

Tyrosinase Related Protein-1 (TYRP1) Monoclonal Mouse Antibody (TA99)



Product Description

This antibody reacts with a 75 kDa melanocyte-specific gene product, identified as Tyrosinase-related protein-1 (TRP-1). It is involved in melanin synthesis. TRP1 is present on the melanosomal membranes of melanoma, normal melanocytes and nevi.Recent evidence suggests that TRP-1 is involved in maintaining stability of tyrosinase protein and modulating its catalytic activity. TRP-1 is also involved in maintenance of melanosome ultrastructure and affects melanocyte proliferation and cell death.

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 0806, Anti-TRP1 (TA99)

Call us: 800-304-5357

Product attributes

Antibody number	#0806		
Antibody reactivity (target)	TRP1		
Antibody type	Primary		
Host species	Mouse		
Clonality	Monoclonal		
Clone	TA99		
Isotype	IgG2a, kappa		
Molecular weight	75 kDa		
Synonyms	5, 6 dihydroxyindole 2 carboxylic acid oxidase, 6-dihydroxyindole-2-carboxylic acid oxidase, Associated with iris pigmentation, CAS2, Catalase B (CATB), DHICA oxidase, Glycoprotein75 (GP75), Melanoma antigen gp75, Tyrosinase-related protein 1 (TYRP1), TYRRP		
Human gene symbol	TYRP1		
Entrez gene ID	7306		
SwissProt	P17643		
Unigene	270279		
Immunogen	SK-MEL-23 cells		
Antibody target cellular localization	Vesicular		
Expected antibody applications	Flow, surface (published for clone), IHC (FFPE) (published for clone), IHC (frozen) (published for clone), IF (published for clone), IP (published for clone)		
Species reactivity	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: 1-2 ug/mL, Immunohistology (frozen only) 0.5-1 ug/mL, Flow Cytometry 0.5-1 ug/million cells/0.1 mL, Optimal dilution for a specific application should be determined by user		
	0.5-1 ug/million cells/0.1 mL, Optimal dilution for a specific		
Positive control	0.5-1 ug/million cells/0.1 mL, Optimal dilution for a specific		
Positive control Shipping condition	0.5-1 ug/million cells/0.1 mL, Optimal dilution for a specific application should be determined by user		
	0.5-1 ug/million cells/0.1 mL, Optimal dilution for a specific application should be determined by user SK-MEL-23, SK-MEL-19, SK-MEL-30 cells or Melanoma.		
Shipping condition	0.5-1 ug/million cells/0.1 mL, Optimal dilution for a specific application should be determined by user SK-MEL-23, SK-MEL-19, SK-MEL-30 cells or Melanoma. Room temperature Store at 2 to 8 °C, Protect fluorescent conjugates from light,		
Shipping condition Storage Conditions	0.5-1 ug/million cells/0.1 mL, Optimal dilution for a specific application should be determined by user SK-MEL-23, SK-MEL-19, SK-MEL-30 cells or Melanoma. Room temperature Store at 2 to 8 °C, Protect fluorescent conjugates from light, Note: store BSA-free antibodies at -10 to -35 °C Guaranteed for at least 24 months from date of receipt when		
Shipping condition Storage Conditions Shelf life	0.5-1 ug/million cells/0.1 mL, Optimal dilution for a specific application should be determined by user SK-MEL-23, SK-MEL-19, SK-MEL-30 cells or Melanoma. Room temperature Store at 2 to 8 °C, Protect fluorescent conjugates from light, Note: store BSA-free antibodies at -10 to -35 °C Guaranteed for at least 24 months from date of receipt when stored as recommended		
Shipping condition Storage Conditions Shelf life Regulatory status Antibody/conjugate	0.5-1 ug/million cells/0.1 mL, Optimal dilution for a specific application should be determined by user SK-MEL-23, SK-MEL-19, SK-MEL-30 cells or Melanoma. Room temperature Store at 2 to 8 °C, Protect fluorescent conjugates from light, Note: store BSA-free antibodies at -10 to -35 °C Guaranteed for at least 24 months from date of receipt when stored as recommended 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/0.05% azide, Purified: 0.2 mg/mL in PBS/0.05% BSA/0.05% azide, Purified: 0.2 mg/mL in PBS/0.05% BSA/0.05% azide, Purified: 0.58A-free: 1 mg/mL in PBS/0.05% BSA/0.05% azide, Purified; 0.58A-free: 1 mg/mL in PBS/0.05% BSA/0.05% azide, Purified; 0.58A-free: 1 mg/mL in PBS/0.05% BSA/0.05% azide, Purified, BSA-free: 1 mg/mL in PBS/0.05% BSA/0.05% BSA/0.		
Shipping condition Storage Conditions Shelf life Regulatory status Antibody/conjugate formulation	0.5-1 ug/million cells/0.1 mL, Optimal dilution for a specific application should be determined by user SK-MEL-23, SK-MEL-19, SK-MEL-30 cells or Melanoma. Room temperature Store at 2 to 8 °C, Protect fluorescent conjugates from light, Note: store BSA-free antibodies at -10 to -35 °C Guaranteed for at least 24 months from date of receipt when stored as recommended 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/0.05% azide, Purified, BSA-free: 1 mg/mL in PBS without azide		
Shipping condition Storage Conditions Shelf life Regulatory status Antibody/conjugate formulation Antibody research areas	0.5-1 ug/million cells/0.1 mL, Optimal dilution for a specific application should be determined by user SK-MEL-23, SK-MEL-19, SK-MEL-30 cells or Melanoma. Room temperature Store at 2 to 8 °C, Protect fluorescent conjugates from light, Note: store BSA-free antibodies at -10 to -35 °C Guaranteed for at least 24 months from date of receipt when stored as recommended 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/0.05% azide, Purified; BSA-free: 1 mg/mL in PBS without azide Cancer 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.

References

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

- J Invest Dermatol (1985) 85:169-174. (IP; IHC, frozen)
 PNAS USA (1987) 84, 4200-4204. (in vivo imaging melanoma)
 J Exp Med (1995) 182: 1609-1614. (functional studies)
 J Cell Biol (2015) 209(4): 563-577. (Flow, surface; IF)
 J Cell Mol Med (2017) 21(4): 802-815. (IHC, FFPE)

This datasheet was generated on November 30, 2025 at 03:46:28 AM. Visit product page to check for updated information before use. Product link: https://biotium.com/product/monoclonal-anti-trp1-tyrosinase-related-protein-1-ta99/