Human IgM Immunoglobulin Recombinant Monoclonal Rabbit Antibody (IGHM/3776R)



Product Description

Recognizes a protein of 75 kDa, identified as mu heavy chain of human immunoglobulins. It does not cross-react with alpha (IgA), gamma (IgG), epsilon (IgE), or delta (IgD), heavy chains, T-cells, monocytes, granulocytes, or erythrocytes. This MAb is useful in the identification of leukemias, plasmacytomas, and certain non-Hodgkin's lymphomas. The most common feature of these malignancies is the restricted expression of a single heavy chain class. Demonstration of clonality in lymphoid infiltrates indicates that the infiltrate is clonal and therefore malignant.

Primary antibodies are available purified, or with a selection of fluorescent CF® dyes and other labels. CF® dyes offer exceptional brightness and photostability. See the CF® Dye Brochure for more information. Note: Conjugates of blue fluorescent dyes like CF®405S and CF®405M are not recommended for detecting low abundance targets, because blue dyes have lower fluorescence and can give higher non-specific background than other dye colors.

Stock status: Because Biotium offers a large number of antibody and conjugation options, primary antibody conjugates may be made to order. Typical lead times are up to one week for CF® dye and biotin conjugates, and up to 2-3 weeks for fluorescent protein and enzyme conjugates. Please email order@biotium.com to inquire about stock status and lead times before placing your order.

Catalog number key for antibody number 3776, Anti-IgM Mu Heavy Chain (IGHM/3776R)

Product attributes

Product attributes			
Antibody number	#3776		
Antibody reactivity (target)	IgM Mu Heavy Chain		
Antibody type	Primary, Anti-Human Immunoglobulin		
Host species	Rabbit		
Clonality	Recombinant Monoclonal		
Clone	IGHM/3776R		
Isotype	IgG		
Molecular weight	50-75 kDa		
Synonyms	AGM1; IGHM; Constant Region of Heavy Chain of IgM; Ig Mu Chain C Region		
Human gene symbol	IGHM		
Entrez gene ID	3507		
SwissProt	P01871 & P20769		
Unigene	510635		
Immunogen	Recombinant full-length human IGHM protein		
Verified antibody applications	IHC (FFPE) (verified)		
Antibody target cellular	Plasma membrane, Secreted (extracellular)		
localization	(
localization Species reactivity	Human		
Species reactivity	Human		
Species reactivity Positive control	Human 293T, Raji or hPBL cells. Human tonsil or spleen (IHC).		
Species reactivity Positive control Shipping condition	Human 293T, Raji or hPBL cells. Human tonsil or spleen (IHC). Room temperature Store at 2 to 8 °C, Protect fluorescent conjugates from light,		
Species reactivity Positive control Shipping condition Storage Conditions	Human 293T, Raji or hPBL cells. Human tonsil or spleen (IHC). Room temperature Store at 2 to 8 °C, Protect fluorescent conjugates from light, Note: store BSA-free antibodies at -10 to -35 °C		
Species reactivity Positive control Shipping condition Storage Conditions Regulatory status Antibody/conjugate	Human 293T, Raji or hPBL cells. Human tonsil or spleen (IHC). Room temperature Store at 2 to 8 °C, Protect fluorescent conjugates from light, Note: store BSA-free antibodies at -10 to -35 °C For research use only (RUO) Conjugates: 0.1 mg/mL in PBS/0.1% BSA/0.05% azide, HRP conjugates: 0.1 mg/mL in PBS/0.05% BSA, Purified: 0.2 mg/mL in PBS/0.05% BSA/0.05% azide, Purified: 8SA-free: 1 mg/mL in		
Species reactivity Positive control Shipping condition Storage Conditions Regulatory status Antibody/conjugate formulation	Human 293T, Raji or hPBL cells. Human tonsil or spleen (IHC). Room temperature Store at 2 to 8 °C, Protect fluorescent conjugates from light, Note: store BSA-free antibodies at -10 to -35 °C For research use only (RUO) Conjugates: 0.1 mg/mL in PBS/0.1% BSA/0.05% azide, HRP conjugates: 0.1 mg/mL in PBS/0.05% BSA, Purified: 0.2 mg/mL in PBS/0.05% BSA/0.05% azide, Purified, BSA-free: 1 mg/mL in PBS without azide Guaranteed for at least 24 months from date of receipt when		
Species reactivity Positive control Shipping condition Storage Conditions Regulatory status Antibody/conjugate formulation Shelf life	Human 293T, Raji or hPBL cells. Human tonsil or spleen (IHC). Room temperature Store at 2 to 8 °C, Protect fluorescent conjugates from light, Note: store BSA-free antibodies at -10 to -35 °C For research use only (RUO) Conjugates: 0.1 mg/mL in PBS/0.1% BSA/0.05% azide, HRP conjugates: 0.1 mg/mL in PBS/0.05% BSA, Purified: 0.2 mg/mL in PBS/0.05% BSA/0.05% azide, Purified; 0.2 mg/mL in PBS without azide Guaranteed for at least 24 months from date of receipt when stored as recommended Product may contain either bovine serum albumin (BSA) from bovine serum (Bos taurus), or recombinant BSA produced in		
Species reactivity Positive control Shipping condition Storage Conditions Regulatory status Antibody/conjugate formulation Shelf life Product origin	Human 293T, Raji or hPBL cells. Human tonsil or spleen (IHC). Room temperature Store at 2 to 8 °C, Protect fluorescent conjugates from light, Note: store BSA-free antibodies at -10 to -35 °C For research use only (RUO) Conjugates: 0.1 mg/mL in PBS/0.1% BSA/0.05% azide, HRP conjugates: 0.1 mg/mL in PBS/0.05% BSA, Purified: 0.2 mg/mL in PBS/0.05% BSA/0.05% azide, Purified: 0.2 mg/mL in PBS without azide Guaranteed for at least 24 months from date of receipt when stored as recommended Product may contain either bovine serum albumin (BSA) from bovine serum (Bos taurus), or recombinant BSA produced in Chinese hamster ovary cells. Inquire for the specific lot.		

Call us: 800-304-5357 Email: btinfo@biotium.com

Antibody # prefix	Conjugation	Ex/Em (nm)	Laser line	Detection channel	Dye Features
BNC04	CF®405S	404/431	405	DAPI (microscopy), AF405	CF®405S Features
BNC88	CF®488A	490/515	488	GFP, FITC	CF®488A Features
BNC68	CF®568	562/583	532, 561	RFP, TRITC	CF®568 Features
BNC94	CF®594	593/614	561	Texas Red®	CF®594 Features
BNC40	CF®640R	642/662	633-640	Cy®5	CF®640R Features
BNC47	CF®647	650/665	633-640	Cy®5	CF®647 Features
BNC74	CF®740	742/767	633-685	775/50	CF®740 Features
BNCB	Biotin	N/A	N/A	N/A	
BNUB	Purified	N/A	N/A	N/A	
BNUM	Purified, BSA-free	N/A	N/A	N/A	

Alexa Fluor, Pacific Blue, Pacific Orange, and Texas Red are trademarks or registered trademarks of Thermo Fisher Scientific; Cy is a registered trademark of Cytiva; IRDye, LI-COR, and Odyssey are registered trademarks of LI-COR Bioscience.

This datasheet was generated on November 4, 2025 at 02:59:29 PM. Visit product page to check for updated information before use. Product link: https://biotium.com/product/human-igm-immunoglobulin-recombinant-monoclonal-rabbit-antibody-ighm-3776r/