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

Citations: 1

View Online »

**Ordering Information** 

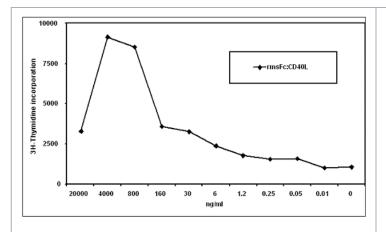
Order Online »

ALX-522-072-C010

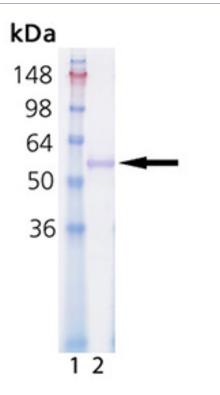
10µg

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.

UniProt ID P27548

Last modified: May 29, 2024

