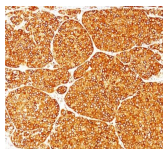


Tyrosinase Monoclonal Mouse Antibody (T311)



Product Description

Recognizes a cluster of proteins between 70-80 kDa, identified as tyrosinase. Occasionally a minor band at 55 kDa is also detected. This MAb shows no cross-reaction with MAGE-1 and tyrosinase-related protein 1, TRP-1/gp75. Tyrosinase is a copper-containing metalloglycoprotein that catalyzes several steps in the melanin pigment biosynthetic pathway; the hydroxylation of tyrosine to L-3,4-dihydroxy-phenylalanine (dopa), and the subsequent oxidation of dopa to dopaquinone. Mutations of the tyrosinase gene occur in various forms of albinism. Tyrosinase is one of the targets for cytotoxic T-cell recognition in melanoma patients. Staining of melanomas with this MAb shows tyrosinase in melanotic as well as amelanotic variants. This MAb is a useful marker for melanocytes and melanomas.

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 0012, Anti-Tyrosinase (T311)

Call us : [800-304-5357](tel:800-304-5357)

Product attributes

| | |
|---------------------------------------|--|
| Antibody number | #0012 |
| Antibody reactivity (target) | Tyrosinase |
| Antibody type | Primary |
| Host species | Mouse |
| Clonality | Monoclonal |
| Clone | T311 |
| Isotype | IgG2a, kappa |
| Molecular weight | 70-80 kDa |
| Synonyms | ATN, CMM8, LB24-AB, Monophenol monooxygenase, OCA1, OCA1A, Oculocutaneous albinism IA, SHEP3, SK29-AB, Tumor rejection antigen AB, TYR |
| Human gene symbol | TYR |
| Entrez gene ID | 7299 |
| SwissProt | P14679 |
| Unigene | 503555 |
| Immunogen | Recombinant tyrosinase protein |
| Antibody target cellular localization | Cytoplasmic |
| Verified antibody applications | IHC (FFPE) (verified) |
| Species reactivity | Human |
| Expected antibody applications | ELISA (published for clone), Flow (intracellular) (published for clone), IP (published for clone), WB (published for clone) |
| 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 formalin-fixed 0.5-1 ug/mL, Staining of formalin-fixed tissues requires boiling tissue sections in 10 mM Tris with 1 mM EDTA, pH 9.0, or 10 mM citrate buffer, pH6.0, for 10-20 min followed by cooling at RT for 20 minutes, 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-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 | 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 Clin Pathol (2001) 54:196-200. (IHC, FFPE)
- PNAS USA (1995) 92: 8125-8129. (WB, IHC, ELISA)
- J Biol Chem (2001) 276(15): 11933-11938. (WB, IP)
- Cytometry Part A (2012) 81A: 374381. (Flow, intracellular)

This datasheet was generated on January 21, 2026 at 03:18:43 AM. Visit product page to check for updated information before use.
Product link: <https://biotium.com/product/monoclonal-anti-tyrosinase-t311/>