

# Anti-hPD-L1-hIgG1

Recombinant monoclonal human IgG1 antibody against human PD-L1

Catalog code: hpd11-mab1, hpd11-mab1-03

<https://www.invivogen.com/anti-human-pdl1-igg1-atezolizumab>

For research use only, not for diagnostic or therapeutic use

Version 23L18-MM

## PRODUCT INFORMATION

**Contents:** Anti-hPD-L1-hIgG1 purified monoclonal antibody (mAb) is provided azide-free and lyophilized. It is available in two quantities:

**hpd11-mab1:** 100 µg Anti-hPD-L1-hIgG1

**hpd11-mab1-03:** 3 x 100 µg Anti-hPD-L1-hIgG1

**Target:** Human Programmed death-ligand 1

**Species reactivity:** Human and mouse

**Source:** CHO cells

**Isotype:** Human IgG1

**Light chain type:** Kappa

**Clonality:** Monoclonal

**Purification:** By affinity chromatography with protein G

**Formulation:** 0.2 µm filtered solution in a sodium phosphate buffer with glycine, saccharose, and stabilizing agents

## Storage and stability

- Product is shipped at room temperature. Upon receipt, store at -20 °C.
- Reconstituted antibody is stable for 1 month when stored at 4 °C and for 1 year when aliquoted and stored at -20 °C. Avoid repeated freeze-thaw cycles.

## Quality control

- Binding of Anti-hPD-L1-hIgG1 to surface expressed human PD-L1 on target cells has been validated using flow cytometry.
- Absence of bacterial contamination (e.g. lipoproteins and endotoxins) has been confirmed using HEK-Blue™ TLR2 and TLR4 cellular assays.

## PRODUCT DESCRIPTION

Anti-hPD-L1-hIgG1 is a recombinant monoclonal antibody (mAb) featuring a fully sequenced variable region equivalent to Atezolizumab, that recognizes human (h)PD-L1, and the constant region of the human (h)IgG1 isotype. Anti-hPD-L1-hIgG1 was generated by recombinant DNA technology, produced in CHO cells, and purified by affinity chromatography with protein G.

## PD-L1 background

Programmed cell death ligand 1 (PD-L1), also known as cluster of differentiation 274 (CD274) or B7 homolog 1 (B7-H1) is a transmembrane protein that can be constitutively expressed or induced in myeloid, lymphoid, and normal epithelial cells, as well as in cancer<sup>1,2</sup>. PD-L1 is the principle ligand for programmed cell death protein 1 (PD-1). This interaction is essential in the development of immune tolerance preventing excessive immune cell activity. However, PD-L1 expression is an immune evasion mechanism exploited by various malignancies<sup>3</sup>. Specifically, over-expressed PD-L1 on tumor cells and tumor infiltrating immune cells, such as macrophages, is able to bind to PD-1 on cytotoxic T cells, and ultimately inhibit the anti-tumor T cell response<sup>2, 4</sup>.

Thus, there are numerous PD-L1 inhibitors in development as promising immuno-oncology therapies. Notably, Atezolizumab (also known as MPDL3280A), a fully humanized IgG1 (N298A) mAb that blocks the interaction of PD-L1 with PD-1 and induces anti-tumor immune reactivation, has been approved by the FDA for combinational use in the treatment of lung and breast cancer<sup>2, 5</sup>.

## IgG1 isotype effector function

Human IgG1 binds with high affinity to the Fc receptor on phagocytic cells. Therefore, Anti-hPD-L1-hIgG1 displays high effector function, including antibody-dependent cell-mediated cytotoxicity (ADCC) and complement-dependent cytotoxicity (CDC).

**1. Juneja V.R. et al. 2017.** PD-L1 on tumor cells is sufficient for immune evasion in immunogenic tumors and inhibits CD8 T cell cytotoxicity. *J Exp Med* 214, 895-904. **2. Kythreotou A. et al. 2018.** PD-L1. *J Clin Pathol* 71, 189-194. **3. Sun C. et al. 2018.** Regulation and Function of the PD-L1 Checkpoint. *Immunity* 48, 434-452. **4. Lau J. et al. 2017.** Tumour and host cell PD-L1 is required to mediate suppression of anti-tumour immunity in mice. *Nat Commun* 8, 14572. **5. Heimes A.S. & Schmidt M. 2019.** Atezolizumab for the treatment of triple-negative breast cancer. *Expert Opin Investig Drugs* 28, 1-5.

## METHODS

### Anti-hPD-L1-hIgG1 resuspension (100 µg/ml)

*Note: Ensure you see the lyophilized pellet before resuspension.*

- Add 1 ml of sterile water to 100 µg and gently pipette until completely resuspended.
- Prepare aliquots and store at -20 °C until required.

## ANTIBODY ISOTYPE COLLECTION

For your research, InvivoGen provides an Anti-hPD-L1 isotype family. This isotype family consists of mAbs comprising a variable region equivalent to Atezolizumab and differing constant regions of both natural and engineered human isotypes (*see related products*). The isotypes differ in their functional and effector functions, such as antibody-dependent cell-mediated cytotoxicity (ADCC) and complement dependent cytotoxicity (CDC).

**Disclaimer:** The terms "Atezolizumab" and "MPDL3280A" are only used as references. Anti-hPD-L1-hIgG1 is not a pharmaceutical biosimilar of Atezolizumab. It has not been developed nor approved by Atezolizumab owner(s), and is not intended for any therapeutic or diagnostic use in human or animal.

## RELATED PRODUCTS

Product	Catalog Code
Anti-hPD-L1-hIgG1 (N298A)	hpd11-mab12
Anti-hPD-L1-hIgG1fut	hpd11-mab13
Raji-hPD-L1 Cells	raji-hpd11

## 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 3622-34-80

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