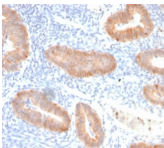


# Interleukin 10 Recombinant Monoclonal Rabbit Antibody (IL10/2651R)



## Product Description

Interleukin 10 (IL10) is a cytokine produced primarily by monocytes and to a lesser extent by lymphocytes. This cytokine has pleiotropic effects in immuno-regulation and inflammation. It down-regulates the expression of Th1 cytokines, MHC class II antigens, and costimulatory molecules on macrophages. It also enhances B cell survival, proliferation, and antibody production. This cytokine can block NF-kappa B activity, and is involved in the regulation of the JAK-STAT signaling pathway. Knockout studies in mice suggested the function of this cytokine as an essential immuno-regulator in the intestinal tract. Mutations in this gene are associated with an increased susceptibility to HIV-1 infection and rheumatoid arthritis. 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 2651, Anti-Interleukin 10 (IL10/2651R)**

| Product attributes                    |                                                                                                                                                                                                                                                                                                                                                                                                                                                        |
|---------------------------------------|--------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------|
| Antibody number                       | #2651                                                                                                                                                                                                                                                                                                                                                                                                                                                  |
| Antibody reactivity (target)          | Interleukin 10                                                                                                                                                                                                                                                                                                                                                                                                                                         |
| Antibody type                         | Primary                                                                                                                                                                                                                                                                                                                                                                                                                                                |
| Host species                          | Rabbit                                                                                                                                                                                                                                                                                                                                                                                                                                                 |
| Clonality                             | Recombinant Monoclonal                                                                                                                                                                                                                                                                                                                                                                                                                                 |
| Clone                                 | IL10/2651R                                                                                                                                                                                                                                                                                                                                                                                                                                             |
| Isotype                               | IgG                                                                                                                                                                                                                                                                                                                                                                                                                                                    |
| Molecular weight                      | 20 kDa (Monomer); 37 kDa (Dimer)                                                                                                                                                                                                                                                                                                                                                                                                                       |
| Synonyms                              | Cytokine Synthesis Inhibitory Factor (CSIF); GVHDS; IL-10; IL10; IL10A; Interleukin-10; T-cell Growth Inhibitory Factor (TGIF)                                                                                                                                                                                                                                                                                                                         |
| Human gene symbol                     | IL10                                                                                                                                                                                                                                                                                                                                                                                                                                                   |
| Entrez gene ID                        | 3586                                                                                                                                                                                                                                                                                                                                                                                                                                                   |
| SwissProt                             | P22301                                                                                                                                                                                                                                                                                                                                                                                                                                                 |
| Unigene                               | 193717                                                                                                                                                                                                                                                                                                                                                                                                                                                 |
| Immunogen                             | Recombinant full-length human Interleukin 10 (IL10) protein                                                                                                                                                                                                                                                                                                                                                                                            |
| Verified antibody applications        | IHC (FFPE) (verified)                                                                                                                                                                                                                                                                                                                                                                                                                                  |
| Antibody target cellular localization | Secreted (extracellular)                                                                                                                                                                                                                                                                                                                                                                                                                               |
| 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. Immunohistology (formalin): 1-2 ug/mL for 30 minutes at RT. Staining of formalin-fixed tissues requires boiling tissue sections in 10 mM citrate buffer, pH 6.0, for 10-20 minutes followed by cooling at RT for 20 minutes. Optimal dilution for a specific application should be determined by user |
| Positive control                      | SK-BR3 Cells, Adrenal, Lymph Node, Bladder, Prostate or Endometrium.                                                                                                                                                                                                                                                                                                                                                                                   |
| 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               | Cytokines, Immunology, Inflammation                                                                                                                                                                                                                                                                                                                                                                                                                    |

| 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>  |
| 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, LI-COR Bioscience.