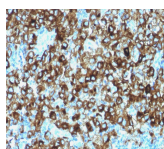


gp100 / Melanosome / PMEL17 / SILV Polyclonal Rabbit Antibody



Product Description

Cytotoxic T lymphocytes (CTLs) recognize melanoma-associated antigens, which belong to three main groups. These groups include tumor-associated testis-specific antigens, melanocyte differentiation antigens and mutated or aberrantly expressed antigens, which are routinely used as markers to identify melanomas based on their binding to specific monoclonal antibodies. gp100, also designated ME20-M, ME20-S and PMEL 17, is classified as a melanocyte differentiation antigen and is expressed at low levels in normal cell lines and tissues, but is upregulated in melanocytes. gp100 is a highly glycosylated protein. It is also the product of proteolytic cleavage, which results in a secreted protein.

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 to inquire about stock status and lead times before placing your order.

Catalog number key for antibody number 1826, gp100 / Melanosome / PMEL17 / SILV Polyclonal Rabbit Antibody

Product attributes

Antibody number	#1826
Antibody reactivity (target)	gp100, Melanosome, Pmel17, SILV
Antibody type	Primary
Host species	Rabbit
Clonality	Polyclonal
Clone	N/A
Isotype	IgG
Molecular weight of antigen	90-100 kDa
Synonyms	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
Human gene symbol	SILV
Entrez gene ID	6490
SwissProt	P40967
Unigene	95972
Immunogen	Recombinant full-length human SILV protein
Antibody target cellular localization	Endoplasmic reticulum, Golgi apparatus
Verified antibody applications	IHC (FFPE) (verified), WB (verified)
Species reactivity	Human
Positive control	SK-MEL-28 cells, Melanoma.
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
Product origin	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.
Cell/tissue expression	Melanocytes
Tumor expression	Melanoma

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 June 4, 2026 at 10:49:41 AM. Visit product page to check for updated information before use. Product link: <https://biotium.com/product/gp100-melanosome-pmel17-silv-polyclonal-rabbit-antibody-n-a/>