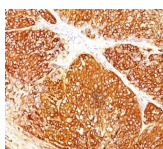


# Melanoma Marker Monoclonal Mouse Antibody (HMB45 + A103 + T311)



## 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](mailto: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

<b>Antibody number</b>	#0702
<b>Antibody reactivity (target)</b>	Melanoma Marker
<b>Antibody type</b>	Primary
<b>Host species</b>	Mouse
<b>Clonality</b>	Monoclonal
<b>Clone</b>	HMB45 + A103 + T311
<b>Isotype</b>	IgG's
<b>Molecular weight</b>	20-22 kDa (doublet) (MART); 70-80 kDa (Tyrosinase); 90-100 kDa (gp100)
<b>Synonyms</b>	Melanoma antigen recognized by T-cells 1 (MART-1), MLAN-A; TYR; PMEL17
<b>Entrez gene ID</b>	2315-MART-1; 7299-Tyrosinase; 6490-gp100
<b>SwissProt</b>	Q16655-MART-1; P14679-Tyrosinase; P40967-gp100
<b>Unigene</b>	154069-MART-1; 503555-Tyrosinase; 95972-gp100
<b>Immunogen</b>	Recombinant hMART-1 protein (A103); Recombinant tyrosinase protein (T311); Extract of pigmented melanoma metastases from lymph nodes (HMB45)
<b>Antibody target cellular localization</b>	Cytoplasmic
<b>Verified antibody applications</b>	IHC (FPPE) (verified), WB (verified)
<b>Species reactivity</b>	Human
<b>Antibody application notes</b>	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
<b>Positive control</b>	SK-MEL-13 and SK-MEL-19 Melanoma cell lines; Melanomas
<b>Shipping condition</b>	Room temperature
<b>Storage Conditions</b>	Store at 2 to 8 °C, Protect fluorescent conjugates from light, Note: store BSA-free antibodies at -10 to -35 °C
<b>Shelf life</b>	Guaranteed for at least 24 months from date of receipt when stored as recommended
<b>Regulatory status</b>	For research use only (RUO)
<b>Antibody/conjugate formulation</b>	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
<b>Antibody research areas</b>	Cancer
<b>Product origin</b>	Product may contain either bovine serum albumin (BSA) from bovine serum (Bos taurus), or recombinant BSA produced in Chinese hamster ovary cells. Inquire for the specific lot.
<b>Cell/tissue expression</b>	Melanocytes
<b>Tumor expression</b>	Melanoma

Antibody # prefix	Conjugation	Ex/Em (nm)	Laser line	Detection channel	Dye Features
BNC04	CF®405S	404/431	405	DAPI (microscopy), AF405	<a href="#">CF®405S Features</a>
BNC88	CF®488A	490/515	488	GFP, FITC	<a href="#">CF®488A Features</a>
BNC68	CF®568	562/583	532, 561	RFP, TRITC	<a href="#">CF®568 Features</a>
BNC94	CF®594	593/614	561	Texas Red®	<a href="#">CF®594 Features</a>
BNC40	CF®640R	642/662	633-640	Cy®5	<a href="#">CF®640R Features</a>
BNC47	CF®647	650/665	633-640	Cy®5	<a href="#">CF®647 Features</a>
BNC74	CF®740	742/767	633-685	775/50	<a href="#">CF®740 Features</a>
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 December 23, 2024 at 12:02:05 PM. Visit product page to check for updated information before use.

Product link: <https://biotium.com/product/monoclonal-anti-melanoma-marker-hmb45-a103-t311/>