## C1QB / Complement C1q B-Chain Monoclonal Mouse Antibody (C1QB/2961)



## **Product Description**

Complement C1q B Chain (C1qb) is a subunit of C1q. C1q is composed of 6 polypeptide chains each of C1qa, C1qb, and C1qc. C1q is a component of the classical complement pathway, and thereby plays an important role in the immune response. Each C1q subunit contains an N-terminal collagen-like region and a C-terminal C1q globular domain. The presence of receptors for C1q on effector cells modulates its activity, which may be antibody-dependent or independent. Macrophages are the primary source of C1q, while anti-inflammatory drugs as well as cytokines differentially regulate expression of the mRNA as well as the protein. C1q deficiency is associated with lupus erythematosus and glomerulonephritis.

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 <a href="mailto:order@biotium.com">order@biotium.com</a> to inquire about stock status and lead times before placing your order.

Catalog number key for antibody number 2961, C1QB / Complement C1q B-Chain Monoclonal Mouse Antibody (C1QB/2961)

## Product attributes

Fiduuci attiibutes			
Antibody number	•		
Antibody reactivity (target)	C1QB, Complement C1q B-Chain		
Antibody type	Primary		
Host species	Mouse		
Clonality	Monoclonal		
Clone	C1QB/2961		
Isotype	IgG1, kappa		
Molecular weight	26-29 kDa		
Synonyms	C1QB; Complement component 1 q subcomponent B chain; Complement component C1q B chain		
Human gene symbol	C1QB		
Entrez gene ID	713		
SwissProt	P02746		
Unigene	8986		
Immunogen	Recombinant fragment (around aa 41-188) of human C1QB protein (exact sequence is proprietary)		
Verified antibody applications	IHC (FFPE) (verified)		
Antibody target cellular localization	Secreted (extracellular)		
Species reactivity	Human		
Positive control	Human liver, kidney or brain tissue (IHC).		
Shipping condition	,,		
Storage Conditions	Store at 2 to 8 °C, Protect fluorescent conjugates from light, Note: store BSA-free antibodies at -10 to -35 °C		
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% BSA/0.05% azide, Purified, BSA-free: 1 mg/mL in PBS without azide		
	F B3 WILLIOUT AZIDE		
Validated in protein array	Monospecific Monospecific		
Validated in protein array Shelf life			
•	Monospecific  Guaranteed for at least 24 months from date of receipt when		
Shelf life	Monospecific  Guaranteed for at least 24 months from date of receipt when stored as recommended		

Call us: 800-304-5357 Email: techsupport@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 September 9, 2025 at 07:38:01 PM. Visit product page to check for updated information before use. Product link: <a href="https://biotium.com/product/c1qb-complement-c1q-b-chain-monoclonal-mouse-antibody-c1qb-2961/">https://biotium.com/product/c1qb-complement-c1q-b-chain-monoclonal-mouse-antibody-c1qb-2961/</a>