

Validation data for hICOS-L-Fc

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

For research use only

Version 24D04-NJ

hICOS-L-Fc is a soluble human ICOS-L chimera protein generated by fusing the N-terminal extracellular domain of human ICOS-L to the N-terminus of a human IgG1 Fc domain with a TEV (Tobacco Etch Virus) sequence linker. hICOS-L-Fc has an apparent molecular weight of ~70 kDa on an SDS-PAGE gel (Figure 1). It has been functionally validated by the detection of cell surface hICOS using flow cytometry (Figure 2), binding of an Anti-hICOS-L monoclonal antibody (mAb) using ELISA (Figure 3), as well as NFAT activation in Jurkat-Lucia™-hICOS reporter cells (Figure 4).

hICOS-L-Fc analysis by SDS-PAGE

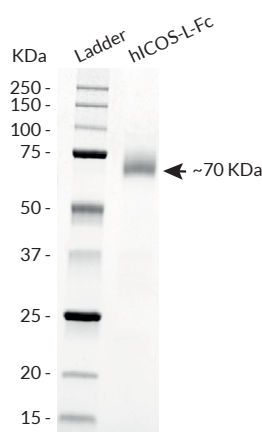


Figure 1: SDS-PAGE analysis of the hICOS-L-Fc protein. 0.5 µg of the fusion protein was loaded on a 12% Mini-PROTEAN® TGX Stain-Free™ Precast Gels (Bio-Rad). Detection was performed as per the manufacturer's instructions.

Cell surface staining using hICOS-L-Fc

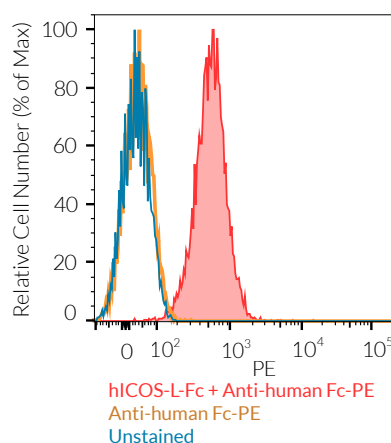


Figure 2: Human ICOS cell surface detection using hICOS-L-Fc. ~5 × 10⁵ Jurkat-Lucia™ hICOS cells were incubated with 500 ng of hICOS-L-Fc for 30 min at 4°C. Cells were then washed and incubated with 1 µl of mouse anti-human IgG Fc antibody coupled to PE for 30 min at 4°C. Cell surface staining was analyzed by flow cytometry.

ELISA detection of hICOS-L-Fc

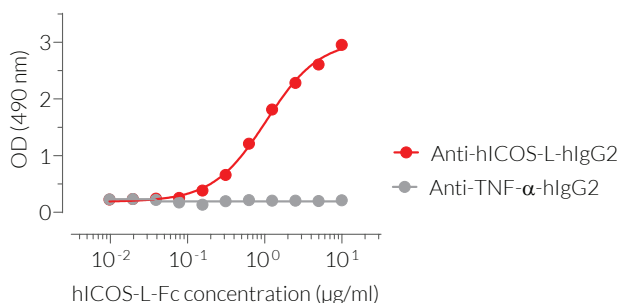


Figure 3: ELISA detection of hICOS-L-Fc with Anti-hICOS-L mAb. A 2-fold serial dilution of the hICOS-L-Fc fusion protein was performed and coated on ELISA plates overnight. Anti-hICOS-L-hlgG2 (red curve) or Anti-TNFα-hlgG2 control mAb (grey curve) at 5 µg/ml was added for the capture step. An HRP-labeled 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.

Activation of Jurkat-Lucia™-hICOS cells

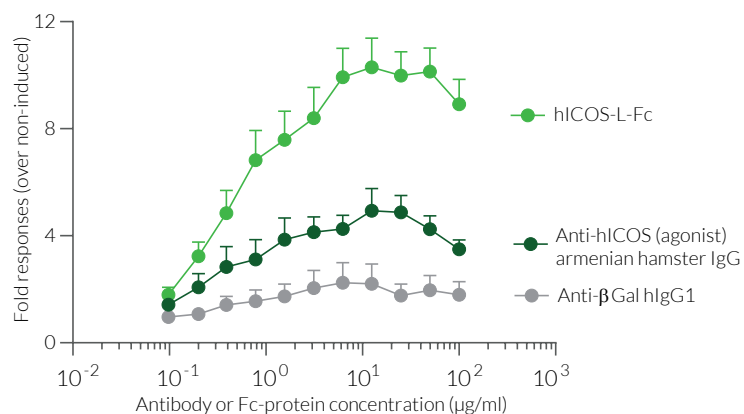


Figure 4: Activation of Jurkat-Lucia™-hICOS cells. Jurkat-Lucia™-hICOS cells were incubated with increasing concentrations of recombinant hICOS-L-Fc fusion protein, agonist Anti-hICOS IgG, or control Anti-βGal hlgG1 mAbs for 6 hours. NFAT activation in the Jurkat-Lucia™-hICOS cells was assessed by determining Lucia luciferase activity in the supernatant using QUANTI-Luc™. Fold responses are shown as mean + 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