

CD80 Hamster Anti-Mouse

Publication Number: MAN0007072

Rev. 1.00

Store at 2°C to 8°C

Catalog No.	Form	Amount	Excitation	Peak Emission
A14896	Pacific Blue™	25 µg	405 nm	455 nm
A14895	PerCP-Cy®5.5	25 µg	482 nm	695 nm

Product Description

CD80 is a highly glycosylated member of the Ig superfamily. It is constitutively expressed on dendritic cells, monocytes, and macrophages, and inducibly expressed on activated B and T cells. CD80 plays a co-stimulatory role with CD28 in T cell activation, proliferation, and cytokine production. It also plays an inhibitory role in T cells, in conjunction with CD152 (CTLA-4).

Product Specifications

Clonality:	Monoclonal
Host/Class:	Armenian Hamster IgG
Reactivity:	Mouse CD80
Immunogen:	Recombinant mouse CD80
Alternate Names:	B7-1, B7, Ly-53
Apparent MW:	60 kDa
Gene ID:	12519
Clone/PAD:	16-10A1
Isotype:	IgG
Lot:	See product label

Product Applications

Applications reported for the CD80 Hamster Anti-Mouse mAb include flow cytometry.¹

It is suggested that investigators titrate reagents to determine optimal conditions for use in their systems.

Stability

When stored as instructed, expires six months from date of receipt unless otherwise indicated on product label.

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.

Storage Buffer

Phosphate-buffered solution, pH 7.2, with 0.09% 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 Sheets (SDSs) and follow the handling instructions. Wear appropriate protective eyewear, clothing, and gloves. SDSs are available at www.lifetechnologies.com/support.

For research use only. Not for human or animal therapeutic or diagnostic use.

7335 Executive Way • Frederick • MD 21704 • E-mail: techsupport@lifetech.com

References

1. Andoniu CE, et al. 2005. *Nature Immunology* 6:1011.

Product Documentation

To obtain a Certificate of Analysis or Safety Data Sheets (SDSs), visit 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
Protein A Agarose	5 mL	15918-014
Recombinant Protein G (rProtein G) Agarose	5 mL	15920-010

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.

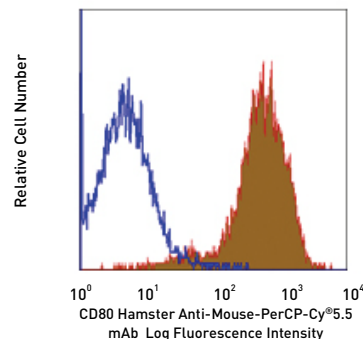


Figure 1 Thioglycolate-elicited Balb/c mouse peritoneal macrophages stained with CD80 Hamster Anti-Mouse-PerCP-Cy[®]5.5 Monoclonal Antibody (Cat. no. A14895).

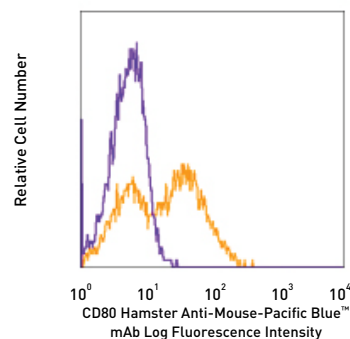


Figure 2 C57BL/6 splenocytes stimulated (4 days) with anti-mouse IgM and CD40, stained with CD80 Hamster Anti-Mouse-Pacific Blue™ Monoclonal Antibody (Cat. no. A14896).

Note: All flow cytometric data shown may not necessarily have been generated using the enclosed lot of reagent. For this reason, and due to differences in flow cytometers and cytometer settings, results may vary from those illustrated above.

Manufactured for Life Technologies by BioLegend, Inc.



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		

LIFE TECHNOLOGIES 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.

©2012 Life Technologies Corporation. All rights reserved. The trademarks mentioned herein are the property of Life Technologies Corporation or their respective owners.

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

www.lifetechnologies.com

15 September 2012

