

MEGACD40L[®] Protein (soluble) (mouse), (recombinant)

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

MEGACD40L[®] is a high activity protein in which two trimeric CD40 ligands are artificially linked via the collagen domain of ACRP30/adiponectin. This protein very effectively simulates the natural membrane-assisted aggregation of CD40L. It provides a simple and equally potent alternative to [CD40L+enhancer] combinations (Prod. No. ALX-850-075).

Citations: 14

[View Online »](#)

Ordering Information

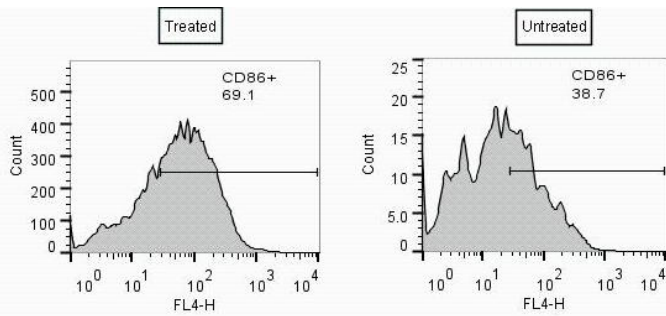
[Order Online »](#)

ALX-522-120-C010	10µg
------------------	------

Manuals, SDS & CofA

[View Online »](#)

- Highly active MEGACD40L[®] 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



B cell lymphocytes activation after titration with MEGACD40L[®] (soluble) (mouse) (rec.). Method: PBMCs (Peripheral Blood Mononuclear Cells) were incubated at 37°C, 5% CO₂ for 48 hours in 48 well plates (1 x 10⁶ cells per well) containing 200µl serum free test media with or without 1000ng/mL MEGACD40L[®] (soluble) (mouse) (rec.) (Prod. No. ALX-522-120). 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 by on a BD FACS Calibur flow cytometer. The data is presented as the percent of CD86 positive cells per CD19 positive B cells with or without treatment with 1000ng/mL mega mouse CD40L.

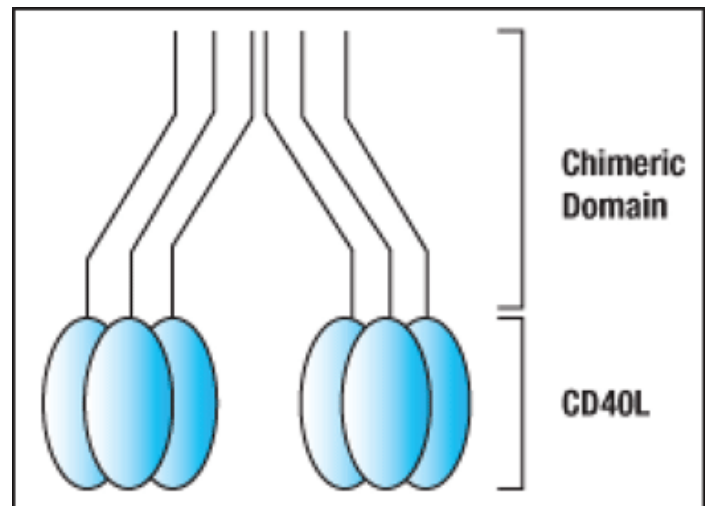
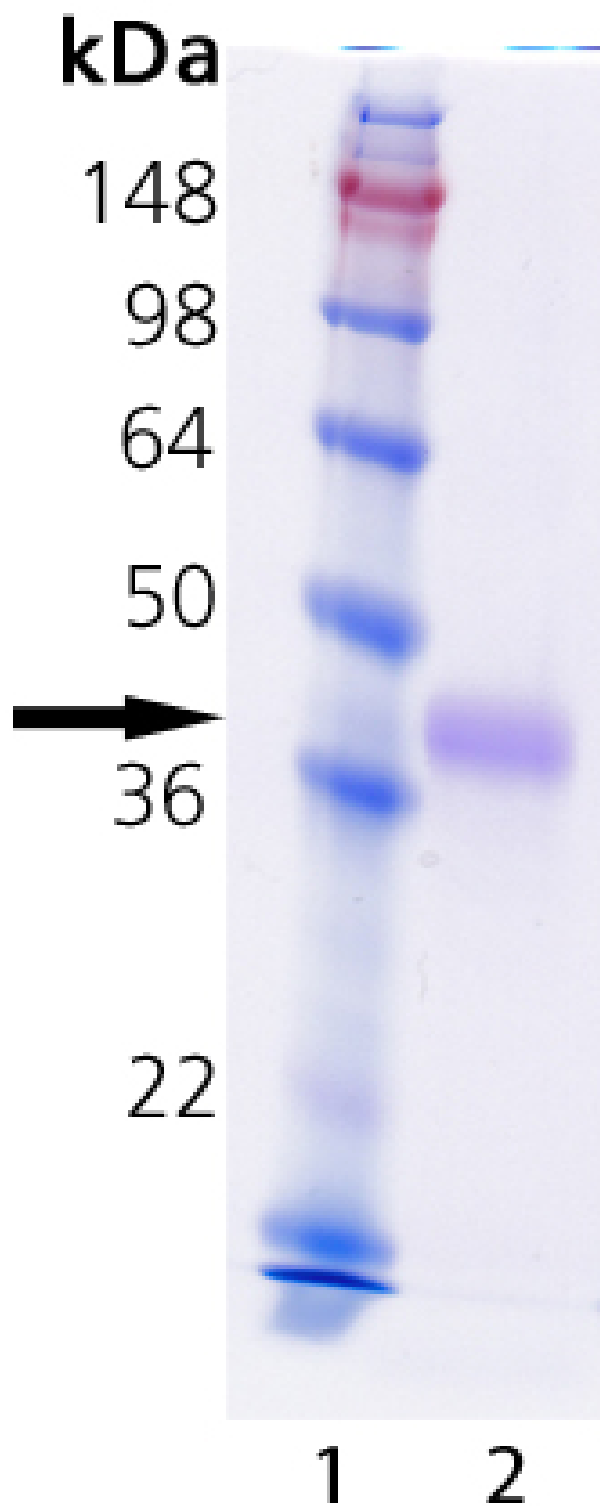


Figure 1: Graphical representation of MEGACD40L[®] (soluble) (mouse) (recombinant) (Prod. No. ALX-522-120).



Coomassie stained SDS-PAGE. Lane 1: Molecular weight marker. Lane 2: 1.0 μg MEGACD40L[®] (soluble) (mouse) (rec.) (Prod. No. ALX-522-120).

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, TNFSF5:ACRP30headless, gp39:ACRP30headless
Biological Activity	Binds to mouse and human CD40. Induces B cells activation (as demonstrated by dose-dependent upregulation of CD86).
Concentration	0.1mg/ml after reconstitution.
Endotoxin Content	<0.1EU/µg purified protein (LAL test; Associates of CAPE COD Inc.).
Formulation	Lyophilized. Contains PBS.
MW	~35-40kDa
Purity	≥95% (SDS-PAGE)
Reconstitution	Reconstitute with 100µl sterile water. Further dilutions should be made with medium containing 5% fetal calf serum.
Source	Produced in CHO cells. The extracellular domain of mouse CD40L (CD154) (aa 115-260) is fused at the N-terminus to mouse ACRP30 <i>headless</i> (aa 18-111) and a FLAG [®] -tag.
Technical Info / Product Notes	FLAG is a registered trademark of Sigma-Aldrich Co.
UniProt ID	P27548 (CD40L), Q60994 (ACRP30)



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)