## CD68 Monoclonal Mouse Antibody (C68/2501)



## **Product Description**

This antibody recognizes a glycoprotein of 110 kDa, which is identified as CD68. It is important for identifying macrophages in tissue sections. It stains macrophages in a wide variety of human tissues, including Kupffer cells and macrophages in the red pulp of the spleen, in lamina propria of the gut, in lung alveoli, and in bone marrow. It reacts with myeloid precursors and peripheral blood granulocytes. It also reacts with plasmacytoid T cells, which are supposed to be of monocyte/macrophage origin. It shows strong granular cytoplasmic staining of chronic and acute myeloid leukemia and also reacts with rare cases of true histiocytic neoplasia. Lymphomas are negative or show few granules.

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 2501, Anti-CD68 (C68/2501)

Product attributes				
Antibody number	#2501			
Antibody reactivity (target)	CD68			
Antibody type	Primary			
Host species	Mouse			
Clonality	Monoclonal			
Clone	C68/2501			
Isotype	IgG2c, kappa			
Molecular weight	~110 kDa			
Synonyms	GP110, LAMP4, Microsialin, Macrosialin, SCARD1, Scavenger Receptor Class D Member-1			
Human gene symbol	CD68			
Entrez gene ID	968			
SwissProt	P34810			
Unigene	647419			
Immunogen	Recombinant fragment of human CD68 protein (around aa 150-221) (exact sequence is proprietary)			
Verified antibody applications	IHC (FFPE) (verified)			
Antibody target cellular localization	Plasma membrane			
Species reactivity	Human			
Antibody application notes	Higher concentration may be required for direct detection using primary antibody conjugates than for indirect detection with secondary antibody. Flow cytometry: 0.5-1 ug/million cells, Immunofluorescence: 0.5-1 ug/mL, Immunofluorescence: 0.5-1 ug/mL, Immunofluorescence: 0.5-1 ug/mL, Immunofluorescence: 0.5-1 ug/mL, Istaining of formalin-fixed tissues is enhanced by boiling tissue sections in 10 mM citrate buffer pH 6.0 for 10-20 minutes followed by cooling at RT for 20 minutes, Western: 0.5-1 ug/mL, Optimal dilution for a specific application should be determined by user			
	buffer pH 6.0 for 10-20 minutes followed by cooling at RT for 20 minutes, Western: 0.5-1 ug/mL, Optimal dilution for a specific			
Positive control	buffer pH 6.0 for 10-20 minutes followed by cooling at RT for 20 minutes, Western: 0.5-1 ug/mL, Optimal dilution for a specific			
Positive control Shipping condition	buffer pH 6.0 for 10-20 minutes followed by cooling at RT for 20 minutes, Western: 0.5-1 ug/mL, Optimal dilution for a specific application should be determined by user			
	buffer pH 6.0 for 10-20 minutes followed by cooling at RT for 20 minutes, Western: 0.5-1 ug/mL, Optimal dilution for a specific application should be determined by user  Tonsil, lymph node, or spleen			
Shipping condition	buffer pH 6.0 for 10-20 minutes followed by cooling at RT for 20 minutes, Western: 0.5-1 ug/mL, Optimal dilution for a specific application should be determined by user  Tonsil, lymph node, or spleen  Room temperature  Store at 2 to 8 °C, Protect fluorescent conjugates from light,			
Shipping condition Storage Conditions	buffer pH 6.0 for 10-20 minutes followed by cooling at RT for 20 minutes, Western: 0.5-1 ug/ml., Optimal dilution for a specific application should be determined by user  Tonsil, lymph node, or spleen  Room temperature  Store at 2 to 8 °C, Protect fluorescent conjugates from light, Note: store BSA-free antibodies at -10 to -35 °C  Guaranteed for at least 24 months from date of receipt when			
Shipping condition Storage Conditions Shelf life	buffer pH 6.0 for 10-20 minutes followed by cooling at RT for 20 minutes, Western: 0.5-1 ug/ml., Optimal dilution for a specific application should be determined by user  Tonsil, lymph node, or spleen  Room temperature  Store at 2 to 8 °C, Protect fluorescent conjugates from light, Note: store BSA-free antibodies at -10 to -35 °C  Guaranteed for at least 24 months from date of receipt when stored as recommended			
Shipping condition Storage Conditions Shelf life Regulatory status Antibody/conjugate	buffer pH 6.0 for 10-20 minutes followed by cooling at RT for 20 minutes, Western 0.5-1 ug/ml., Optimal dilution for a specific application should be determined by user  Tonsil, lymph node, or spleen  Room temperature  Store at 2 to 8 °C, Protect fluorescent conjugates from light, Note: store BSA-free antibodies at -10 to -35 °C  Guaranteed for at least 24 months from date of receipt when stored as recommended  For research use only (RUO)  Conjugates: 0.1 mg/mL in PBS/0.05% BSA, Purified: 0.2 mg/mL in PBS/0.05% BSA, Purified; 0.2 mg/mL in PBS/0.05% BSA, Puri			
Shipping condition Storage Conditions Shelf life Regulatory status Antibody/conjugate formulation	buffer pH 6.0 for 10-20 minutes followed by cooling at RT for 20 minutes, Western: 0.5-1 ug/ml., Optimal dilution for a specific application should be determined by user  Tonsil, lymph node, or spleen  Room temperature  Store at 2 to 8 °C, Protect fluorescent conjugates from light, Note: store BSA-free antibodies at -10 to -35 °C  Guaranteed for at least 24 months from date of receipt when stored as recommended  For research use only (RUO)  Conjugates: 0.1 mg/mL in PBS/0.05% BSA, Purified: 0.2 mg/mL in PBS/0.05% BSA, Purified; 0.2 mg/mL in PBS without azide			
Shipping condition Storage Conditions Shelf life Regulatory status Antibody/conjugate formulation Validated in protein array	buffer pH 6.0 for 10-20 minutes followed by cooling at RT for 20 minutes, Western: 0.5-1 ug/ml., Optimal dilution for a specific application should be determined by user  Tonsil, lymph node, or spleen  Room temperature  Store at 2 to 8 °C, Protect fluorescent conjugates from light, Note: store BSA-free antibodies at -10 to -35 °C  Guaranteed for at least 24 months from date of receipt when stored as recommended  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% BSA/0.05% azide, Purified, BSA-free: 1 mg/mL in PBS without azide  Monospecific			
Shipping condition Storage Conditions Shelf life Regulatory status Antibody/conjugate formulation  Validated in protein array Antibody research areas	buffer pH 6.0 for 10-20 minutes followed by cooling at RT for 20 minutes, Western: 0.5-1 ug/mL, Optimal dilution for a specific application should be determined by user  Tonsil, lymph node, or spleen  Room temperature  Store at 2 to 8 °C, Protect fluorescent conjugates from light, Note: store BSA-free antibodies at -10 to -35 °C  Guaranteed for at least 24 months from date of receipt when stored as recommended  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  Monospecific  Immunology  Product may contain either bovine serum albumin (BSA) from bovine serum (Bos taurus), or recombinant BSA produced in			

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 2, 2025 at 05:53:37 PM. Visit product page to check for updated information before use. Product link: <a href="https://biotium.com/product/cd68-monoclonal-mouse-antibody-c68-2501/">https://biotium.com/product/cd68-monoclonal-mouse-antibody-c68-2501/</a>