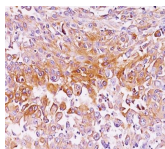


# gp100 / Melanosome / PMEL17 / SILV Monoclonal Mouse Antibody (NKI-beteb)



## Product Description

By immunohistochemistry, this antibody specifically recognizes a protein in melanocytes and melanomas. It reacts with junctional and blue nevus cells and variably with fetal and neonatal melanocytes. Intra-dermal nevi, normal adult melanocytes, and non-melanocytic cells are negative. It does not stain tumor cells of epithelial, lymphoid, glial, or mesenchymal origin. This Mab labels 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 0226, Anti-gp100 (NKI-beteb)**

Call us : [800-304-5357](tel:800-304-5357)

## Product attributes

<b>Antibody number</b>	#0226
<b>Antibody reactivity (target)</b>	gp100, Pmel17, SILV
<b>Antibody type</b>	Primary
<b>Host species</b>	Mouse
<b>Clonality</b>	Monoclonal
<b>Clone</b>	NKI-beteb
<b>Isotype</b>	IgG2b, kappa
<b>Molecular weight of antigen</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>	PMEL
<b>Entrez gene ID</b>	6490
<b>SwissProt</b>	P40967
<b>Unigene</b>	95972
<b>Immunogen</b>	Membranes from a human melanoma metastasis
<b>Verified antibody applications</b>	IHC (FFPE) (verified)
<b>Antibody target cellular localization</b>	Endoplasmic reticulum, Golgi apparatus
<b>Expected antibody applications</b>	ELISA (published for clone), Flow (intracellular) (published for clone), IHC (frozen) (published for clone), IF (published for clone), IP (published for clone)
<b>Species reactivity</b>	Horse, Human
<b>Antibody application notes</b>	Immunohistology formalin-fixed 1-2 ug/mL, Staining of formalin-fixed tissues requires boiling tissue sections in 10 mM citrate buffer, pH 6.0, for 10-20 min followed by cooling at RT for 20 minutes, Flow Cytometry 0.5-1 ug/million cells/0.1 mL, Immunofluorescence 1-2 ug/mL, Does not react with rat, others not tested, Optimal dilution for a specific application should be determined by user
<b>Positive control</b>	SK-MEL-28 cells or 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>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.

## References

Note: References for this clone sold by other suppliers may be listed for expected applications.

1. Am J Pathol (1988) 130(1): 179. (mAb characterization; IHC, FFPE; IF; ELISA; IP; WB)
2. Am J Pathol (1993) 143(6): 1579. (IHC, frozen; IF)
3. J Biol Chem (2004) 279(27): 28330-28338. (epitope mapping; IP; reports that NK1-beteb does not work well for WB)
4. Exp Dermatol (2010) 19(8): e282-e285. (IF; Flow, intracellular)

This datasheet was generated on June 4, 2026 at 12:09:54 PM. Visit product page to check for updated information before use.

Product link: <https://biotium.com/product/monoclonal-anti-pmel17-gp100-silv-nki-beteb/>