

CD86 (B7-2) Rat Anti-Mouse mAb FITC Conjugate

Store at 2°C to 8°C

Catalog Number A18658

Pub. No. MAN0009008 Rev. 2.0

Catalog No.	Form	Amount	Excitation	Peak Emission
A18658	FITC	25 µg	488 nm	519 nm
A27076	FITC	100 µg	488 nm	519 nm
A27077	FITC	500 µg	488 nm	519 nm

Product description

The CD86 (B7-2) Rat Anti-Mouse Monoclonal Antibody (mAb) reacts with mouse CD86, also known as B7-2, an 80 kDa cell surface protein which is a ligand for CD28, a co-stimulatory receptor for the T cell receptor (TCR). CD28 can also bind a second B7 ligand known as CD80 (B7-1). Both CD80 and CD86 are expressed on activated B cells and antigen-presenting cells. These ligands trigger CD28 signaling in concert with TCR activation to drive T cell proliferation, induce high-level expression of IL-2, impart resistance to apoptosis, and enhance T cell cytotoxicity. The interaction between the B7 ligands and CD28 provides crucial communication between T cells and B cells or APCs to coordinate the adaptive immune response. The GL-1 antibody may be used as a marker for CD86 expression on B cells, macrophages, and dendritic cells.

Product specifications

Clonality:	Monoclonal
Host/Class:	Rat IgG
Reactivity:	Mouse CD86
Clone/PAD:	GL-1 (GL1)
Isotype:	IgG2ak
Lot:	See product label

Product applications

Applications reported for the CD86 (B7-2) Rat Anti-Mouse mAb include Flow Cytometry.

Storage and handling

Store reagents at 2°C to 8°C. If the reagent is being diluted, it is recommended that only the quantity to be used within one week be diluted. Cells should be analyzed within 18 hours of staining for best results.

Avoid light exposure with fluorochrome-conjugated antibodies. Use dim light during handling, incubation with cells, and prior to analysis.

Stability

When stored as instructed, expires six months from date of receipt unless otherwise indicated on the Certificate of Analysis.

Storage buffer

Phosphate buffered saline (PBS) with 0.1% sodium azide.



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.

Product documentation

To obtain a Certificate of Analysis or Safety Data Sheet (SDS), visit <http://www.lifetechnologies.com/support>.

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.

Related products

Product Name	Quantity	Catalog No.
AbC™ Anti-Mouse Bead Kit	1 kit	A10344
AbC™ Anti-Rat/Hamster Bead Kit	1 kit	A10389
FIX & PERM® Reagents (200 tests)	1 kit	GAS004

Product Name	Quantity	Catalog No.
Blue (UV excitation)	1 kit	L23105
Violet (405 nm excitation)	(200 assays)	L34955
Aqua (405 nm excitation)		L34957
Yellow (405 nm excitation)		L34959
Green (488 nm excitation)		L23101
Red (488 nm excitation)		L23102
Far-red (633/635 nm excitation)		L10210
Near-IR (633/635 nm excitation)		L10119

References

1. Liu Z, Geboes K, Hellings P, Maerten P, Heremans H, Vandenberghe P, Boon L, van Kooten P, Rutgeerts P, and Ceuppens JL. 2011. J. Immunol. 167: 1830-1838. (in vivo blocking, Immunohistochemistry – OCT embedded frozen tissue)
2. Kastenmuller W, Gasteiger G, Subramanian N, Sparwasser T, Busch DH, Belkaid Y, Drexler I, and Germain RN, 2011. J. Immunol. 187: 3186-3197. (in vivo blocking)
3. Zheng SG, Wang JH, Stohl, W, Kim KS, Gray JD, and Horwitz DA. 2006. J. Immunol. 176:3321-3329. (in vitro blocking)
4. Leithauser F, Meinhardt-Krajina T, Fink K, Wotschke B, Moller P and Reimann J. 2006. Am. J. Pathol. 168(6): 1898-1909. (Immunohistochemistry – frozen tissue)
5. Odobasic D, Kitching AR, Semple TJ, Timoshanko JR, Tipping PG, and Holdsworth SR. 2005. J. Am. Soc. Nephrol. 16: 2012-2022. (in vivo activation, Immunofluorescence microscopy – frozen tissue, Immunohistochemistry – frozen tissue)
6. Lenschow DJ, Ho SC, Sattar H, Rhee L, Gray G, Nabavi N, Herold KC, and Bluestone JA. 1995. J. Exp. Med. 181:1145-155. (in vitro blocking)
7. Blazar BR, Taylor PA, Panoskaltsis-Mortari A, Gray GS, and Vallera DA. 1995. Blood. 85: 2607-2618. (Immunohistochemistry – OCT embedded frozen tissue)

Explanation of symbols

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

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.

©2014 Thermo Fisher Scientific Inc. All rights reserved. All trademarks are the property of Thermo Fisher Scientific and its subsidiaries unless otherwise specified. FIX & PERM is a registered trademark of An Der Grub Bio Research GmbH.

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

lifetechnologies.com

12 November 2014

