

Validation data for Anti-PD-L1-mIgG1e3 (Atezo) InvivoFit™

www.invivogen.com/recombinant-anti-mouse-pdl1-atezolizumab-10F.9G2-d265a

For research use only

Version 24C21-AK

Anti-PD-L1-mIgG1e3 (Atezo) InvivoFit™ is a recombinant monoclonal antibody (mAb) designed for *in vivo* studies in mice. This mAb features the variable region of the previously described anti-PD-L1 atezolizumab and the engineered murine IgG1e3 effectorless constant region. Atezolizumab (formerly known as MPDL3280A) is a therapeutic mAb that targets programmed cell death ligand 1 (PD-L1), blocking its interaction with the receptor, PD-1. This mAb binds both mouse and human PD-L1. The binding capacity of Anti-PD-L1-mIgG1e3 (Atezo) InvivoFit™ was compared to the 10F.9G2-derived Anti-mPD-L1 mAb using ELISA and flow cytometry (Figures 1 & 2). Anti-PD-L1-mIgG1e3 (Atezo) InvivoFit™ recognizes both mouse and human PD-L1, unlike Anti-mPD-L1-mIgG1e3 (10F.9G2) InvivoFit™, which is mouse PD-L1-specific (Figure 1).

Evaluation of binding specificity by ELISA

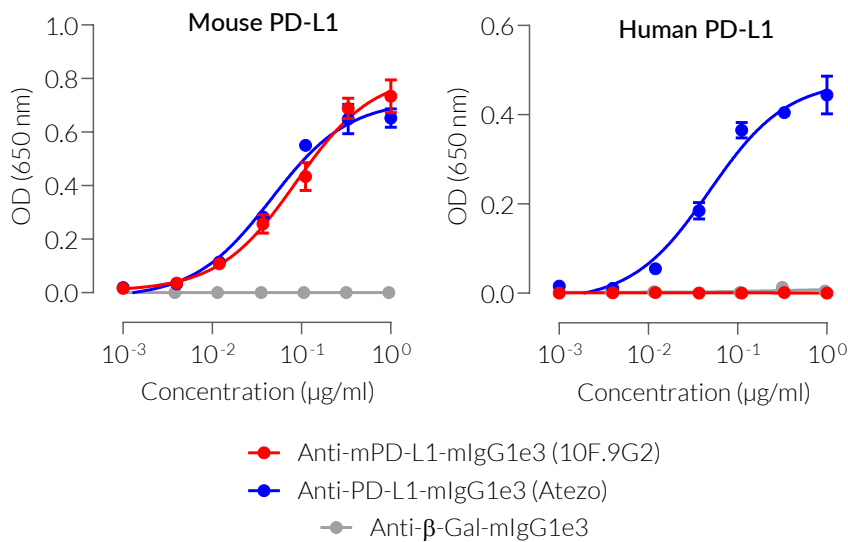


Figure 1. Binding specificity to recombinant mouse or human PD-L1.

Mouse or human PD-L1 (2 µg/ml) was coated on ELISA plates overnight. A 3-fold serial dilution of the Anti-PD-L1-mIgG1e3 (Atezo) InvivoFit™ (blue curve), Anti-mPD-L1-mIgG1e3 (F10.9G2) InvivoFit™ (red curve), or of the Anti-β-Gal-mIgG1e3 control antibody (grey curve) was performed for the capture step. An HRP-labelled anti-mIgG antibody (1/1000 dilution) and the HRP substrate OPD (o-phenylenediamine dihydrochloride) were used for the detection step. Absorbance was read at 650 nm.

Evaluation of binding specificity by flow cytometry

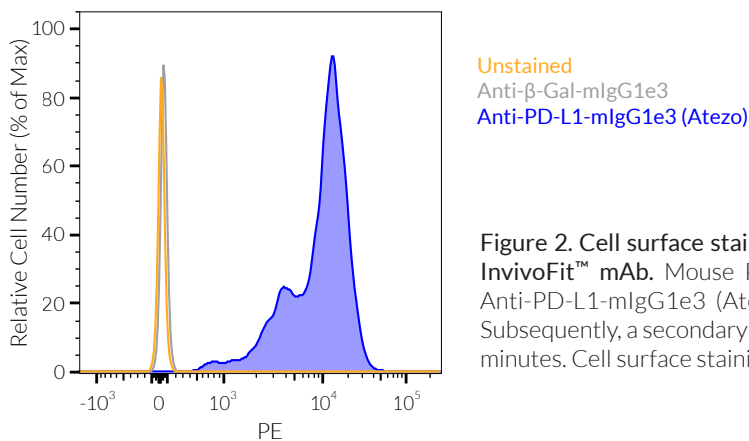


Figure 2. Cell surface staining of mouse PD-L1 using Anti-PD-L1-mIgG1e3 (Atezo) InvivoFit™ mAb. Mouse PD-L1-expressing EL4 cells were incubated with 2 µg of Anti-PD-L1-mIgG1e3 (Atezo) InvivoFit™ mAb or an isotype control for 1h at 4°C. Subsequently, a secondary PE-labeled antibody was added and incubated at 4°C for 30 minutes. Cell surface staining was analyzed by flow cytometry.

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