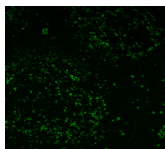


## PD1 / PDCD1 / CD279 Monoclonal Mouse Antibody (PDCD1/922)



### Product Description

PD-1 (also known as Programmed Death 1, CD279, and PDCD1) is a negative regulatory T cell surface receptor. PD-1 is a Type I transmembrane protein expressed on the plasma membrane of T-cells. It is a T-cell checkpoint protein, involved in preventing autoimmunity. Because binding of PD-1 by its ligands PD-L1 or PD-L2 has a suppressive effect on the immune system, inhibition of the PD-1/PD-L1 pathway is a major strategy of immune-oncology therapies. PD-L1 is upregulated in certain cancers, and blocking PD-1 or PD-L1 can inhibit cancer growth.

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](mailto:order@biotium.com) to inquire about stock status and lead times before placing your order.

**Catalog number key for antibody number 0922, Anti-CD279|PD1|PDCD1 (PDCD1/922)**

Call us : [800-304-5357](tel:800-304-5357)

### Product attributes

Antibody number	#0922
Antibody reactivity (target)	CD279, PD1, PDCD1
Antibody type	Primary
Host species	Mouse
Clonality	Monoclonal
Clone	PDCD1/922
Isotype	IgG1, kappa
Molecular weight	55 kDa
Synonyms	CD279; hPD-1; hSLE1; PD1; PDCD1; Programmed Cell Death Protein 1; Protein PD-1; SLEB2; Systemic lupus erythematosus susceptibility 2
Human gene symbol	PDCD1
Entrez gene ID	5133
SwissProt	Q15116
Unigene	158297
Immunogen	Recombinant full-length human PDCD1 protein
Antibody target cellular localization	Plasma membrane
Verified antibody applications	Flow (verified), IF (verified), IHC (FFPE) (verified)
Species reactivity	Human
Antibody application notes	Higher concentration may be required for direct detection using primary antibody conjugates than for indirect detection with secondary antibody. Immunofluorescence: 1-2 ug/mL. Immunohistology formalin-fixed 0.5-1 ug/mL. Staining of formalin-fixed tissues requires boiling tissue sections in 10 mM Tris with 1 mM EDTA, pH 9.0, for 10-20 min followed by cooling at RT for 20 minutes. Flow Cytometry 0.5-1 ug/million cells/0.1 mL. Optimal dilution for a specific application should be determined by user
Positive control	TY cells or Tonsil.
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
Shelf life	Guaranteed for at least 24 months from date of receipt when stored as recommended
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
Cell/tissue expression	B-cells, Myeloid cells, T-cells
Antibody research areas	Apoptosis, Cancer, Cancer immunotherapy/immune checkpoint
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.
Tumor expression	Leukemia/lymphoma

Antibody # prefix	Conjugation	Ex/Em (nm)	Laser line	Detection channel	Dye Features
BNC04	CF®405S	404/431	405	DAPI (microscopy), AF405	<a href="#">CF®405S Features</a>
BNC88	CF®488A	490/515	488	GFP, FITC	<a href="#">CF®488A Features</a>
BNC68	CF®568	562/583	532, 561	RFP, TRITC	<a href="#">CF®568 Features</a>
BNC94	CF®594	593/614	561	Texas Red®	<a href="#">CF®594 Features</a>
BNC40	CF®640R	642/662	633-640	Cy®5	<a href="#">CF®640R Features</a>
BNC47	CF®647	650/665	633-640	Cy®5	<a href="#">CF®647 Features</a>
BNC74	CF®740	742/767	633-685	775/50	<a href="#">CF®740 Features</a>
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 December 13, 2025 at 06:20:28 AM. Visit product page to check for updated information before use.

Product link: <https://biotium.com/product/monoclonal-anti-pd1-pdcd1-cd279-pdcd1922/>