus to mouse ACRP30 <i>headless</i> (aa 18-111) and a FLAG <sup>®</sup> -tag.
<b>Technical Info / Product Notes</b>	FLAG is a registered trademark of Sigma-Aldrich Co.
<b>UniProt ID</b>	P29965 (human CD40L), Q60994 (mouse Adiponectin)

**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)