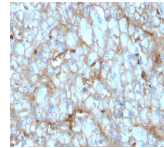


C1QA / Complement C1q A-Chain Monoclonal Mouse Antibody (C1QA/2954)



Product Description

Complement C1q A Chain (C1qa) 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 order@biotium.com to inquire about stock status and lead times before placing your order. **Catalog number key for antibody number 2954, C1QA / Complement C1q A-Chain Monoclonal Mouse Antibody (C1QA/2954)**

Product attributes

Antibody number	#2954
Antibody reactivity (target)	C1QA, Complement C1q A-Chain
Antibody type	Primary
Host species	Mouse
Clonality	Monoclonal
Clone	C1QA/2954
Isotype	IgG2b, kappa
Molecular weight	29 kDa
Synonyms	C1QA; Complement C1q subcomponent subunit A; Complement component 1 q subcomponent A chain; Complement component 1 q subcomponent alpha polypeptide; Complement component C1q A chain
Human gene symbol	C1QA
Entrez gene ID	712
SwissProt	P02745
Unigene	632379
Immunogen	Recombinant fragment (around aa 104-237) of human C1QA protein (exact sequence is proprietary)
Verified antibody applications	IHC (FFPE) (verified), WB (verified)
Antibody target cellular localization	Secreted (extracellular)
Species reactivity	Human
Positive control	Liver or Kidney.
Shipping condition	Room temperature
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% azide, Purified, BSA-free: 1 mg/mL in PBS without azide
Validated in protein array	Monospecific
Shelf life	Guaranteed for at least 24 months from date of receipt when stored as recommended
Antibody research areas	Immunology
Cell/tissue expression	Macrophage/monocyte

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
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, LI-COR Bioscience.

This datasheet was generated on June 29, 2024 at 04:37:58 PM. Visit product page to check for updated information before use.

Product link: <https://biotium.com/product/c1qa-complement-c1q-a-chain-monoclonal-mouse-antibody-c1qa-2954/>