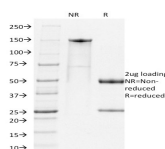


# PD1 / PDCD1 / CD279 Mouse Monoclonal Rat Antibody (RMP1-14)



## 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](mailto: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

<b>Antibody number</b>	#2002
<b>Antibody reactivity (target)</b>	PD1
<b>Antibody type</b>	Primary
<b>Host species</b>	Rat
<b>Clonality</b>	Monoclonal
<b>Clone</b>	RMP1-14
<b>Isotype</b>	IgG2a, kappa
<b>Molecular weight of antigen</b>	~50-55 kDa
<b>Synonyms</b>	CD279; CD279 antigen; PD1; PDCD1; Programmed cell death 1; Programmed cell death protein 1; Systemic lupus erythematosus susceptibility 2 (SLEB2)
<b>Entrez gene ID</b>	18566
<b>SwissProt</b>	Q02242
<b>Unigene</b>	5024
<b>Immunogen</b>	PDCD1 transfected BHK cells
<b>Antibody target cellular localization</b>	Plasma membrane
<b>Expected antibody applications</b>	Flow, surface (published for clone), Functional studies (published for clone)
<b>Species reactivity</b>	Mouse
<b>Positive control</b>	Spleen, Thymus or Lymph Node.
<b>Shipping condition</b>	Room temperature
<b>Storage Conditions</b>	Note: store BSA-free antibodies at -10 to -35 °C, Store at 2 to 8 °C, Protect fluorescent conjugates from light
<b>Regulatory status</b>	For research use only (RUO)
<b>Antibody/conjugate formulation</b>	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
<b>Shelf life</b>	Guaranteed for at least 24 months from date of receipt when stored as recommended
<b>Product origin</b>	Product may contain either bovine serum albumin (BSA) from bovine serum ( <i>Bos taurus</i> ), or recombinant BSA produced in Chinese hamster ovary cells. Inquire for the specific lot.
<b>Cell/tissue expression</b>	B-cells, Myeloid cells, T-cells
<b>Tumor expression</b>	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.

## References

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

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

This datasheet was generated on June 4, 2026 at 05:54:25 PM. Visit product page to check for updated information before use.

Product link: <https://biotium.com/product/pd1-pdcd1-cd279-mouse-monoclonal-rat-antibody-rmp1-14/>