

PDL-2 antibody (pAb)

Rabbit Anti-Human/Mouse/Rat PDL-2

Instruction Manual

Catalog Number	PK-AB718-4063
Synonyms	PDL-2 Antibody: PDL-2, Programmed cell death 1 ligand-2, programmed death ligand 1, PD-L2, B7DC
Description	Cell-mediated immune responses are initiated by T lymphocytes that are themselves stimulated by cognate peptides bound to MHC molecules on antigen-presenting cells (APC). T-cell activation is generally self-limited as activated T cells express receptors such as PD-1 (also known as PDCD-1) that mediate inhibitory signals from the APC. PD-1 can bind two different but related ligands, PDL-1 and PDL-2, both of which are thought to act as a negative regulator of T cell activation. However, it has been suggested that PDL-2 can act to stimulate an immunogenic response through an alternative receptor from PD-1. At least three isoforms of PDL-2 are known to exist; this antibody is specific to the largest isoform. PDL-2 antibody has no cross-reactivity to PDL-1.
Quantity	100 µg
Source / Host	Rabbit
Immunogen	PDL-2 antibody was raised in rabbits against a 16 amino acid peptide from near the center of human PDL-2.
Purification Method	Affinity chromatography purified via peptide column.
Clone / IgG Subtype	Polyclonal antibody
Species Reactivity	Human, Mouse, Rat
Specificity	
Formulation	Antibody is supplied in PBS containing 0.02% sodium azide.
Reconstitution	During shipment, small volumes of antibody will occasionally become entrapped in the seal of the product vial. For products with volumes of 200 µl or less, we recommend gently tapping the vial on a hard surface or briefly centrifuging the vial in a tabletop centrifuge to dislodge any liquid in the container's cap.
Storage & Stability	Antibody can be stored at 4°C for three months and at -20°C for up to one year. As with all antibodies care should be taken to avoid repeated freeze thaw cycles. Antibodies should not be exposed to prolonged high temperatures.
Applications	E, WB, IHC, IF INote: Antibody might be suitable for other applications not tested so far. Optimal concentrations for each application have to be determined individually. PDL-2 antibody can be used for detection of PDL-2 by Western blot at 0.5 - 2 µg/mL. Antibody can also be used for immunohistochemistry starting at 2.5 µg/mL. For immunofluorescence start at 20 µg/mL.
Images	Available upon request.
References	Holling TM, Schooten E, and van Den Elsing PJ. Function and regulation of MHC class II molecules in T-lymphocytes: of mice and men. Hum. Immunol. 2004; 65:282-90. Ishida Y, Agata Y, Shibahara K, et al. Induced expression of PD-1, a novel member of the immunoglobulin gene superfamily, upon programmed cell death. EMBO J. 1992; 11:3887-95. LaGier J and Pober JS. Immune accessory functions of human endothelial cells are modulated by overexpression of B7-H1 (PDL1). Hum. Immunol. 2006; 67:568-78. Zhang Y, Chung Y, Bishop C, et al. Regulation of T cell activation and tolerance by PDL2. Proc. Natl. Acad. Sci. USA 2006; 103:11695-700.
Images	Available upon request.
Related Products	Cat.No. PK-AB718-4063P; PDL-2 Peptide Cat.No. PK-AB718-1207; Raji Cell Lysate

FOR IN VITRO RESEARCH USE ONLY. NOT FOR DIAGNOSTIC OR THERAPEUTIC PROCEDURES.