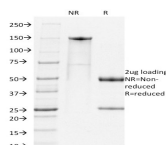


PD1 / PDCD1 / CD279 Mouse Monoclonal Rat Antibody (RMP1-14)



Product Description

PD-1 is expressed on a subset of CD4-CD8-thymocytes, and on activated T and B cells. PD-1 is thought to be involved in lymphocyte clonal selection and peripheral tolerance. The PD-1 ligands, PD-L1 (also known as B7-H1) and PD-L2 (B7-DC), are members of the B7 immunoglobulin superfamily. PD-1 is transiently expressed on CD4 and CD8 thymocytes as well as activated T and B-lymphocytes and myeloid cells. Additionally, Pdc1mRNA is expressed in developing B-lymphocytes during the pro-B-cell stage. The RMP1-14 antibody has been reported to block the binding of PD-1 to its ligands (B7-H1 and B7-DC) and to inhibit T cell proliferation and cytokine production co-stimulated by macrophages (but not by dendritic cells and B cells).

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 2002, Anti-PD1 (RMP1-14)

Product attributes

Antibody number	#2002
Antibody reactivity (target)	PD1
Antibody type	Primary
Host species	Rat
Clonality	Monoclonal
Clone	RMP1-14
Isotype	IgG2a, kappa
Molecular weight	~50-55 kDa
Synonyms	CD279; CD279 antigen; PD1; PDCD1; Programmed cell death 1; Programmed cell death protein 1; Systemic lupus erythematosus susceptibility 2 (SLEB2)
Entrez gene ID	18566
SwissProt	Q02242
Unigene	5024
Immunogen	PDCD1 transfected BHK cells
Antibody target cellular localization	Plasma membrane
Expected antibody applications	Flow, surface (published for clone), Functional studies (published for clone)
Species reactivity	Mouse
Positive control	Spleen, Thymus or Lymph Node.
Shipping condition	Room temperature
Storage Conditions	Note: store BSA-free antibodies at -10 to -35 °C, Store at 2 to 8 °C, Protect fluorescent conjugates from light
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.
Cell/tissue expression	B-cells, Myeloid cells, T-cells
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	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.

References

Note: References for this clone sold by other suppliers may be listed for expected applications.

1. J Immunol (2005) 175:1586-1592. (Flow; functional studies)
2. Cancer Res (2014) 74(14): 3652-8. (functional studies)
3. Cancer Discov (2017) 7(7): 694-703. (functional studies)

This datasheet was generated on July 28, 2025 at 05:29:49 PM. Visit product page to check for updated information before use.

Product link: <https://biotium.com/product/pd1-pdc1-cd279-mouse-monoclonal-rat-antibody-rmp1-14/>