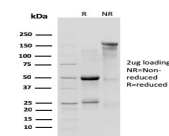


IL-10R1, Mouse Monoclonal Rat Antibody (1B1.3a)



Product Description

This monoclonal antibody recognizes Mouse Interleukin-10 receptor 1 (IL-10R1), a cell surface glycoprotein of approximately 110 kDa. Its epitope is localized in the ligand-binding domain in the extracellular region of Mouse IL-10R1. This antibody blocks the binding of IL-10 to its receptor. IL-10R1 is a class II cytokine receptor expressed on thymocytes, T cells, B cells, NK cells, monocytes and macrophages. Binding of IL-10 to its receptor initiates JAK/STAT signaling that leads to B cell proliferation and inhibition of macrophage activation.

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 2001, Anti-IL-10R1, (1B1.3a)

Product attributes

| | |
|---------------------------------------|---|
| Antibody number | #2001 |
| Antibody reactivity (target) | IL-10R1, |
| Antibody type | Primary |
| Host species | Rat |
| Clonality | Monoclonal |
| Clone | 1B1.3a |
| Isotype | IgG1, kappa |
| Molecular weight | 90-110 kDa |
| Synonyms | CD210 antigen; CDw210a; IBD28; IL10R alpha; IL-10R subunit 1; IL-10R1; IL10RA IL10RIL-10R subunit alpha; Interleukin 10 receptor; alpha; interleukin-10 receptor alpha chain; Interleukin-10 receptor subunit 1 |
| Entrez gene ID | 16154 |
| SwissProt | Q61727 |
| Unigene | 379327 |
| Immunogen | Purified recombinant extracellular region of Mouse IL10RA |
| Antibody target cellular localization | Plasma membrane |
| Species reactivity | Mouse |
| Expected antibody applications | Flow, surface (published for clone), Functional studies (published for clone) |
| Positive control | MPBMC's. Spleen, Thymus, Lymph Node, Placenta, Lung or Liver. |
| 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 |
| 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 |
| Shelf life | Guaranteed for at least 24 months from date of receipt when stored as recommended |
| 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 | Lymphocytes, Monocytes/macrophages, NK cells |

| 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.

1. J Immunol (2008) 181:5545-5550. (neutralization)
2. PLoS One (2017) 12(10): e0186317. (Flow, surface)

This datasheet was generated on July 16, 2025 at 11:34:16 AM. Visit product page to check for updated information before use.
Product link: <https://biotium.com/product/il-10r1-mouse-monoclonal-rat-antibody-1b1-3a/>