

CD44 Rat Anti-Human/Mouse mAb PE-Cyanine7 Conjugate

Store at 2°C to 8°C

Catalog Number A18709

Pub. No. MAN0009074 Rev. 2.0

Catalog No.	Form	Amount	Excitation	Peak Emission
A18709	PE-Cyanine7	25 µg	488 nm	767 nm

Product description

The CD44 Rat Anti-Human/Mouse Monoclonal Antibody (mAb) recognizes CD44, a ubiquitously expressed cell surface receptor which is important for extracellular matrix organization, cell-cell and cell-matrix adhesion and migration. CD44 may be expressed in a number of different isoforms (splice variants) from the most typical or "standard" form, known as CD44s, to variants designated CD44v, e.g. CD44v1 or CD44v6. These receptors interact with several ligands, but most often associate with an extracellular matrix component hyaluronate, through which it mediates adhesion. The IM7 clone may be used for detection of all isoforms of CD44, as it recognizes constant epitopes near the extracellular proximal domain. It has been reported to be cross-reactive with many non-human species including Baboon, Chimpanzee, Cynomolgus, Rhesus, Horse, Cow, Pig, Dog and Cat CD44.

Product specifications

Clonality:	Monoclonal
Host/Class:	Rat IgG
Reactivity:	Human/Mouse CD44
Clone/PAD:	IM7
Isotype:	IgG2bk
Lot:	See product label

Product applications

Applications reported for the CD44 Rat Anti-Human/Mouse mAb include flow cytometry^{2,5}, functional assays^{1,8,9}, immunoprecipitation³, immunohistochemistry (frozen tissue)^{4,6}, immunohistochemistry (paraffin embedded tissue)⁷, and western blot⁹.

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. Chandler HL, Haeussler DJ, Gemensky-Metzler AJ, Wilkie DA, and Lutz EA. 2012. *Invest. Ophthalmol. Vis. Sci.* 53:1835-1845.
2. Lee L-F, Logronio K, Tu GH, Zhai W, Ni I, Mei L, Dilley J, Yu J, et al. 2012. *Proc. Natl. Acad. Sci.* 10.1073.
3. Ruffell B, Poon GFT, Lee SSM, Brown KL, Tjew S-L, Cooper J, and Johnson P. 2011. *J. Biol. Chem.* 286:19179-19190.
4. Miyake Y, Matsumoto H, Yokoo M, Miyazawa K et al. 2006. *Biol. Reprod.* 74: 501-510.
5. Veir JK, Lappin MR, and Dow SW. 2006. *Journal of Feline Medicine and Surgery.* 8:400-411.
6. Frank NY, Margaryan A, Huang Y, Schatton T, Waaga-Gasser AM, Gasser M, Sayegh MH, Sadee W, and Frank MH. 2005. *Cancer Res.* 65:4320-4333.
7. Fischer A, Schumacher N, Maier M, Sendtner M, and Gessler M. 2004. *Genes and Dev.* 18:901-911.
8. Xu H, Manivannan A, Liversidge J, Sharp PF, Forrester JV, and Crane IJ. 2002. *J. Leukoc. Biol.* 72:1133-1141.
9. Si-Tahar M, Sitaraman S, Shibahara T, and Madara JL. 2001. *Am. J. Physiol. Cell Physiol.* 280:C423-C432.

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