

Validation data for hPD1-Fc

<https://www.invivogen.com/pd1-fc>

For research use only

Version 23L08-NJ

hPD1-Fc is a soluble human PD-1 chimera protein generated by fusing the N-terminal extracellular domain of human PD-1 to the N-terminus of a human IgG1 Fc domain with a TEV (Tobacco Etch Virus) sequence linker. hPD1-Fc has an apparent molecular weight of ~62 kDa on an SDS-PAGE gel (Figure 1). It has been functionally validated by the detection of cell surface hPD-L1 using flow cytometry (Figure 2) and by binding of Anti-hPD-1 monoclonal antibodies (mAbs) using ELISA (Figure 3).

hPD1-Fc analysis by SDS-PAGE

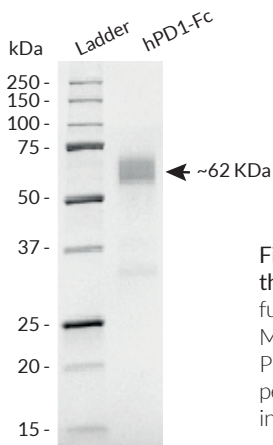


Figure 1: SDS-PAGE analysis of the hPD1-Fc protein. 0.5 μ g of the fusion protein was loaded on a 12% Mini-PROTEAN® TGX Stain-Free™ Precast Gels (Bio-Rad). Detection was performed as per the manufacturer's instructions.

Cell surface staining using hPD1-Fc

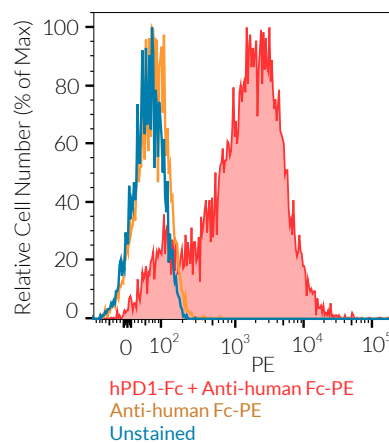


Figure 2: Human PD-L1 cell surface detection using hPD1-Fc. $\sim 5 \times 10^5$ Raji-APC-hPD-L1 cells were incubated with 2 μ g of hPD1-Fc for 45 min at 4°C. Cells were then washed and incubated with 1 μ g of mouse anti-human IgG Fc antibody coupled to PE for 1h at 4°C. Cell surface staining was analyzed by flow cytometry.

ELISA detection of hPD1-Fc

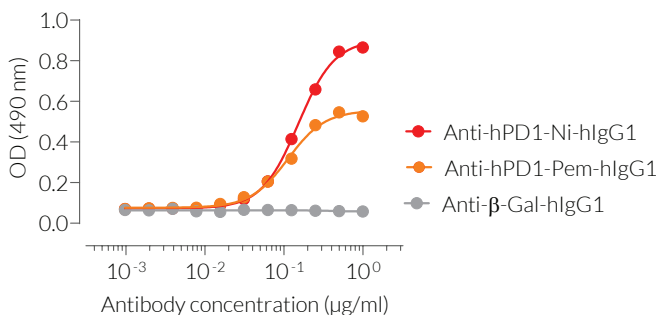


Figure 3: ELISA detection of hPD1-Fc with Anti-hPD1 mAbs. hPD1-Fc fusion protein (1 μ g/ml) was coated on ELISA plates overnight. A 2-fold serial dilution of Anti-hPD1-hlgG1 Nivolumab biosimilar (red curve) or Pembrolizumab biosimilar (orange curve), or Anti- β -Gal-hlgG1 control mAb (grey curve) was performed 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