

CD273 (B7-DC) Rat Anti-Mouse mAb (clone TY25), PE Conjugate

Store at 2°C to 8°C

Pub. No. MAN0009524 Rev. 1.00

Catalog No.	Form	Amount	Excitation	Peak Emission
A18421	PE	50 µg (0.2 mg/mL)	496 nm	578 nm

Clone	TY25
Host/Class	Rat IgG2aκ
Description	CD273 (B7-DC) is a ~25 kDa, a member of the B7 subfamily of the Ig superfamily. Mouse CD273 has a short cytoplasmic tail (4 amino acids), and does not bind CD28/CTLA-4, though it does bind PD-1. The interaction between PD-1 and CD273 have been reported to be involved in co-stimulation or suppression of T cell proliferation depending on state of cellular activation. CD273 is primarily expressed by sub-populations of dendritic cells, monocytes and macrophages.
Alternate Names	B7DC, PD-L2, PDL-2, PDL2
Applications*	FC (mouse splenocyte suspensions and B7-DC-transfected cells), FUNC ⁸
Storage Buffer	The reagent is provided in aqueous buffer with 0.09% sodium azide, and may contain carrier protein/stabilizer. CAUTION! Sodium azide is extremely toxic and may react with lead and copper plumbing to form highly explosive metal azides. Properly dispose of solutions containing sodium azide. Read the Safety Data Sheet (SDS) and follow the handling instructions. Wear appropriate protective eyewear, clothing, and gloves. SDSs are available at www.lifetechnologies.com/support .
Storage	Store reagents in the dark at 2° to 8°C. Do not freeze. If the reagent is being diluted, it is recommended that only the quantity to be used within one week be diluted. Avoid prolonged light exposure with fluorochrome-conjugated antibodies. Use dim light during handling, incubation with cells, and prior to analysis.
Stability	When stored as instructed, expires one year from date of receipt unless otherwise indicated on Certificate of Analysis.
Lot Number	See product label.
References	<ol style="list-style-type: none"> 1. Aramaki O, Shirasugi N, Takayama T, Shimazu M, Kitajima M, Ikeda Y, Azuma M, Okumura K, Yagita H, Niimi M. 2004. Programmed death-1-programmed death-L1 interaction is essential for induction of regulatory cells by intratracheal delivery of alloantigen. <i>Transplantation</i>. Jan 15;77(1):6-12. 2. Tseng SY, Otsuji M, et al. 2001. B7-DC, a new dendritic cell molecule with potent costimulatory properties for T cells. <i>J Exp Med</i>. 193(7): 839-46. 3. Carter L, Fouser L, et al. 2002. PD-1:PD-L inhibitory pathway affects both CD4(+) and CD8(+) T cells and is overcome by IL-2. <i>Eur J Immunol</i>. 32(3): 634-43. 4. Latchman Y, Wood CR, et al. 2001. PD-L2 is a second ligand for PD-1 and inhibits T cell activation. <i>Nat Immunol</i>. 2(3): 261-8. 5. Yamazaki T, Akiba H, et al. Expression of programmed death 1 ligands by murine T cells and APC. <i>J Immunol</i>. 2002 Nov 15;169(10):5538-45.

* Because conditions may vary, it is recommended that each investigator determine the optimal amount of antibody to be used for each application.

FC = flow cytometry; FUNC = functional assay; ICC = immunocytochemistry; IHC(F) = immunohistochemistry (frozen sample); IHC(P) = immunohistochemistry (paraffin embedded sample); IP = immunoprecipitation; RIA = radioimmunoassay; WB = western blot








For Research Use Only. Not for use in diagnostic procedures.

Manufacturing site: 7335 Executive Way | Frederick, MD 21704 | Toll Free in USA 800.955.6288

References, continued	6. Kanai T, Totsuka T, Uraushihara K, Makita S, Nakamura T, Koganei K, Fukushima T, Akiba H, Yagita H, Okumura K, Machida U, Iwai H, Azuma M, Chen L, Watanabe M. 2003. Blockade of B7-H1 suppresses the development of chronic intestinal inflammation. <i>J Immunol.</i> 171(8):4156-63.
	7. Mohammed Javeed I. Ansari, Alan D. Salama, Tanuja Chitnis, R. Neal Smith, Hideo Yagita, Hisaya Akiba, Tomohide Yamazaki, Miyuki Azuma, Hideyuki Iwai, Samia J. Khoury, Hugh Auchincloss, Jr. and Mohamed H. Sayegh. 2003. The Programmed Death-1 (PD-1) Pathway Regulates Autoimmune Diabetes in Nonobese Diabetic (NOD) Mice. <i>J Exp Med.</i> 198 (1): 63-69.
	8. Tanaka K., M. Albin et al. 2007. PDL1 is required for peripheral transplantation tolerance and protection from chronic allograft rejection. <i>J Immunol.</i> 179(8):5204-5210.

Explanation of Symbols

The symbols present on the product label are explained below:

Symbol	Description	Symbol	Description	Symbol	Description
	Manufacturer		Catalog number		Batch code
	Use by		Temperature limitation		
	Consult instructions for use		Caution, consult accompanying documents		

Important Licensing Information

This product may be covered by one or more Limited Use Label Licenses. By use of this product, you accept the terms and conditions of all applicable Limited Use Label Licenses.

Limited Product Warranty

Life Technologies Corporation and/or its affiliate(s) warrant their products as set forth in the Life Technologies' General Terms and Conditions of Sale found on Life Technologies' website at www.lifetechnologies.com/termsandconditions. If you have any questions, please contact Life Technologies at www.lifetechnologies.com/support.

© 2013 Life Technologies Corporation. All rights reserved. The trademarks mentioned herein are the property of Life Technologies Corporation and/or its affiliate(s) or their respective owners.

DISCLAIMER: LIFE TECHNOLOGIES CORPORATION AND/OR ITS AFFILIATE(S) DISCLAIM ALL WARRANTIES WITH RESPECT TO THIS DOCUMENT, EXPRESSED OR IMPLIED, INCLUDING BUT NOT LIMITED TO THOSE OF MERCHANTABILITY, FITNESS FOR A PARTICULAR PURPOSE, OR NON-INFRINGEMENT. TO THE EXTENT ALLOWED BY LAW, IN NO EVENT SHALL LIFE TECHNOLOGIES AND/OR ITS AFFILIATE(S) BE LIABLE, WHETHER IN CONTRACT, TORT, WARRANTY, OR UNDER ANY STATUTE OR ON ANY OTHER BASIS FOR SPECIAL, INCIDENTAL, INDIRECT, PUNITIVE, MULTIPLE OR CONSEQUENTIAL DAMAGES IN CONNECTION WITH OR ARISING FROM THIS DOCUMENT, INCLUDING BUT NOT LIMITED TO THE USE THEREOF.

For support visit www.lifetechnologies.com/support or email techsupport@lifetech.com

www.lifetechnologies.com

15 November 2013

