## Myeloid Associated Differentiation Marker Monoclonal Mouse Antibody (MYADM/971)



## **Product Description**

Recognizes a myeloid associated differentiation antigen in the cytoplasm of mature granulocytes. It shows no reactivity with any other cell type in human tissues. Markers of myeloid cells are useful in the identification of different levels of cellular differentiation. It reacts with early precursor and mature forms of human and monkey myeloid cells. This MAb is useful for the detection of myeloid leukemias and granulocytic sarcomas. It can be used as a marker of granulocytes in normal tissues or inflammatory processes.

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 0971, Anti-Myeloid Associated Differentiation Marker (MYADM/971)

## Product attributes

Product attributes				
Antibody number	#0971			
Antibody reactivity (target)	Myeloid Associated Differentiation Marker			
Antibody type	Primary			
Host species	Mouse			
Clonality	Monoclonal			
Clone	MYADM/971			
Isotype	lgG1			
Molecular weight	Unknown			
Synonyms	MYADM; myeloid associated differentiation marker; Myeloid upregulated protein; Protein SB135			
Entrez gene ID	91663			
SwissProt	Q96S97			
Unigene	380906			
Immunogen	Recombinant human MYADM protein			
Verified antibody applications	IHC (FFPE) (verified)			
Antibody target cellular localization	Cytoplasmic			
Species reactivity	Human, Macaque			
Antibody application notes	Higher concentration may be required for direct detection using primary antibody conjugates than for indirect detection with secondary antibody, Immunofluorescence: 1-2 ug/mL, Immunohistology formalin-fixed 0.5-1 ug/mL, Staining of formalin-fixed tissues requires boiling tissue sections in 10 mM citrate buffer, pH 6.0, for 10-20 min followed by cooling at RT for 20 minutes, Flow Cytometry 0.5-1 ug/million cells/0.1 mL, Optimal dilution for a specific application should be determined by user			
	secondary antibody, İmmunofluorescence: 1-2 ug/mL, Immunohistology formalin-fixed 0.5-1 ug/mL, Staining of formalin-fixed tissues requires boiling tissue sections in 10 mM citrate buffer, pH 6.0, for 10-20 min followed by cooling at RT for 20 minutes, Flow Cytometry 0.5-1 ug/million cells/0.1 mL, Optimal dilution for a specific application should be determined			
Positive control	secondary antibody, İmmunofluorescence: 1-2 ug/mL, Immunohistology formalin-fixed 0.5-1 ug/mL, Staining of formalin-fixed tissues requires boiling tissue sections in 10 mM citrate buffer, pH 6.0, for 10-20 min followed by cooling at RT for 20 minutes, Flow Cytometry 0.5-1 ug/million cells/0.1 mL, Optimal dilution for a specific application should be determined			
Positive control Shipping condition	secondary antibody, İmmunofluorescence: 1-2 ug/mL, Immunohistology formalin-fixed 0.5-1 ug/mL, Staining of formalin-fixed tissues requires boiling tissue sections in 10 mM citrate buffer, pH 6.0, for 10-20 min followed by cooling at RT for 20 minutes, Flow Cytometry 0.5-1 ug/million cells/0.1 mL, Optimal dilution for a specific application should be determined by user			
	secondary antibody, İmmunofluorescence: 1-2 ug/ml., Immunohistology formalin-fixed 0.5-1 ug/ml., Staining of formalin-fixed tissues requires boiling tissue sections in 10 mM citrate buffer, pH 6.0, for 10-20 min followed by cooling at RT for 20 minutes, Flow Cytometry 0.5-1 ug/million cells/0.1 mL, Optimal dilution for a specific application should be determined by user  HL60 cells. Tonsil or lymph node			
Shipping condition	secondary antibody, İmmunofluorescence: 1-2 ug/mL, Immunohistology formalin-fixed 0.5-1 ug/mL, Staining of formalin-fixed tissues sequires boiling tissue sections in 10 mM citrate buffer, pH 6.0, for 10-20 min followed by cooling at RT for 20 minutes, Flow Cytometry 0.5-1 ug/million cells/0.1 mL, Optimal dilution for a specific application should be determined by user  HL60 cells. Tonsil or lymph node  Room temperature  Store at 2 to 8 °C, Protect fluorescent conjugates from light,			
