

CD181 (CXCR1) Mouse Anti-Human mAb (clone eBio8F1-1-4), FITC Conjugate

Store at 2°C to 8°C

Pub. No. MAN0009447 Rev. 1.00

Catalog No.	Form	Amount	Excitation	Peak Emission
A18342	FITC	25 tests (5 µL/test; 0.1 µg/µL)	488 nm	519 nm

Clone	eBio8F1-1-4 (8F1-1-4)
Host/Class	Mouse IgG2bk
Description	The CD181 (CXCR1) Mouse Anti-Human Monoclonal Antibody (mAb) reacts with human CD181 (CXCR1). CD181 is a 67–70 kDa member of the 7-transmembrane spanning G-protein coupled receptor (GPCR) family. It is involved in cell activation, chemotaxis, proliferation and angiogenesis. Its ligands, include IL-8, NAP-2, GCP-2 and GRO-alpha. CD181 is expressed as a homodimer, or a heterodimer with CD182 (CXCR2, IL-8R beta) and is expressed on granulocytes, NK cells, a subset of T cells, mast cells, monocytes, endothelial cells, megakaryocytes, and oligodendrocytes.
Alternate Names	Interleukin-8 receptor A, IL-8Ra, IL8RA, IL-8R alpha
Applications*	FC (normal human peripheral blood)
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. Wolf M, Delgado MB, Jones SA, Dewald B, Clark-Lewis I, Baggiolini M. Granulocyte chemotactic protein 2 acts via both IL-8 receptors, CXCR1 and CXCR2. <i>Eur J Immunol.</i> 1998 Jan;28(1):164-70. 2. Lippert U, Artuc M, Grutzkau A, Moller A, Kenderessy-Szabo A, Schadendorf D, Norgauer J, Hartmann K, Schweitzer-Stenner R, Zuberbier T, Henz BM, Kruger-Krasagakes S. Expression and functional activity of the IL-8 receptor type CXCR1 and CXCR2 on human mast cells. <i>J Immunol.</i> 1998 Sep 1;161(5):2600-8. 3. Li F, Zhang X, Gordon JR. CXCL8((3-73))K11R/G31P antagonizes ligand binding to the neutrophil CXCR1 and CXCR2 receptors and cellular responses to CXCL8/IL-8. <i>Biochem Biophys Res Commun.</i> 2002 May 10;293(3):939-44. 4. Sprenger H, Lloyd AR, Meyer RG, Johnston JA, Kelvin DJ. Genomic structure, characterization, and identification of the promoter of the human IL-8 receptor A gene. <i>J Immunol.</i> 1994 Sep 15;153(6):2524-32. 5. Chuntharapai A, Lee J, Hebert CA, Kim KJ. Monoclonal antibodies detect different distribution patterns of IL-8 receptor A and IL-8 receptor B on human peripheral blood leukocytes. <i>J Immunol.</i> 1994 Dec 15;153(12):5682-8.

* 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

