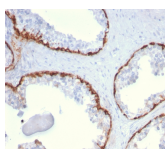


# Cytokeratin 15 Monoclonal Mouse Antibody (KRT15/2959)



## Product Description

Keratin 15 is a type I keratin which is expressed only in basal keratinocytes in stratified epithelia and does not appear to have a natural type II expression partner. Keratin 15 is down regulated in activated keratinocytes. Cytokeratin 15 is a specific marker of stem cells of the hair-follicle bulge and may be a useful marker for diagnosis between basal cell carcinoma (BCC) and trichoepithelioma. Trichoblastoma are benign neoplasms of follicular differentiation frequently found in nevus sebaceous. Many morphologic features are shared with nodular basal cell carcinoma, sometimes rendering a diagnosis difficult. Trichoblastoma and BCC show variable expression of Cytokeratin 15 and Cytokeratin 19, and absence of hair keratins.

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 2959, Anti-Cytokeratin 15 (KRT15/2959)**

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

## Product attributes

|  |   |
|--|---|
| <b>Antibody number</b>                       | #2959   |
| <b>Antibody reactivity (target)</b>          | Cytokeratin 15  |
| <b>Antibody type</b>                         | Primary   |
| <b>Host species</b>                          | Mouse   |
| <b>Clonality</b>                             | Monoclonal  |
| <b>Clone</b>                                 | KRT15/2959  |
| <b>Isotype</b>                               | IgG2c, kappa  |
| <b>Molecular weight of antigen</b>           | 52 kDa  |
| <b>Synonyms</b>                              | CK15; Cytokeratin 15; K1CO; Ka15; Keratin 15 basic; Keratin 15 beta; Keratin complex 1 acidic gene 15; Keratin type I cytoskeletal 15; KRT15; KRTB; KRTL15; Type I cytoskeletal 15; Type I keratin Ka15 |
| <b>Human gene symbol</b>                     | KRT15   |
| <b>Entrez gene ID</b>                        | 3866  |
| <b>SwissProt</b>                             | P19012  |
| <b>Unigene</b>                               | 654570  |
| <b>Immunogen</b>                             | Recombinant full-length human KRT15 protein   |
| <b>Antibody target cellular localization</b> | Cytoskeleton  |
| <b>Verified antibody applications</b>        | IF (verified), IHC (FFPE) (verified), WB (verified)   |
| <b>Species reactivity</b>                    | Human   |
| <b>Positive control</b>                      | MCF-7, A431 cells. Skin or Colon or Thymus.   |
| <b>Shipping condition</b>                    | Room temperature  |
| <b>Storage Conditions</b>                    | Store at 2 to 8 °C, Protect fluorescent conjugates from light, Note: store BSA-free antibodies at -10 to -35 °C   |
| <b>Regulatory status</b>                     | For research use only (RUO)   |
| <b>Antibody/conjugate formulation</b>        | 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       |
| <b>Validated in protein array</b>            | Monospecific  |
| <b>Shelf life</b>                            | Guaranteed for at least 24 months from date of receipt when stored as recommended   |
| <b>Product origin</b>                        | 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.                         |
| <b>Cell/tissue expression</b>                | Epidermal cells   |
| <b>Antibody research areas</b>               | Cancer, Cytoskeleton  |

| 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 June 26, 2026 at 02:29:48 AM. Visit product page to check for updated information before use.

Product link: <https://biotium.com/product/cytokeratin-15-monoclonal-mouse-antibody-krt15-2959/>