Shipping condition Storage Conditions	secondary antibody, İmmunofluorescence: 1-2 ug/mL, Immunohistology formalin-fixed 0.5-1 ug/mL, Staining of formalin-fixed tissues requires boiling tissue sections in 10 mM citrate buffer, pH 6.0, for 10-20 min followed by cooling at RT for 20 minutes, Flow Cytometry 0.5-1 ug/million cells/0.1 mL, Optimal dilution for a specific application should be determined by user  HL60 cells. Tonsil or lymph node  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	secondary antibody, İmmunofluorescence: 1-2 ug/mL, Immunohistology formalin-fixed 0.5-1 ug/mL, Staining of formalin-fixed tissues requires boiling tissue sections in 10 mM citrate buffer, pH 6.0, for 10-20 min followed by cooling at RT for 20 minutes, Flow Cytometry 0.5-1 ug/million cells/0.1 mL, Optimal dilution for a specific application should be determined by user  HL60 cells. Tonsil or lymph node  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	secondary antibody, İmmunofluorescence: 1-2 ug/ml., Immunohistology formalin-fixed 0.5-1 ug/ml., Staining of formalin-fixed tissues requires boiling tissue sections in 10 mM citrate buffer, pH 6.0, for 10-20 min followed by cooling at RT for 20 minutes, Flow Cytometry 0.5-1 ug/million cells/0.1 mL. Optimal dilution for a specific application should be determined by user  HL60 cells. Tonsil or lymph node  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, Purified; 0.2 mg/mL in PBS/0.05% BSA, Purified; BSA-free: 1 mg/mL in PBS/0.05% BSA, Purified; mg/mL in PBS/0.05% BSA, Purified; mg/mL in PBS/0.05% BSA, Purified; mg/mL in PBS/0.05% BSA, Purified; mg/mL in PBS/0.05% BSA, Purified; mg/mL in PBS/0.05% BSA, Purified; mg/mL in PBS/0.05% BSA, Purified; mg/mL in PBS/0.05% BSA, Purified; mg/mL in PBS/0.05% BSA, Purified; mg/mL in PBS/0.05% BSA, Purified; mg/mL in PBS/0.05% BSA/0.05% acide, Purified; mg/mL in PBS/0.05% BSA/0.05% acide, Purified; mg/mL in PBS/0.05% BSA/0.05% acide, Purified; mg/mL in PBS/0.05% BSA/0.05% acide, Purified; mg/mL in PBS/0.05% BSA/0.05% acide, Purified; mg/mL in PBS/0.05% BSA/0.05% acide, Purified; mg/mL in PBS/0.05% BSA/0.05% acide, Purified; mg/mL in PBS/0.05% BSA/0.05% acide, Purified; mg/mL in PBS/0.05% BSA/0.05% acide, Purified; mg/mL in PBS/0.05% BSA/0.05% acide, Purified; mg/mL in PBS/0.05% BSA/0.05% acide, Purified; mg/mL in PBS/0.05% BSA/0.05% acide, Purified; mg/mL in PBS/0.05% BSA/0.05% acide, Purified; mg/mL in PBS/0.05% BSA/0.05% acide, Purified; mg/mL in PBS/0.05% BSA/0.05% acide, Purified; mg/mL in PBS/0.05% BSA/0.05% acide, Purified; mg/mL in PBS/0.05% BSA/0.05% acide, Purified; mg/mL in PBS/0.05% BSA/0.05% acide, Purified; mg/mL in PBS/0.05% BSA/0.05% acide, Purified; mg/mL			
Shipping condition Storage Conditions Shelf life Regulatory status Antibody/conjugate formulation	secondary antibody, İmmunofluorescence: 1-2 ug/ml., Immunohistology formalin-fixed 0.5-1 ug/ml., Staining of formalin-fixed tissues requires boiling tissue sections in 10 mM citrate buffer, pH 6.0, for 10-20 min followed by cooling at RT for 20 minutes, Flow Cytometry 0.5-1 ug/million cells/0.1 mL. Optimal dilution for a specific application should be determined by user  HL60 cells. Tonsil or lymph node  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/0.05% azide, HRP conjugates: 0.1 mg/mL in PBS/0.05% BSA/Preie: 1 mg/mL in PBS/0.05% BSA/Preie: 1 mg/mL in PBS without azide			

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 5, 2025 at 11:05:21 PM. Visit product page to check for updated information before use. Product link: <a href="https://biotium.com/product/monoclonal-anti-myeloid-associated-differentiation-marker-myadm971/">https://biotium.com/product/monoclonal-anti-myeloid-associated-differentiation-marker-myadm971/</a>