

CD8a Mouse Anti-Human Monoclonal Antibody (clone RPA-T8) PE Conjugate

Store at 2°C to 8°C

Pub. No. MAN0008180 Rev. 1.00

Catalog No.	Form	Amount	Excitation	Peak Emission
A15827	PE	25 tests (5 µL/test, 0.025 µg/µL)	496 nm	578 nm

Clone	RPA-T8
Host/Class	Mouse IgG _{1κ}
Description	The CD8a Mouse Anti-Human Monoclonal Antibody reacts with human CD8a, a cell surface receptor of 32–34 kDa. CD8a can exist as an α/β heterodimer with the CD8 beta chain, or as an α/α homodimer. CD8 binds MHC class I antigen, is involved in T-cell development, and activation of mature T cells. CD8a is expressed on most thymocytes and a subpopulation of mature T cells and NK cells.
Alternate Names	CD8 alpha, leu-2a
Applications*	FC (human peripheral blood leukocytes)
Storage Buffer	The reagent is provided in aqueous buffer with 0.09% sodium azide, and may contain carrier protein/stabilizer.
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"> Knapp, W., B. Dorken, et al. eds. (1989). <i>Leucocyte Typing IV: White Cell Differentiation Antigens</i>. Oxford University Press. New York. Schlossman, S., L. Bloumsell, et al. eds. (1995). <i>Leucocyte Typing V: White Cell Differentiation Antigens</i>. Oxford University Press. New York. Kishimoto, T., A.E.G., von dem Borne, et al. eds. (1998). <i>Leucocyte Typing VI: White Cell Differentiation Antigens</i>. Garland Publishing, Inc. London.

* 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

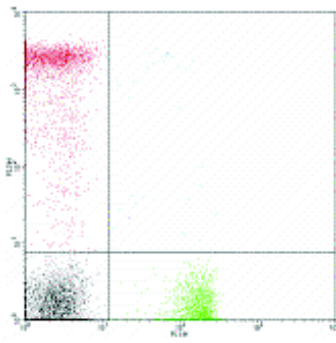


Figure 1 Staining of normal human peripheral blood cells with CD8a Mouse Anti-Human PE Conjugate (Cat. no. A15827). Staining of normal human peripheral blood cells with CD8a Mouse Anti-Human PE Conjugate and a human CD4 FITC conjugate.

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 affiliates 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

24 April 2013

