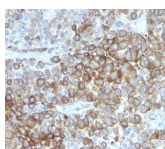


Melanoma Marker Monoclonal Mouse Antibody (PNL2)



Product Description

Anti-PNL2 is a novel monoclonal antibody, which has recently been introduced as an immunohistochemical reagent to stain melanocytes and tumors derived therefrom. The antigen recognized by PNL2 is different from Melan A and gp100. Its epitope is not destroyed by digestion with neuraminidase i.e. its epitope is not glycosylated. Anti-PNL2 may be most useful because of its high sensitivity for metastatic melanoma (87%), as opposed to 76% for anti-HMB45 and 82% for anti-MART-1. Anti-PNL2 labels intra-epidermal nevi while the dermal component of compound nevi are largely non-reactive with anti-PNL2. Antibodies against PNL2, MART-1 (Melan A) and HMB45 stain most clear cell sarcoma cells and a few cells in angio-myolipomas and lymphangioid leiomyomatosis. Anti-PNL2 is a useful antibody for the identification of melanomas and clear cell sarcomas. Differential diagnosis is aided by the results from a panel of antibodies, including antibodies against HMB45, MART-1, tyrosinase, and MITF.

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 0894, Anti-Melanoma Marker (PNL2)

Product attributes

Antibody number	#0894
Antibody reactivity (target)	Melanoma Marker
Antibody type	Primary
Host species	Mouse
Clonality	Monoclonal
Clone	PNL2
Isotype	IgG1, kappa
Molecular weight of antigen	Multiple (160 kDa, 100 kDa and
Synonyms	Human Melanoma Associated Antigen; PNL2
Entrez gene ID	Not Known
SwissProt	Not Known
Unigene	Not Known
Immunogen	Melanocyte antigen
Antibody target cellular localization	Cytoplasmic
Verified antibody applications	IHC (FFPE) (verified)
Species reactivity	Dog, Human, Mouse
Antibody application notes	Higher concentration may be required for direct detection using primary antibody conjugates than for indirect detection with secondary antibody. Immunofluorescence: 0.5-1 µg/mL. Immunohistochemistry (formalin). 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. Optimal dilution for a specific application should be determined by user
Positive control	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
Shelf life	Guaranteed for at least 24 months from date of receipt when stored as recommended
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
Antibody research areas	Cancer
Product origin	Product may contain either bovine serum albumin (BSA) from bovine serum (<i>Bos taurus</i>), 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 24, 2026 at 08:34:35 PM. Visit product page to check for updated information before use. Product link: <https://biotium.com/product/monoclonal-anti-melanoma-marker-pnl2/>