

# Validation data for Fc-hCD70

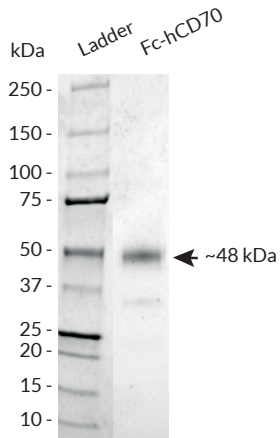
<https://www.invivogen.com/cd70-fc>

For research use only

Version 24D29-NJ

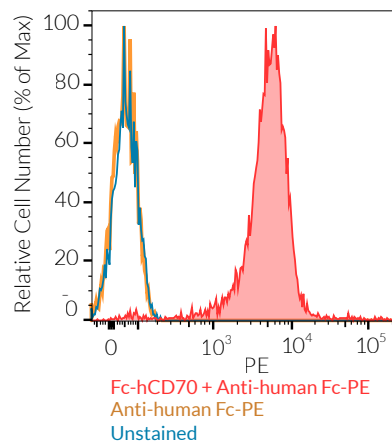
Fc-hCD70 is a soluble human CD70 chimera protein generated by fusing the N-terminal extracellular domain of human CD70 to the C-terminus of a human IgG1 Fc domain with a TEV (Tobacco Etch Virus) sequence linker. Fc-hCD70 has an apparent molecular weight of ~48 kDa on an SDS-PAGE gel (Figure 1). It has been functionally validated by the detection of cell surface hCD27 using flow cytometry (Figure 2), binding of an anti-hCD70 monoclonal antibody (mAb) using ELISA (Figure 3), as well as NF- $\kappa$ B activation in Jurkat-Lucia™ hCD27 reporter cells (Figure 4).

## Fc-hCD70 analysis by SDS-PAGE



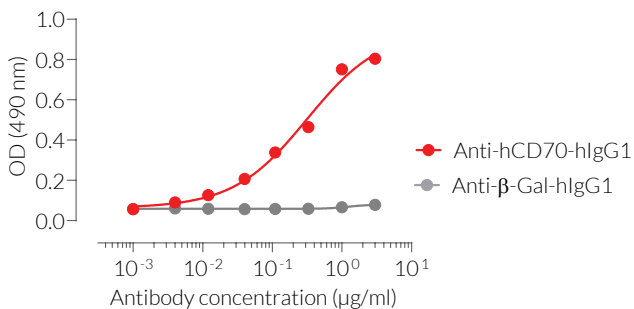
**Figure 1: SDS-PAGE analysis of the Fc-hCD70 protein.** 0.5  $\mu$ g of the fusion protein was loaded on a 12% Mini-PROTEAN® TGX Stain-Free™ Precast Gels (Bio-Rad). Detection was performed as per the manufacturer's instructions.

## Cell surface staining using Fc-hCD70



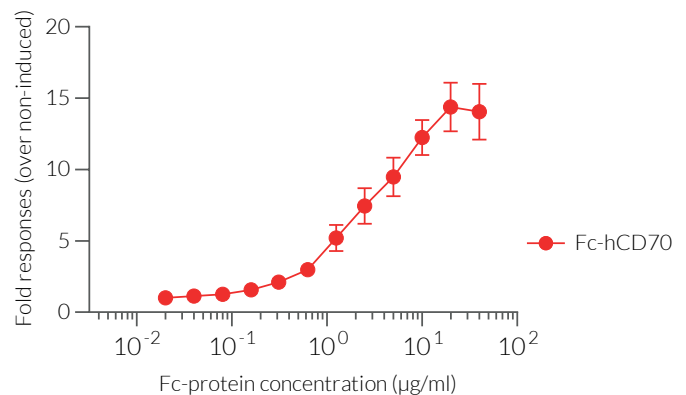
**Figure 2: Human CD27 cell surface detection using Fc-hCD70.**  $\sim 5 \times 10^5$  Jurkat-Lucia™ hCD27 cells were incubated with 2  $\mu$ g of Fc-hCD70 for 30 min at 4°C. Cells were then washed and incubated with 1  $\mu$ l of mouse anti-human IgG Fc antibody coupled to PE for 30 min at 4°C. Cell surface staining was analyzed by flow cytometry.

## ELISA detection of Fc-hCD70



**Figure 3: ELISA detection of Fc-hCD70 with Anti-hCD70-hIgG1 mAb.** The Fc-hCD70 fusion protein was performed and coated at 2  $\mu$ g/ml on ELISA plates overnight. A 3-fold serial dilution of Anti-hCD70-hIgG1 (red curve) or Anti- $\beta$ -Gal-hIgG1 control mAb (grey curve) was added for the capture step. An HRP-labeled anti-human  $\kappa$  light chain antibody (1/1000 dilution) and the HRP substrate OPD (o-phenylenediamine dihydrochloride) were used for the detection step. Absorbance was read at 490 nm.

## Activation of Jurkat-Lucia™ hCD27 cells



**Figure 4: Activation of Jurkat-Lucia™ hCD27 cells.** Jurkat-Lucia™ hCD27 cells were incubated with increasing concentrations of recombinant human Fc-CD70 fusion protein for 24 hours. The NF- $\kappa$ B activation in Jurkat-Lucia™ hCD27 cells was assessed by determining Lucia luciferase activity in the supernatant using QUANTI-Luc™ 4. Fold responses are shown as 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 3-622-34-80  
E-mail: [info@invivogen.com](mailto:info@invivogen.com)

 **InvivoGen**  
[www.invivogen.com](http://www.invivogen.com)