



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

## Product Description

This antibody cocktail recognizes two melanoma-specific proteins, which include MART-1 and gp100. MART-1 is a newly identified melanocyte differentiation antigen recognized by autologous cytotoxic T lymphocytes. Function of gp100 is not known but it is reported to be a useful marker for melanocytes and melanomas. A cocktail of these two markers is sensitive for labeling formalin-fixed, paraffin-embedded melanomas and other tumors showing melanocytic differentiation.

### Product attributes

Antibody number	#0703
Antibody reactivity (target)	Melanoma cells
Antibody type	Primary
Host species	Mouse
Clonality	Monoclonal
Clone	DT101+ BC199 + HMB45
Isotype	IgG1, kappa
Molecular weight	20-22 kDa (doublet) (MART-1); 90-100 kDa (gp100)
Synonyms	Melanoma antigen recognized by T-cells 1 (MART-1), MLAN-A; PMEL17
Human gene symbol	MLANA & SILV
Entrez gene ID	2315; 6490
SwissProt	Q16655 ; P40967
Unigene	154069 ; 95972
Immunogen	Recombinant hMART-1 protein (DT101 & BC199); Extract of pigmented melanoma metastases from lymph nodes (HMB45)
Antibody target cellular	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
Positive control	SK-MEL-13 and SK-MEL-19 melanoma cell lines. Human melanoma tissues.
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
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
Shelf life	Guaranteed for at least 24 months from date of receipt when stored as recommended
Cell/tissue expression	Melanocytes
Tumor expression	Melanoma
Antibody research areas	Cancer

This datasheet was generated on August 30, 2024 at 04:13:38 AM. Visit product page to check for updated information before use.  
 Product link: <https://biotium.com/product/melanoma-marker-monoclonal-mouse-antibody-dt101-bc199-hmb45/>