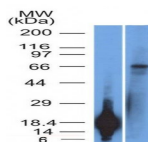


TRIM29 Monoclonal Mouse Antibody (TRIM29/1042)



Product Description

This antibody recognizes a 66 kDa protein, which is identified as Tripartite motif-containing protein 29 (TRIM29). It interacts with the intermediate filament protein vimentin, a substrate for the PKC family of protein kinases, and with hPKC δ -1, an inhibitor of the PKCs. TRIM29 protein contains both zinc finger and leucine zipper motifs, suggesting that it may form homodimers and possibly associate with DNA. High expression of TRIM29 has been reported in gastric cancer and pancreatic cancer, and correlates with enhanced tumor growth and lymph node metastasis. TRIM29 is also able to distinguish lung squamous cell carcinoma from lung adenocarcinoma with ~90% positive accuracy, when used in a panel with TTF-1, p63, CK5/6, and Napsin-A antibodies.

Primary antibodies are available purified, or with a selection of fluorescent CF $\text{\textcircled{R}}$ dyes and other labels. CF $\text{\textcircled{R}}$ dyes offer exceptional brightness and photostability. See the [CF \$\text{\textcircled{R}}\$ Dye Brochure](#) for more information. Note: Conjugates of blue fluorescent dyes like CF $\text{\textcircled{R}}$ 405S and CF $\text{\textcircled{R}}$ 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 $\text{\textcircled{R}}$ 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 1042, Anti-TRIM29 (TRIM29/1042)

Product attributes

| | |
|---------------------------------------|---|
| Antibody number | #1042 |
| Antibody reactivity (target) | TRIM29 |
| Antibody type | Primary |
| Host species | Mouse |
| Clonality | Monoclonal |
| Clone | TRIM29/1042 |
| Isotype | IgG2b, kappa |
| Molecular weight | 66 kDa |
| Synonyms | Ataxia telangiectasia group D complementing gene (ATDC); Tripartite motif-containing protein 29 (TRIM29); |
| Human gene symbol | TRIM29 |
| Entrez gene ID | 23650 |
| SwissProt | Q14134 |
| Unigene | 504115 |
| Immunogen | Recombinant fragment (126 Amino acid residues between aa 1-200) of human TRIM29 protein |
| Verified antibody applications | WB (verified) |
| Antibody target cellular localization | Cytoplasmic, Cytoskeleton, Nucleus |
| Species reactivity | Human |
| Antibody application notes | Higher concentration may be required for direct detection using primary antibody conjugates than for indirect detection with secondary antibody, Immunofluorescence: 0.5-1 μ g/mL, Flow Cytometry 0.5-1 μ g/million cells/0.1 mL, Western blotting 0.5-1 μ g/mL, Optimal dilution for a specific application should be determined by user |
| Positive control | A431 cells, Tonsil or Squamous cell carcinoma. |
| Shipping condition | Room temperature |
| Storage Conditions | Store at 2 to 8 $^{\circ}$ C, Protect fluorescent conjugates from light, Note: store BSA-free antibodies at -10 to -35 $^{\circ}$ 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, Cytoskeleton, Signal transduction |
| 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. |
| Tumor expression | Lung cancer, Squamous cell carcinoma |

| Antibody # prefix | Conjugation | Ex/Em (nm) | Laser line | Detection channel | Dye Features |
|-------------------|----------------------------------|------------|------------|------------------------------------|--|
| BNC04 | CF $\text{\textcircled{R}}$ 405S | 404/431 | 405 | DAPI (microscopy), AF405 | CF$\text{\textcircled{R}}$405S Features |
| BNC88 | CF $\text{\textcircled{R}}$ 488A | 490/515 | 488 | GFP, FITC | CF$\text{\textcircled{R}}$488A Features |
| BNC68 | CF $\text{\textcircled{R}}$ 568 | 562/583 | 532, 561 | RFP, TRITC | CF$\text{\textcircled{R}}$568 Features |
| BNC94 | CF $\text{\textcircled{R}}$ 594 | 593/614 | 561 | Texas Red $\text{\textcircled{R}}$ | CF$\text{\textcircled{R}}$594 Features |
| BNC40 | CF $\text{\textcircled{R}}$ 640R | 642/662 | 633-640 | Cy $\text{\textcircled{R}}$ 5 | CF$\text{\textcircled{R}}$640R Features |
| BNC47 | CF $\text{\textcircled{R}}$ 647 | 650/665 | 633-640 | Cy $\text{\textcircled{R}}$ 5 | CF$\text{\textcircled{R}}$647 Features |
| BNC74 | CF $\text{\textcircled{R}}$ 740 | 742/767 | 633-685 | 775/50 | CF$\text{\textcircled{R}}$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 September 15, 2025 at 01:24:09 PM. Visit product page to check for updated information before use.
Product link: <https://biotium.com/product/monoclonal-anti-trim29-trim291042/>