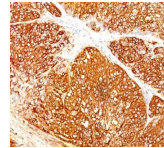


Melanoma Marker Monoclonal Mouse Antibody (HMB45 + A103 + T311)



Product Description

This antibody cocktail recognizes three melanoma-specific proteins, which include MART-1, Tyrosinase and gp100. MART-1 is a newly identified melanocyte differentiation antigen recognized by autologous cytotoxic T lymphocytes. Tyrosinase is one of the targets for cytotoxic T-cell recognition in melanoma patients. The function of gp100 is not known but it is reported to be a useful marker for melanocytes and melanomas. This cocktail of three markers is designed for extremely sensitive labeling of formalin-fixed, paraffin-embedded melanomas and other tumors showing melanocytic differentiation.

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 0702, Anti-Melanoma Marker (HMB45 A103 T311)

Product attributes

Antibody number	#0702
Antibody reactivity (target)	Melanoma Marker
Antibody type	Primary
Host species	Mouse
Clonality	Monoclonal
Clone	HMB45 + A103 + T311
Isotype	IgG's
Molecular weight	20-22 kDa (doublet) (MART); 70-80 kDa (Tyrosinase); 90-100 kDa (gp100)
Synonyms	Melanoma antigen recognized by T-cells 1 (MART-1), MLAN-A; TYR; PMEL17
Entrez gene ID	2315-MART-1; 7299-Tyrosinase; 6490-gp100
SwissProt	Q16655-MART-1; P14679-Tyrosinase; P40967-gp100
Unigene	154069-MART-1; 503555-Tyrosinase; 95972-gp100
Immunogen	Recombinant hMART-1 protein (A103); Recombinant tyrosinase protein (T311); Extract of pigmented melanoma metastases from lymph nodes (HMB45)
Antibody target cellular localization	Cytoplasmic
Verified antibody applications	IHC (FFPE) (verified), WB (verified)
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-fixed) 0.5-1.0 ug/mL, Staining of formalin-fixed tissues is enhanced by boiling tissue sections in 10 mM citrate buffer, pH 6.0, for 10-20 min followed by cooling at RT for 20 minutes, Optimal dilution for a specific application should be determined by user
Positive control	SK-MEL-13 and SK-MEL-19 Melanoma cell lines; Melanomas
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	Cancer
Cell/tissue expression	Melanocytes
Tumor expression	Melanoma

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
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, of LI-COR Bioscience.