

gp100 / Melanosome / PMEL17 / SILV Monoclonal Mouse Antibody (HMB45 + PMEL/783)

Product Description

By immunohistochemistry, this antibody specifically recognizes a protein in melanocytes and melanomas. This MAb reacts with junctional and blue nevus cells and variably with fetal and neonatal melanocytes. Intradermal nevi, normal adult melanocytes, and non-melanocytic cells are negative. It does not stain tumor cells of epithelial, lymphoid, glial, or mesenchymal origin. Metastatic amelanotic melanoma can often be confused with a variety of poorly differentiated carcinomas, large cell lymphomas, and sarcomas using H E stains alone. It is also difficult to differentiate melanoma from spindle cell carcinomas and various types of mesenchymal neoplasms. This MAb stains fetal and neonatal melanocytes, junctional and blue nevus cells, and malignant melanoma. This MAb also stains Angiomyolipoma (PEComa). Catalog number key for antibody number 0951, Anti-gp100|Melanosome|Pmel17|SILV (HMB45 PMEL/783)

Call us: 800-304-5357

Product attributes

Antibody number				
Antibody number	#0951			
Antibody reactivity (target)	gp100, Melanosome, Pmel17, SILV			
Antibody type	Primary			
Host species	Mouse			
Clonality	Monoclonal			
Clone	HMB45 + PMEL/783			
Isotype	IgG1, kappa			
Molecular weight	90-100 kDa			
Synonyms	95kDa melanocyte-specific secreted glycoprotein, M-beta, Melanocyte lineage specific antigen GP100, Melanocyte protei Pmel 17, Melanoma associated ME20 antigen, Melanosomal matrix protein17, p100, p26, PMEL17, Premelanosome protein Secreted melanoma-associated ME20 antigen, SILV, Silver homolog			
Human gene symbol	SILV			
Entrez gene ID	6490			
SwissProt	P40967			
Unigene	95972			
Immunogen	Extract of pigmented melanoma metastases from lymph nodes (HMB45); Recombinant human SILV protein (PMEL/783)			
Antibody target cellular localization	Endoplasmic reticulum, Golgi apparatus			
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, Immunohistochemistry (formalin-fixed): 0.5-1.0 ug/mL for 30 minutes at RT, Staining of formalin-fixed tissues requires 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			
	minutes, Optimal dilution for a specific application should be			
Positive control	minutes, Optimal dilution for a specific application should be			
Positive control Shipping condition	minutes, Optimal dilution for a specific application should be determined by user			
	minutes, Optimal dilution for a specific application should be determined by user SK-MEL-28 cells. Melanoma.			
Shipping condition	minutes, Optimal dilution for a specific application should be determined by user SK-MEL-28 cells. Melanoma. Room temperature Store at 2 to 8 °C, Protect fluorescent conjugates from light,			
Shipping condition Storage Conditions	minutes, Optimal dilution for a specific application should be determined by user SK-MEL-28 cells. Melanoma. 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			
Shipping condition Storage Conditions Regulatory status Antibody/conjugate	minutes, Optimal dilution for a specific application should be determined by user SK-MEL-28 cells. Melanoma. 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, Purified: 0.2 mg/mL in PBS/0.05% BSA, Purified: 0.2 mg/mL in PBS/0.05% BSA/0.05% azide, Puri			
Shipping condition Storage Conditions Regulatory status Antibody/conjugate formulation	minutes, Optimal dilution for a specific application should be determined by user SK-MEL-28 cells. Melanoma. 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/0.05% azide, HRP conjugates: 0.1 mg/mL in PBS/0.05% 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	minutes, Optimal dilution for a specific application should be determined by user SK-MEL-28 cells. Melanoma. 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/0.05% BSA/0.05% azide, HRP conjugates: 0.1 mg/mL in PBS/0.05% BSA/0.05% azide, Purified; 0.2 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	minutes, Optimal dilution for a specific application should be determined by user SK-MEL-28 cells. Melanoma. 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, 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 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, 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 14, 2025 at 09:04:36 PM. Visit product page to check for updated information before use. Product link: https://biotium.com/product/gp100-melanosome-pmel17-silv-monoclonal-mouse-antibody-hmb45-pmel-783/