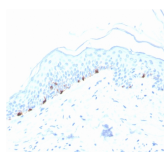


# Tyrosinase Related Protein-1 (TYRP1) Monoclonal Mouse Antibody (TYRP1/3280)



## Product Description

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](mailto:order@biotium.com) to inquire about stock status and lead times before placing your order.

**Catalog number key for antibody number 3280, Anti-Tyrosinase Related Protein-1 (TYRP1/3280)**

## Product attributes

Antibody number	#3280
Antibody reactivity (target)	Tyrosinase Related Protein-1
Antibody type	Primary
Host species	Mouse
Clonality	Monoclonal
Clone	TYRP1/3280
Isotype	IgG2b, 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
Human gene symbol	TYRP1
Entrez gene ID	7306
SwissProt	P17643
Unigene	270279
Immunogen	A recombinant fragment (around aa 257-377) of human TYRP1 protein (exact sequence is proprietary)
Verified antibody applications	IHC (FFPE) (verified)
Antibody target cellular localization	Vesicular
Species reactivity	Human
Positive control	SK-MEL-23, SK-MEL-19, SK-MEL-30, SK-MEL-37 cells. Human skin or melanoma tissues (IHC).
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
Validated in protein array	Monospecific
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	<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.

This datasheet was generated on August 7, 2025 at 07:26:13 PM. Visit product page to check for updated information before use.  
Product link: <https://biotium.com/product/tyrosinase-related-protein-1-tyrp1-monoclonal-mouse-antibody-tyrp1-3280/>