

# CD40 (human):Fc (human), (recombinant)

Citations: 6

[View Online »](#)

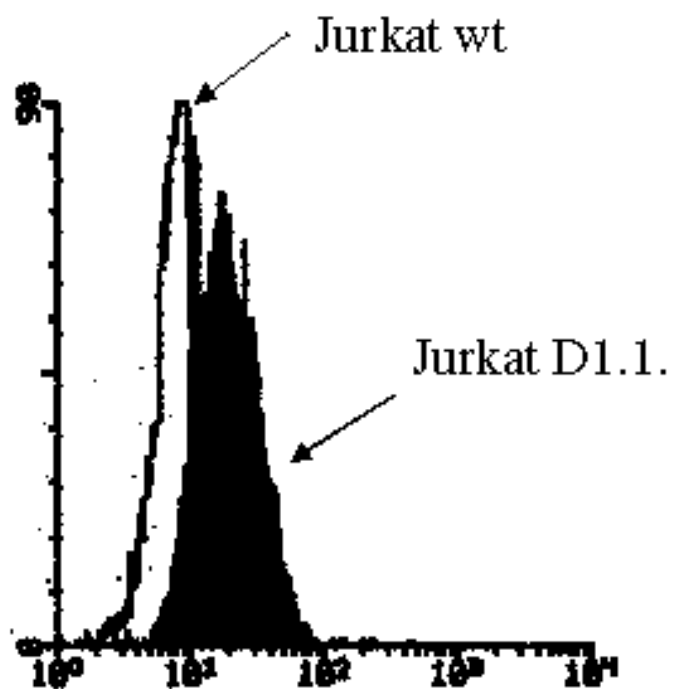
Ordering Information

[Order Online »](#)

ALX-522-016-C050	50µg
------------------	------

Manuals, SDS & CofA

[View Online »](#)



**Figure:** Jurkat T cells clone D1.1 expresses high surface levels of CD40L. Cells were immunostained at 4°C for 30 minutes with 10µg/ml of CD40:Fc (Prod. No [ALX-522-016](#)). After washing, cells were incubated for 30 min. at 4°C with 10µg/ml goat anti-human RPE (Prod. No ALX-211-501). After washing

# 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	TNFRSF 5
Application Notes	ELISA: binds to CD40L at 1-10 ng/ml. Can be used together with anti-human IgG1 PAb to detect membrane-bound human or mouse CD40L. CD40:Fc should be used at concentrations between 1 and 100µg/ml depending on the level of CD40L surface expression.
Concentration	1mg/ml after reconstitution.
Endotoxin Content	<0.1EU/µg purified protein (LAL test; Associates of CAPE COD Inc).
Formulation	Lyophilized. Contains PBS.
MW	~54kDa (SDS-PAGE).
Purity	≥95% (SDS-PAGE)
Reconstitution	Reconstitute with 50µl 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 cysteine-rich region (aa 21-193) of human CD40 is fused to the Fc portion of human IgG1.
Specificity	Binds human and mouse CD40L (CD154).
UniProt ID	P25942



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)