

Fc (human):CD40L (soluble) (mouse), (recombinant)

Citations: 1

[View Online »](#)

Ordering Information

[Order Online »](#)

ALX-522-072-C010	10µg
------------------	------

Manuals, SDS & CofA

[View Online »](#)

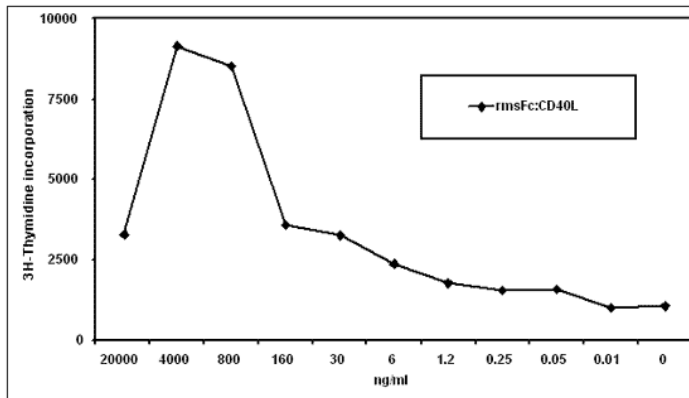
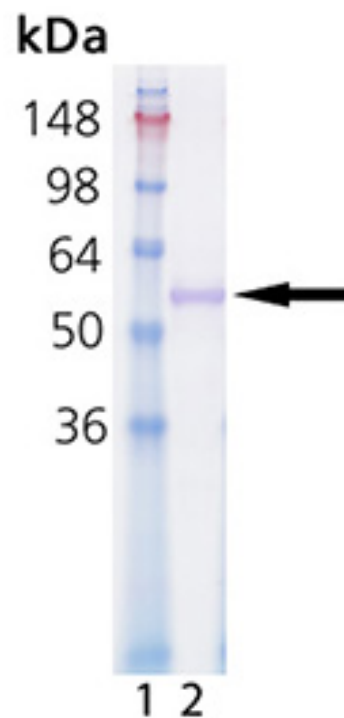


Figure: Proliferation of mouse B cells induced by Fc (human):CD40L, Soluble (mouse) (rec.) (Prod. No. [ALX-522-072](#)). **Method:** Splenocytes from Balb/c mice were purified on B220 magnetic beads using the MACS system. The resulting B cells were put in culture at a cell density of 70'000 cells/well in a 96-well plate. Cells were activated with the indicated concentrations of Fc (human):CD40L, Soluble (mouse) (rec.). After 36 hours, 0.5µCi of 3H-Thymidine/well was added. Cells were pulsed for 10 hours, then freeze-thawed, harvested and counted."",Fc (human):CD40L (soluble) (mouse)



SDS-PAGE analysis of Fc (human):CD40L (soluble) (mouse), (recombinant) (Prod. no. ALX-522-072): Lane 1: MW Marker, Lane 2: 1 µg.

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 human CD40:COMP at 10-100 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	~58kDa (SDS-PAGE).
Purity	≥90% (SDS-PAGE)
Reconstitution	Reconstitute with 100μl sterile water. Further dilutions should be made with medium containing 5% fetal calf serum or a carrier protein.
Source	Produced in HEK 293 cells. The extracellular domain of mouse CD40L (CD154) (aa 115-260) is fused at the N-terminus to the Fc portion of human IgG1 and a linker peptide (20 aa).
Specificity	Binds to human and mouse CD40.
Technical Info / Product Notes	Historical data has shown that Fc (human):CD40L (mouse) stimulates the proliferation of mouse B cells and dendritic cells. Note: Does not require an enhancer.



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

P27548
European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
info-eu@enzolifesciences.com

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
info-be@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