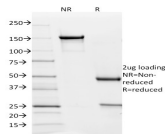


PD-L2 / PDCD1LG2 / CD273 Monoclonal Mouse Antibody (PDL2/1850)



Product Description

Recognizes a protein of about 31 kDa, which is identified as PD-L2 (same as PDCD1LG2). Engagement of CD28 by B7-1 (CD80) or B7-2 (CD86) in the presence of antigen promotes T cell proliferation, cytokine production, differentiation of effector T cells and the induction of Bcl-x, a promoter of T cell survival. Conversely, engagement of CTLA4 by B7-1 or B7-2 may inhibit proliferation and IL-2 production. PD-L2 does not bind CD28, cytotoxic T lymphocyte A4 or ICOS (inducible co-stimulator). The constitutive expression of PD-L1 and PD-L2 on parenchymal cells of heart, lung and kidney suggests that the Pdccl-1-Pdccl-L system could provide unique negative signaling to help prevent autoimmune disease.

Primary antibodies are available purified, or with a selection of fluorescent CF® dyes and other labels. CF® dyes offer exceptional brightness and photostability. See the [CF® Dye Brochure](#) for more information. Note: Conjugates of blue fluorescent dyes like CF®405S and CF®405M are not recommended for detecting low abundance targets, because blue dyes have lower fluorescence and can give higher non-specific background than other dye colors.

Stock status: Because Biotium offers a large number of antibody and conjugation options, primary antibody conjugates may be made to order. Typical lead times are up to one week for CF® dye and biotin conjugates, and up to 2-3 weeks for fluorescent protein and enzyme conjugates. Please email order@biotium.com to inquire about stock status and lead times before placing your order.

Catalog number key for antibody number 1850, Anti-CD273|PD-L2|PDCD1LG2 (PDL2/1850)

Product attributes

Antibody number	#1850
Antibody reactivity (target)	CD273, PD-L2, PDCD1LG2
Antibody type	Primary
Host species	Mouse
Clonality	Monoclonal
Clone	PDL2/1850
Isotype	IgG1, kappa
Molecular weight	31 kDa
Synonyms	B7 dendritic cell molecule; B7-DC; B7DC; Btdc; Butyrophilin B7DC; CD273; PD-L2; PDCD1L2; PDCD1LG2; PDL2; Programmed death ligand 2
Human gene symbol	PDCD1LG2
Entrez gene ID	80380
SwissProt	Q9Q51
Unigene	532279; 617872
Immunogen	Recombinant fragment (around aa 27-220) of human PD-L2 protein (exact sequence is proprietary)
Verified antibody applications	Flow (verified)
Antibody target cellular localization	Plasma membrane
Species reactivity	Human
Antibody application notes	Flow cytometry: 0.5-1 ug/million cells; Optimal dilution for a specific application should be determined., Higher concentration may be required for direct detection using primary antibody conjugates than for indirect detection with secondary antibody
Positive control	Placenta or Pancreas. Thyroid Carcinoma.
Shipping condition	Room temperature
Storage Conditions	Store at 2 to 8 °C, Protect fluorescent conjugates from light, Note: store BSA-free antibodies at -10 to -35 °C
Regulatory status	For research use only (RUO)
Antibody/conjugate formulation	Conjugates: 0.1 mg/mL in PBS/0.1% BSA/0.05% azide, HRP conjugates: 0.1 mg/mL in PBS/0.05% BSA, Purified: 0.2 mg/mL in PBS/0.05% BSA/0.05% azide, Purified, BSA-free: 1 mg/mL in PBS without azide
Shelf life	Guaranteed for at least 24 months from date of receipt when stored as recommended
Product origin	Product may contain either bovine serum albumin (BSA) from bovine serum (Bos taurus), or recombinant BSA produced in Chinese hamster ovary cells. Inquire for the specific lot.

Antibody # prefix	Conjugation	Ex/Em (nm)	Laser line	Detection channel	Dye Features
BNC04	CF®405S	404/431	405	DAPI (microscopy), AF405	CF®405S Features
BNC88	CF®488A	490/515	488	GFP, FITC	CF®488A Features
BNC68	CF®568	562/583	532, 561	RFP, TRITC	CF®568 Features
BNC94	CF®594	593/614	561	Texas Red®	CF®594 Features
BNC40	CF®640R	642/662	633-640	Cy®5	CF®640R Features
BNC47	CF®647	650/665	633-640	Cy®5	CF®647 Features
BNC74	CF®740	742/767	633-685	775/50	CF®740 Features
BNCB	Biotin	N/A	N/A	N/A	
BNUB	Purified	N/A	N/A	N/A	
BNUM	Purified, BSA-free	N/A	N/A	N/A	

Alexa Fluor, Pacific Blue, Pacific Orange, and Texas Red are trademarks or registered trademarks of Thermo Fisher Scientific; Cy is a registered trademark of Cytiva; IRDye, LI-COR, and Odyssey are registered trademarks of LI-COR Bioscience.

This datasheet was generated on August 7, 2025 at 06:23:45 PM. Visit product page to check for updated information before use.

Product link: <https://biotium.com/product/pd-l2-pdccl1lg2-cd273-monoclonal-mouse-antibody-pdl2-1850/>