## CD14 Monoclonal Mouse Antibody (LPSR/2408)



## **Product Description**

This antibody recognizes a protein of 55 kDa, identified as CD14 (also known lipopolysaccharide receptor). CD14 is expressed strongly on monocytes and macrophage and weakly on the surface of neutrophils. CD14 is anchored to cells by linkage to glycosylphosphatidylinositol (GPI) and functions as a high affinity receptor for complexes of LPS and LPS binding protein (LBP). Soluble CD14, also binding to LPS, acts at physiological concentration as an LPS agonist and has, at higher concentrations, an LPS antagonizing effect in 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 <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 2408, Anti-CD14 (LPSR/2408)

## Product attributes

Call us: 800-304-5357

Product attributes				
Antibody number	#2408			
Antibody reactivity (target)	CD14			
Antibody type	Primary			
Host species	Mouse			
Clonality	Monoclonal			
Clone	LPSR/2408			
Isotype	lgG1, kappa			
Molecular weight	ight 55 kDa			
Synonyms	Lipopolysaccharide receptor; LPS-R; LPSR; Mo2; Monocyte Differentiation Antigen 14; Myeloid cell-specific leucine-rich glycoprotein			
Human gene symbol	CD14			
Entrez gene ID	929			
SwissProt	P08571			
Unigene	163867			
Immunogen	Recombinant fragment of human CD14 protein (around aa 25-148) (exact sequence is proprietary)			
Verified antibody applications	IHC (FFPE) (verified), WB (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, Immunohistology (formalin): 1-2 ug/mL for 30 minutes at RT, Staining of formalin-fixed tissues is enhanced by boiling tissue sections in 10 mM Tris with 1 mM EDTA pH 9.0			
	for 10-20 minutes followed by cooling at RT for 20 minutes, Optimal dilution for a specific application should be determined by user			
Positive control	Optimal dilution for a specific application should be determined			
Positive control Shipping condition	Optimal dilution for a specific application should be determined by user			
	Optimal dilution for a specific application should be determined by user  THP-1 cells and human monocytes. Lymph nodes and tonsils.			
Shipping condition	Optimal dilution for a specific application should be determined by user  THP-1 cells and human monocytes. Lymph nodes and tonsils.  Room temperature  Store at 2 to 8 °C, Protect fluorescent conjugates from light,			
Shipping condition Storage Conditions	Optimal dilution for a specific application should be determined by user  THP-1 cells and human monocytes. Lymph nodes and tonsils.  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	Optimal dilution for a specific application should be determined by user  THP-1 cells and human monocytes. Lymph nodes and tonsils.  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	Optimal dilution for a specific application should be determined by user  THP-1 cells and human monocytes. Lymph nodes and tonsils.  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/0.05% azide, Purified; BSA-free: 1 mg/mL in			
Shipping condition Storage Conditions Shelf life Regulatory status Antibody/conjugate formulation	Optimal dilution for a specific application should be determined by user  THP-1 cells and human monocytes. Lymph nodes and tonsils.  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			
Shipping condition Storage Conditions Shelf life Regulatory status Antibody/conjugate formulation Validated in protein array	Optimal dilution for a specific application should be determined by user  THP-1 cells and human monocytes. Lymph nodes and tonsils.  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  Monospecific			

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,	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 3, 2025 at 11:20:52 PM. Visit product page to check for updated information before use. Product link: <a href="https://biotium.com/product/cd14-monoclonal-mouse-antibody-lpsr-2408/">https://biotium.com/product/cd14-monoclonal-mouse-antibody-lpsr-2408/</a>