

Validation data for recombinant human CD40L

<https://www.invivogen.com/human-cd40l>

For research use only

Version 23J16-AK

CD40 ligand, CD40L (also known as CD154, TRAP or gp39), is a type II transmembrane glycoprotein belonging to the TNF family. Its size and purity was assessed using SDS PAGE (**Figure 1**). The biological activity of CD40L has been confirmed using InvivoGen's HEK-Blue™ CD40L cells (**Figure 2**). Binding of CD40L to its receptor CD40 on the surface of HEK-Blue™ CD40L cells triggers a signaling cascade leading to the activation NF- κ B and the subsequent production of SEAP. Levels of SEAP can be easily measured using QUANTI-Blue™, a detection reagent.

Detection by SDS-PAGE

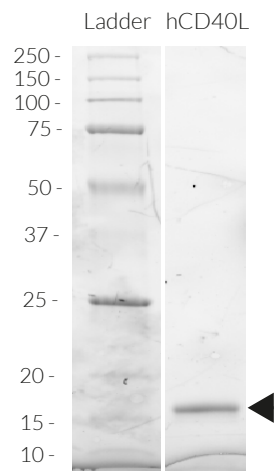


Figure 1. SDS PAGE of the recombinant human (h)CD40L protein. 1.5 μ g of hCD40L was loaded on a 12% Mini-PROTEAN® TGX Stain-Free™ Precast Gel (Bio-Rad). Detection was performed as per the manufacturer's instructions. A band was detected at ~16 kDa.

Dose-response in HEK cells to human CD40L

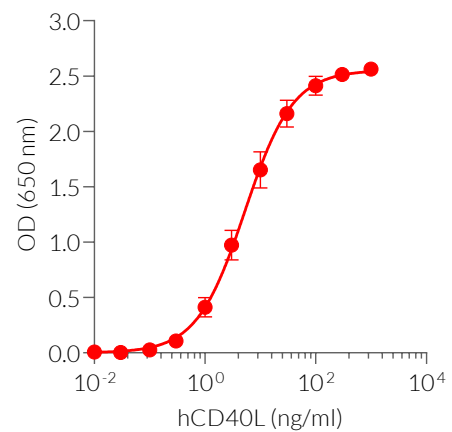


Figure 2. Dose-response in HEK-Blue™ CD40 cells to recombinant hCD40L protein. Cells were stimulated with increasing concentrations of recombinant human (h)CD40L. After overnight incubation, the NF- κ B-induced SEAP activity was determined using QUANTI-Blue™, a SEAP detection reagent. Data are shown as optical density (OD) at 650 nm (mean \pm SEM).

TECHNICAL SUPPORT

InvivoGen USA (Toll-Free): 888-457-5873
InvivoGen USA (International): +1 (858) 457-5873
InvivoGen Europe: +33 (0) 5-62-71-69-39
InvivoGen Asia: +852 3622-3480
E-mail: info@invivogen.com