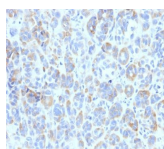


# Tyrosinase Related Protein-1 (TYRP1) Monoclonal Mouse Antibody (TA99)



## 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](mailto: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)**

## 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), TYRPP   |
| 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 |
| Positive control                      | SK-MEL-23, SK-MEL-19, SK-MEL-30 cells or 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 (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,<br>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. J Invest Dermatol (1985) 85:169-174. (IP; IHC, frozen)
2. PNAS USA (1987) 84, 4200-4204. (in vivo imaging melanoma)
3. J Exp Med (1995) 182: 1609-1614. (functional studies)
4. J Cell Biol (2015) 209(4): 563-577. (Flow, surface; IF)
5. J Cell Mol Med (2017) 21(4): 802-815. (IHC, FFPE)

This datasheet was generated on August 18, 2025 at 10:48:30 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/>