CD21 Monoclonal Mouse Antibody (CR2/1953)



Product Description

This antibody recognizes a protein of 140 kDa, which is identified as the complement receptor 2 (CR2) / CD21. This protein is expressed strongly on mature B cells, follicular dendritic cells and weakly on immature thymocytes and T lymphocytes. In B-cell ontogeny, CD21 appears after the pre-B-stage, is maintained during peripheral B-cell development and is lost upon terminal differentiation into plasma cells. CD21 expression is also gradually lost after stimulation of B cells in vitro. CD21 functions as receptor for C3d, C3dg and iC3b Complement components, for EBV and for IFNalpha. CD21 binds to CD23 and associates with CD19, CD81 and Leu13 to form a large signal-transduction complex involved in B cell activation.

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 1953, Anti-CD21 (CR2/1953)

Product attributes

Call us: 800-304-5357

Product attributes			
Antibody number	#1953		
Antibody reactivity (target)	CD21		
Antibody type	Primary		
Host species	Mouse		
Clonality	Monoclonal		
Clone	CR2/1953		
Isotype	IgG1, kappa		
Molecular weight	140 kDa		
Synonyms	CD21; Complement C3d receptor (C3DR); Complement Receptor type 2 (CR2); CVID7; EBV receptor; EBV-R; Epstein-Barr virus receptor; EVBR; SLEB9		
Human gene symbol	CR2		
Entrez gene ID	1380		
SwissProt	P20023		
Unigene	445757		
Immunogen	Recombinant human CR2 (CD21) protein fragment (around aa 44-196) (exact sequence is proprietary)		
Verified antibody applications	IHC (FFPE) (verified)		
Antibody target cellular localization	Plasma membrane		
Species reactivity	Human		
Positive control	WEHI-231 cells. Human tonsil tissue.		
Shipping condition	Room temperature		
Storage Conditions	Store at 2 to 8 $^{\circ}$ C, Protect fluorescent conjugates from light, Note: store BSA-free antibodies at -10 to -35 $^{\circ}$ C		
Shelf life	Guaranteed for at least 24 months from date of receipt when stored as recommended		
Regulatory status	For research use only (RUO)		
Antibody/conjugate formulation	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		
Validated in protein array	Monospecific		
Antibody research areas	Immunology		
Product origin	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.		
Cell/tissue expression	Antigen presenting cells, B-cells, Lymphocytes		

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,	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 December 12, 2025 at 03:18:53 AM. Visit product page to check for updated information before use. Product link: https://biotium.com/product/cd21-monoclonal-mouse-antibody-cr2-1953/