

Melanoma Marker Monoclonal Mouse Antibody (DT101+ BC199 + T311 + HMB45)

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. 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. Catalog number key for antibody number 0704, Anti-Melanoma cells (DT101 BC199 T311 HMB45)

Call us: 800-304-5357

Product attributes

Product attributes				
Antibody number	#0704			
Antibody reactivity (target)	Melanoma cells			
Antibody type	Primary			
Host species	Mouse			
Clonality	Monoclonal			
Clone	DT101+ BC199 + T311 + HMB45			
Isotype	IgG2a, kappa			
Molecular weight	20-22 kDa (doublet) (MART-1); 90-100 kDa (gp100); 70-80 kDa (Tyrosinase)			
Synonyms	Melanoma antigen recognized by T-cells 1 (MART-1), MLAN-A; PMEL17 $$			
Entrez gene ID	2315; 6490			
Immunogen	Recombinant hMART-1 protein (DT101 & BC199); Extract of pigmented melanoma metastases from lymph nodes (HMB45); Recombinant tyrosinase protein (T311)			
Antibody target cellular localization	Cytoplasmic			
Verified antibody applications	IHC (FFPE) (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, Immunohistochemistry (formalin-fixed): 0.5-1.0 ug/mL for 30 minutes at RT, Staining of formalin-fixed tissues is enhanced by boiling tissue sections in 10 mM citrate buffer pH 6.0 for 10-20 minutes followed by cooling at RT for 20 minutes, Optimal dilution for a specific application should be determined by user			
5	determined by user			
Positive control	SK-MEL-13 and SK-MEL-19 melanoma cell lines. Human melanoma tissues.			
Shipping condition	SK-MEL-13 and SK-MEL-19 melanoma cell lines. Human			
	SK-MEL-13 and SK-MEL-19 melanoma cell lines. Human melanoma tissues.			
Shipping condition	SK-MEL-13 and SK-MEL-19 melanoma cell lines. Human melanoma tissues. Room temperature Store at 2 to 8 °C, Protect fluorescent conjugates from light,			
Shipping condition Storage Conditions	SK-MEL-13 and SK-MEL-19 melanoma cell lines. Human melanoma tissues. Room temperature Store at 2 to 8 °C, Protect fluorescent conjugates from light, Note: store BSA-free antibodies at -10 to -35 °C			
Shipping condition Storage Conditions Regulatory status Antibody/conjugate	SK-MEL-13 and SK-MEL-19 melanoma cell lines. Human melanoma tissues. Room temperature Store at 2 to 8 °C, Protect fluorescent conjugates from light, Note: store BSA-free antibodies at -10 to -35 °C 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/0.05% acide, Purified, BSA-free: 1 mg/mL in			
Shipping condition Storage Conditions Regulatory status Antibody/conjugate formulation	SK-MEL-13 and SK-MEL-19 melanoma cell lines. Human melanoma tissues. Room temperature Store at 2 to 8 °C, Protect fluorescent conjugates from light, Note: store BSA-free antibodies at -10 to -35 °C For research use only (RUO) 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 Guaranteed for at least 24 months from date of receipt when			
Shipping condition Storage Conditions Regulatory status Antibody/conjugate formulation Shelf life	SK-MEL-13 and SK-MEL-19 melanoma cell lines. Human melanoma tissues. Room temperature Store at 2 to 8 °C, Protect fluorescent conjugates from light, Note: store BSA-free antibodies at -10 to -35 °C For research use only (RUO) 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 Guaranteed for at least 24 months from date of receipt when stored as recommended			
Shipping condition Storage Conditions Regulatory status Antibody/conjugate formulation Shelf life Cell/tissue expression	SK-MEL-13 and SK-MEL-19 melanoma cell lines. Human melanoma tissues. Room temperature Store at 2 to 8 °C, Protect fluorescent conjugates from light, Note: store BSA-free antibodies at -10 to -35 °C 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 without azide Guaranteed for at least 24 months from date of receipt when stored as recommended Melanocytes Product may contain either bovine serum albumin (BSA) from bovine serum (Bos taurus), or recombinant BSA produced in			

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,	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 26, 2025 at 05:57:00 PM. Visit product page to check for updated information before use. Product link: https://biotium.com/product/melanoma-marker-monoclonal-mouse-antibody-dt101-bc199-t311-hmb45/