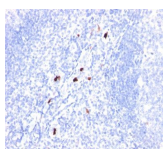


Granulocyte Marker Monoclonal Mouse Antibody (BM-1)



Product Description

Recognizes 183 kDa protein with DNA-binding characteristics, which is identified as a myeloid specific antigen. It reacts with myeloid precursor cells and granulocytes in bone marrow. Its antigen appears to be restricted to M2 and M3 acute myelogenous leukemia (AML) subtypes. Markers of myeloid cells are useful in the identification of different levels of cellular differentiation. This MAb reacts with early precursor and mature forms of human myeloid cells. It is useful in the identification of myelogenous leukemias, distinguishing granulocytic sarcomas from lymphoid malignancies and also in the study of differentiation and transformation of human myeloid cells. The biological function of this antigen is not clear, although it has been proposed that it may play a role in the differentiation of myeloid cells.

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 0061, Anti-Granulocyte Marker (BM-1)

Product attributes

Antibody number	#0061
Antibody reactivity (target)	Granulocyte Marker
Antibody type	Primary
Host species	Mouse
Clonality	Monoclonal
Clone	BM-1
Isotype	IgG1
Molecular weight of antigen	183 kDa
Synonyms	Myeloid specific protein; p183 DNA-binding cytoplasmic protein
Entrez gene ID	Not Known
SwissProt	Not Known
Unigene	Not Known
Immunogen	Human peripheral blood mononuclear cells
Antibody target cellular localization	Nucleus & cytoplasm
Verified antibody applications	IHC (FFPE) (verified)
Species reactivity	Human
Expected antibody applications	RIA (published for clone), IHC (frozen) (published for clone), IF (published for clone), WB (published for clone)
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, No special pretreatment is required for staining of formalin/paraffin tissues, Flow Cytometry 0.5-1 ug/million cells/0.1 mL, Optimal dilution for a specific application should be determined by user
Positive control	HL60 cells, Bone marrow, lymph node or tonsil
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
Shelf life	Guaranteed for at least 24 months from date of receipt when stored as recommended
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
Antibody research areas	Immunology
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	Granulocytes, Myeloid cells
Tumor expression	Leukemia/lymphoma

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.

References

Note: References for this clone sold by other suppliers may be listed for expected applications.

Blood (1987) 70(4): 1124-1130. (IHC, FFPE; IHC, frozen; IF; WB; RIA)

This datasheet was generated on June 3, 2026 at 07:21:07 PM. Visit product page to check for updated information before use.

Product link: <https://biotium.com/product/monoclonal-anti-granulocyte-marker-bm-1/>