

CD147 Rat Anti-Mouse mAb (clone RL73), PE Conjugate

Store at 2°C to 8°C

Pub. No. MAN0009459 Rev. 1.00

Catalog No.	Form	Amount	Excitation	Peak Emission
A18354	PE	50 µg (0.2 mg/mL)	496 nm	578 nm

Clone	RL73
Host/Class	Rat IgG2aκ
Description	The CD147 Rat Anti-Mouse Monoclonal Antibody (mAb) recognizes mouse CD147, which is expressed by a majority of thymocytes, activated peripheral T cells, erythroblasts, and erythrocytes. CD147 expression correlates with cycling of immature thymocytes and inhibition of mature T cell development on immature fetal thymocytes.
Alternate Names	Basigin, EMMPRIN
Applications*	FC (mouse thymocyte cell suspensions)
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"> MacDonald HR, Lees RK, Bron C. 1985. Cell surface glycoproteins involved in the stimulation of interleukin 1-dependent interleukin 2 production by a subline of EL4 thymoma cells. I. Functional characterization by monoclonal antibodies. <i>J Immunol.</i> 135(6):3944-50. Renno T, Wilson A, Dunkel C, Coste I, Maisnier-Patin K, Benoit de Coignac A, Aubry JP, Lees RK, Bonnefoy JY, MacDonald HR, Gauchat JF. 2002. A role for CD147 in thymic development. <i>J Immunol.</i> 168(10):4946-50. Coste I, Gauchat JF, Wilson A, Izui S, Jeannin P, Delneste Y, MacDonald HR, Bonnefoy JY, Renno T. 2001. Unavailability of CD147 leads to selective erythrocyte trapping in the spleen. <i>Blood.</i> 97(12):3984-8.

* 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		

Limited Use Label License: Research Use Only

The purchase of this product conveys to the purchaser the limited, non-transferable right to use the purchased amount of the product only to perform internal research for the sole benefit of the purchaser. No right to resell this product or any of its components is conveyed expressly, by implication, or by estoppel. This product is for internal research purposes only and is not for use in commercial applications of any kind, including, without limitation, quality control and commercial services such as reporting the results of purchaser's activities for a fee or other form of consideration. For information on obtaining additional rights, please contact outlicensing@lifetech.com or Out Licensing, Life Technologies, 5791 Van Allen Way, Carlsbad, California 92008.

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 October 2013

