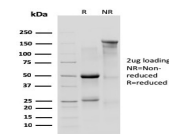


IL-10R1, Mouse Monoclonal Rat Antibody (1B1.3a)



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 (<i>Bos taurus</i>), 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.

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

This datasheet was generated on November 27, 2024 at 11:56:13 PM. Visit product page to check for updated information before use.

Product link: <https://biotium.com/product/il-10r1-mouse-monoclonal-rat-antibody-1b1-3a/>