

Validation data for hCD27-Fc

<https://www.invivogen.com/cd27-fc>

For research use only

Version 24D23-NJ

hCD27-Fc is a soluble human CD27 chimera protein generated by fusing the N-terminal extracellular domain of human CD27 to the N-terminus of a human IgG1 Fc domain with a TEV (Tobacco Etch Virus) sequence linker. hCD27-Fc has an apparent molecular weight of ~55 kDa on an SDS-PAGE gel (Figure 1). It has been functionally validated by the detection of cell surface hCD70 using flow cytometry (Figure 2) and binding of an anti-hCD27 monoclonal antibody (mAb) using ELISA (Figure 3).

hCD27-Fc analysis by SDS-PAGE

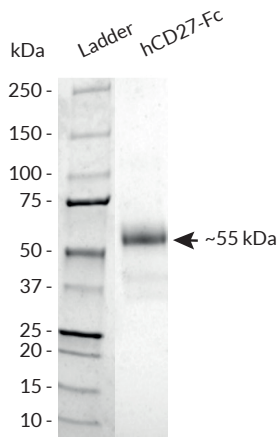


Figure 1: SDS-PAGE analysis of the hCD27-Fc protein. 0.5 μ g of the fusion protein was loaded on a 12% Mini-PROTEAN® TGX Stain-Free™ Precast Gels (Bio-Rad). Detection was performed as per the manufacturer's instructions.

Cell surface staining using hCD27-Fc

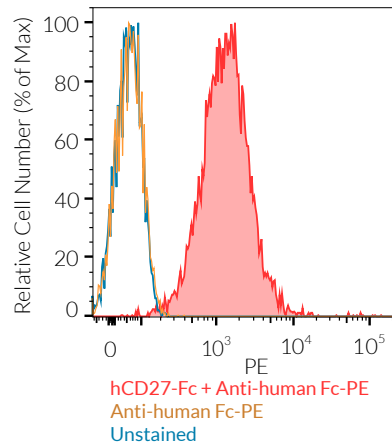


Figure 2: Human CD70 cell surface detection using hCD27-Fc. $\sim 5 \times 10^5$ Raji-APC-Null cells were incubated with 2 μ g of hCD27-Fc for 30 min at 4°C. Cells were then washed and incubated with 1 μ l of mouse anti-human IgG Fc antibody coupled to PE for 30 min at 4°C. Cell surface staining was analyzed by flow cytometry.

ELISA detection of hCD27-Fc

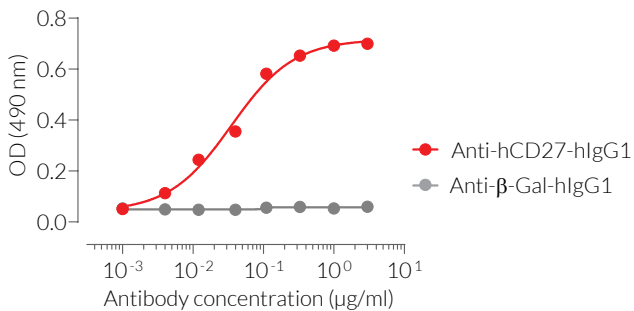


Figure 3: ELISA detection of hCD27-Fc with Anti-hCD27-hIgG1 mAb. The hCD27-Fc fusion protein was performed and coated at 2 μ g/ml on ELISA plates overnight. A 3-fold serial dilution of Anti-hCD27-hIgG1 (red curve) or Anti- β -Gal-hIgG1 control mAb (grey curve) was added for the capture step. An HRP-labeled anti-human κ 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.

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