

## Wilm's Tumor 1 / WT1 Recombinant Monoclonal Rabbit Antibody (WT1/1434R)

## **Product Description**

This antibody recognizes a 47-55 kDa-tumor suppressor protein, identified as Wilm's Tumor (WT1) protein. The antibody reacts with all isoforms of the full-length WT1 and also identifies WT1 lacking exon 2-encoded amino acids, frequently found in subsets of sporadic Wilm's tumors. WT1, a sporadic and familial pediatric kidney tumor, is genetically heterogeneous. Wilm's tumor is associated with mutations of WT1, a zinc-finger transcription factor that is essential for the development of the metanephric kidney and the urogenital system. The WT1 gene is normally expressed in fetal kidney and mesothelium, and its expression has been suggested as a marker for Wilm's tumor and mesothelioma. WT1 protein has been identified in proliferative mesothelial cells, malignant mesothelioma, ovarian carcinoma, gonadoblastoma, nephroblastoma, and desmoplastic small round cell tumor. Lung adenocarcinomas rarely stain positive with this antibody. WT1 protein expression in mesothelial cells has become a reliable marker for the diagnosis of mesotheliomas.

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 <a href="mailto:order@biotium.com">order@biotium.com</a> to inquire about stock status and lead times before placing your order.

Catalog number key for antibody number 1434, Anti-Wilm's Tumor 1 (WT1/1434R)

## Product attributes

Call us: 800-304-5357

| Product attributes                       |  |  |  |
|--|--|--|--|
| Antibody number                          | #1434  |  |  |
| Antibody reactivity (target)             | Wilm's Tumor 1, WT1  |  |  |
| Antibody type                            | Primary  |  |  |
| Host species                             | Rabbit   |  |  |
| Clonality                                | Recombinant Monoclonal   |  |  |
| Clone                                    | WT1/1434R  |  |  |
| Isotype                                  | IgG, kappa   |  |  |
| Molecular weight                         | 47-55 kDa  |  |  |
| Synonyms                                 | WT1; AWT1; FWT1; GUD; NPHS4; WAGR; Wilms tumor 1   |  |  |
| Human gene symbol                        | WT1  |  |  |
| Entrez gene ID                           | 7490   |  |  |
| SwissProt                                | P19544   |  |  |
| Unigene                                  | 591980   |  |  |
| Immunogen                                | Recombinant human full-length WT1 protein  |  |  |
| Antibody target cellular<br>localization | Nucleus & cytoplasm  |  |  |
| Verified antibody applications           | IHC (FFPE) (verified)  |  |  |
| 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 ug/mL, Immunohistology (formalin): 0.5-1 ug/mL, Staining of formalin-fixed tissues requires boiling tissue sections in 10 mM citrate buffer, pH 6.0, for 10-20 min followed by cooling at RT for 20 min, Flow Cytometry 0.5-1 ug/million cells/0.1 mL, Optimal dilution for a specific application should be determined by user |  |  |
| Positive control                         | K562 cells. Wilm s Tumor, mesothelioma or fetal kidney.  |  |  |
| 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<br>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, Developmental biology, Transcription factors   |  |  |
| 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                         | Mesothelioma   |  |  |
|  |  |  |  |

| Antibody # prefix | Conjugation           | Ex/Em (nm) | Laser line | Detection channel           | Dye Features     |
|-------------------|-----------------------|------------|------------|-----------------------------|------------------|
| BNC04             | CF®405S               | 404/431    | 405        | DAPI (microscopy),<br>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,<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.

This datasheet was generated on November 23, 2025 at 02:16:46 PM. Visit product page to check for updated information before use. Product link: <a href="https://biotium.com/product/recombinant-rabbit-monoclonal-anti-wilms-tumor-1-wt11434r/">https://biotium.com/product/recombinant-rabbit-monoclonal-anti-wilms-tumor-1-wt11434r/</a>