

## CD278 (ICOS) Hamster Anti-Mouse mAb (clone 15F9), PE Conjugate

Store at 2°C to 8°C

Pub. No. MAN0009538 Rev. 1.00

Catalog No.	Form	Amount	Excitation	Peak Emission
A18435	PE	50 µg (0.2 mg/mL)	496 nm	578 nm

Clone	15F9
Host/Class	Golden Syrian Hamster IgG
Description	The CD278 (ICOS) Hamster Anti-Mouse Monoclonal Antibody (mAb) recognizes mouse CD278, a T cell specific molecule and CD28/CTLA-4 family member. CD278 forms a homodimer of 47-57 kDa and is expressed on activated T cells. It is required for the humoral immune response and has potent costimulatory activity for T cell activation and proliferation. ICOS binds to its ligand on activated APC and function in inflammatory autoimmune response. The mAb has functional activity and may be involved in Th2 cell development.
Alternate Names	ICOS, Inducible CoStimulatory molecule
Applications*	FC (mouse ConA-activated splenocytes), FUNC <sup>2</sup>
Storage Buffer	The reagent is provided in aqueous buffer with 0.09% sodium azide, and may contain carrier protein/stabilizer. <b>CAUTION!</b> 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 <a href="http://www.lifetechnologies.com/support">www.lifetechnologies.com/support</a> .
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"> <li>1. Lucy S. K. Walker, Helen E. Wiggett, Fabrina M. C. Gaspal, Chandra R. Raykundalia, Margaret D. Goodall, Kai-Michael Toellner, and Peter J. L. Lane. 2003. Established T Cell-Driven Germinal Center B Cell Proliferation Is Independent of CD28 Signaling but Is Tightly Regulated Through CTLA-4. <i>J Immunol.</i> 170:91-98.</li> <li>2. Kohyama M, Sugahara D, Sugiyama S, Yagita H, Okumura K, Hozumi N. Inducible costimulator-dependent IL-10 production by regulatory T cells specific for self-antigen. <i>Proc Natl Acad Sci U S A.</i> 2004 Mar 23;101(12):4192-7.</li> <li>3. McAdam AJ, Chang TT, Lumelsky AE, Greenfield EA, Boussiotis VA, Duke-Cohan JS, Chernova T, Malenkovich N, Jabs C, Kuchroo VK, Ling V, Collins M, Sharpe AH, Freeman GJ. Mouse inducible costimulatory molecule (ICOS) expression is enhanced by CD28 costimulation and regulates differentiation of CD4+ T cells. <i>J Immunol.</i> 2000 Nov 1;165(9):5035-40.</li> </ol>

\* 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

## 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](http://www.lifetechnologies.com/termsandconditions). If you have any questions, please contact Life Technologies at [www.lifetechnologies.com/support](http://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](http://www.lifetechnologies.com/support) or email [techsupport@lifetech.com](mailto:techsupport@lifetech.com)

[www.lifetechnologies.com](http://www.lifetechnologies.com)

15 November 2013

