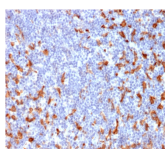


CD163 Monoclonal Mouse Antibody (M130/2162)



Product Description

This MAb recognizes a protein of 140 kDa, identified as CD163. It has been identified as an acute phase-regulated transmembrane protein whose function is to mediate the endocytosis of haptoglobin-hemoglobin complexes. This receptor is expressed on the surface of monocytes with low expression and on tissue macrophages, histiocytes with high expression. Staining with anti-CD163 has been helpful to distinguish synovial macrophages from synovial intimal fibroblasts in rheumatoid arthritis, where its specificity for macrophages was found to be superior to that of anti-CD68. Increased levels of CD163 were also detected in patients with microbial infections and myelomonocytic leukemias.

Anti-CD163 is of considerable value for selective identification of monocytes and macrophages at a certain stage of differentiation and is suitable for diagnosing myelomonocytic or monocytic leukaemia and neoplasms of true histiocytic origin. CD163 is positive in skin (histiocytes), gut, Kupffer cells, a few alveolar macrophages, macrophages in the placenta, and in macrophages in inflamed tissues including tumor tissue.

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 2162, Anti-CD163 (M130/2162)

Product attributes

Antibody number	#2162
Antibody reactivity (target)	CD163
Antibody type	Primary
Host species	Mouse
Clonality	Monoclonal
Clone	M130/2162
Isotype	IgG2b, kappa
Molecular weight	140 kDa
Synonyms	CD163; CD163 antigen; Macrophage-associated antigen; M130; CD163 molecule; Hemoglobin scavenger receptor; MM130; Scavenger receptor cysteine rich type 1 protein M130
Human gene symbol	CD163
Entrez gene ID	9332
SwissProt	Q86V7
Unigene	504641
Immunogen	Recombinant fragment (around aa 43-196) of human CD163 (exact sequence is proprietary)
Verified antibody applications	IHC (FFPE) (verified)
Antibody target cellular localization	Plasma membrane
Species reactivity	Human
Positive control	hPBL Cells. Skin, Lung, Placenta or Histiocytoma.
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
Product origin	Product may contain either bovine serum albumin (BSA) from bovine serum (<i>Bos taurus</i>), or recombinant BSA produced in Chinese hamster ovary cells. Inquire for the specific lot.
Cell/tissue expression	Monocytes/macrophages

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 27, 2024 at 10:29:04 AM. Visit [product page](#) to check for updated information before use.

Product link: <https://biotium.com/product/cd163-monoclonal-mouse-antibody-m130-2162/>