

CD40L (soluble) (mouse), (recombinant)

Citations: 2

[View Online »](#)

Ordering Information

[Order Online »](#)

ALX-522-070-C010	10µg
ALX-522-070-2010	2x10µg

Manuals, SDS & CofA

[View Online »](#)

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 at 10-500 ng/ml.
Concentration	0.1mg/ml after reconstitution.
Endotoxin Content	<50EU/mg purified protein (LAL test; Associates of Cape Cod).
Formulation	Lyophilized. Contains PBS.
MW	~26kDa (SDS-PAGE).
Purity	≥90% (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 HEK 293 cells. The extracellular domain of mouse CD40L (CD154) (aa 115-260) is fused at the N-terminus to a linker peptide (8 aa) and a FLAG [®] -tag.
Specificity	Binds to human and mouse CD40.
Technical Info / Product Notes	<p>Historical data has shown that mouse CD40L stimulates the proliferation of mouse B cells and dendritic cells. The activity of rmsCD40L increases 1'000-fold (stimulation in the ng/ml range) in the presence of a cross-linking enhancer (see Set Prod. No. ALX-850-075).</p> <p>FLAG is a registered trademark of Sigma-Aldrich Co.</p>



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)