

# Validation data for Raji-hPD-L1 Cells

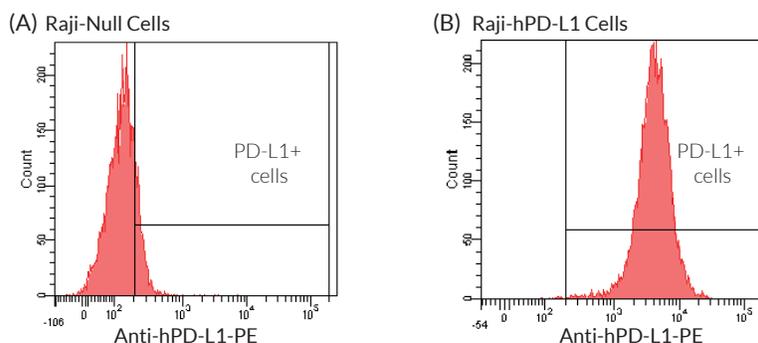
<https://www.invivogen.com/raji-hpd1>

For research use only

Version 20C16-ED

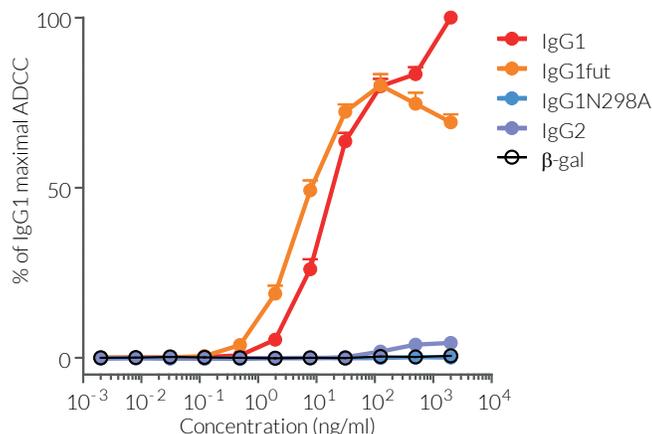
Raji-hPD-L1 cells were developed from the Raji cell line to overexpress the human programmed cell death ligand 1 (PD-L1, also known as CD274 or B7-H1) gene. Raji-hPD-L1 cells were designed as target cells in InvivoGen's antibody-dependent cellular cytotoxicity (ADCC) assay using clinically-relevant anti-human PD-L1 monoclonal antibodies (mAbs). Human PD-L1 expression by Raji-hPD-L1 cells has been verified by flow-cytometry (**Figure 1**), and induction of ADCC has been validated using InvivoGen's combinations of anti-human PD-L1 antibody isotypes and Jurkat-NFAT Lucia™ CD16 reporter cell line (**Figure 2**). The level of ADCC induction is measured as a bioluminescent signal produced by an NFAT-dependent Lucia luciferase reporter protein. Antibodies displaying lower EC<sub>50</sub> have higher ADCC potency.

## Validation of PD-L1 expression



**Figure 1:** Validation of the expression of human PD-L1 by Raji-hPD-L1 cells. Raji-Null (A) and Raji-hPD-L1 (B) cells were incubated with a PE-conjugated Anti-hPD-L1 mAb for 30 minutes. The binding affinity was then measured using flow cytometry.

## ADCC assay using various anti-human PD-L1 (Atezolizumab) antibody isotypes and Raji-hPD-L1 target cells



**Figure 2:** Comparison of ADCC potency for native and engineered anti-human PD-L1 antibody isotypes. Raji-hPD-L1 cells were incubated with gradient concentrations of Anti-hPD-L1 or Anti- $\beta$ -galactosidase ( $\beta$ -gal) mAbs for 1 hour. Jurkat-Lucia™ NFAT-CD16 effector cells were then co-incubated with targets cells for 6 hours. NFAT activation, reflecting the induced ADCC response, was assessed by determining Lucia luciferase activity in the supernatant using QUANTI-Luc™. Percentages of the maximal response normalized to the IgG1 isotype are shown.

### 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 Hong Kong: +852 3622-3480

E-mail: [info@invivogen.com](mailto:info@invivogen.com)

 **InvivoGen**  
www.invivogen.com