

# 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).

### Product attributes

<b>Antibody number</b>	#0951
<b>Antibody reactivity (target)</b>	gp100, Melanosome, Pmel17, SILV
<b>Antibody type</b>	Primary
<b>Host species</b>	Mouse
<b>Clonality</b>	Monoclonal
<b>Clone</b>	HMB45 + PMEL/783
<b>Isotype</b>	IgG1, kappa
<b>Molecular weight</b>	90-100 kDa
<b>Synonyms</b>	95kDa melanocyte-specific secreted glycoprotein, M-beta, Melanocyte lineage specific antigen GP100, Melanocyte protein Pmel 17, Melanoma associated ME20 antigen, Melanosomal matrix protein17, p100, p26, PMEL17, Premelanosome protein, Secreted melanoma-associated ME20 antigen, SILV, Silver homolog
<b>Human gene symbol</b>	SILV
<b>Entrez gene ID</b>	6490
<b>SwissProt</b>	P40967
<b>Unigene</b>	95972
<b>Immunogen</b>	Extract of pigmented melanoma metastases from lymph nodes (HMB45); Recombinant human SILV protein (PMEL/783)
<b>Antibody target cellular localization</b>	Endoplasmic reticulum, Golgi apparatus
<b>Verified antibody applications</b>	IHC (FFPE) (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. 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
<b>Positive control</b>	SK-MEL-28 cells. Melanoma.
<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>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>Shelf life</b>	Guaranteed for at least 24 months from date of receipt when stored as recommended
<b>Cell/tissue expression</b>	Melanocytes
<b>Tumor expression</b>	Melanoma
<b>Antibody research areas</b>	Cancer