

IFN gamma Rat Anti-Mouse mAb PE-Cyanine7 Conjugate

Store at 2°C to 8°C

Catalog Number A18713

Pub. No. MAN0009080 Rev. 2.0

Catalog No.	Form	Amount	Excitation	Peak Emission
A18713	PE-Cyanine7	25 µg	488 nm	767 nm

Product description

The IFN gamma Rat Anti-Mouse Monoclonal Antibody (mAb) is specific for mouse Interferon-gamma (IFN-g), a 20 kDa type II cytokine known for its central roles in protection against bacterial or viral pathogens and for its anti-tumor properties. IFN-g is secreted by several types of immune cells, which allow the cytokine to modulate innate immunity, when secreted by NK and NKT cells, and to function in support adaptive immunity when secreted by Th1 and CD8+ T cells (CTLs). The XMG1.2 clone is suitable for detection of intracellular IFN-g protein, e.g. by flow cytometry, as well as for quantitative analysis of the secreted protein by ELISA. This antibody is also widely used for neutralization of the functional activity of IFN-g in a variety of assays.

Product specifications

Clonality:	Monoclonal
Host/Class:	Rat IgG
Reactivity:	Mouse IFN
Clone/PAD:	XMG1.2
Isotype:	IgG1k
Lot:	See product label

Product applications

Applications reported for the IFN gamma Rat Anti-Mouse mAb include flow cytometry⁶, functional assays^{1, 2, 5}, ELISA^{3, 4}, immunohistochemistry (frozen tissue)⁷, and immunofluorescence microscopy⁸.

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. Choudhry N, Petry F, van Rooijen N, and McDonald V. 2012. *J. of Infect. Disease.* 206: 117-124.
2. Cobb D and Smeltz RB. 2012. *J. Immunol.* 188: 3766-3773.
3. Brown DM, Lee S, Garcia-Hernandez M, and Swain SL. 2012. *J. Virol.* 86: 6792-6803.
4. Yu H, Karunakaran KP, Jiang X, Shen C, Andersen P, and Brunham RC. 2012. *Infect. Immun.* 80: 1510-1518.
5. Kwon M-J, Ma J, Ding Y, Wang R, and Sun Z. 2012. *J. Immunol.* 188: 5887-5897.
6. Barr TA, Shen P, Brown S, Lampropoulou V, Roch T, Lawrie S, Fan B, O'Connor RA, Anderton SM, Bar-Or Am Fillatreau S, and Gray D. 2012. *J. Exp. Med.* 209: 1001-1010.
7. Cardona AE, Restrepo BI, Jaramillo JM, and Teale JM. 1999. *J. Immunol.* 162: 995-1002.
8. Kupfer A, Mosmann TR, and Kupfer H. 1991. *Proc. Natl. Acad. Sci.* 88: 775-779.

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. Cy is a registered trademark of GE Healthcare. 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