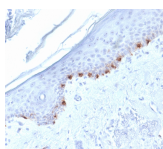


Tyrosinase Related Protein-1 (TYRP1) Monoclonal Mouse Antibody (TYRP1/1986)



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 1986, Anti-Tyrosinase Related Protein-1 (TYRP1/1986)

Product attributes

| | |
|---------------------------------------|--|
| Antibody number | #1986 |
| Antibody reactivity (target) | Tyrosinase Related Protein-1 |
| Antibody type | Primary |
| Host species | Mouse |
| Clonality | Monoclonal |
| Clone | TYRP1/1986 |
| 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) |
| Antibody target cellular localization | Vesicular |
| Verified antibody applications | IHC (FFPE) (verified) |
| Species reactivity | Human |
| Positive control | SK-MEL-23, SK-MEL-19, SK-MEL-30, SK-MEL-37 cells. Human Skin 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 |
| Validated in protein array | Monospecific |
| 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 | 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 August 26, 2025 at 11:15:44 AM. Visit product page to check for updated information before use.
Product link: <https://biotium.com/product/tyrosinase-related-protein-1-monoclonal-mouse-antibody-tyrp1-1986/>