## **HLA-DRB Monoclonal Mouse Antibody** (Bra30)

## **Product Description**

This MAb reacts with the HLA-DRB3 antigen, a member of MHC class II molecules. It does not cross react with HLA-DP and HLA-DQ. HLA-DR is a heterodimeric cell surface glycoprotein comprised of a 36 kDa alpha (heavy) chain and a 28 kDa beta (light) chain. It is expressed on B-cells, activated T-cells, monocytes/macrophages, dendritic cells and other non-professional APCs. In conjunction with the CD3/TCR complex and CD4 molecules, HLA-DR is critical for efficient peptide presentation to CD4 T cells. HLA-DR antigens also occur on a variety of epithelial cells and their corresponding neoplastic counterparts.

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 0407, Anti-HLA-DRB (Bra30)

## Product attributes

Call us: 800-304-5357

Product attributes			
Antibody number	#0407		
Antibody reactivity (target)	HLA-DRB		
Antibody type	Primary		
Host species	Mouse		
Clonality	Monoclonal		
Clone	Bra30		
Isotype	IgG2a, kappa		
Molecular weight	28 kDa (beta chain)		
Synonyms	HLA class II histocompatibility antigen DR beta 3 chain; HLADR3B; Major histocompatibility complex; class II DR beta 3		
Human gene symbol	HLA-DRB3		
Entrez gene ID	3123		
SwissProt	P01911		
Unigene	534322		
Immunogen	Human REH cells (non-T, non-B leukemia cell line)		
Antibody target cellular localization	Plasma membrane		
Species reactivity	Baboon, Chimpanzee, Dog, Human, Monkey, Squirrel		
Expected antibody applications	Flow, surface (published for clone), Functional studies (published for clone), WB (published for clone)		
Antibody application notes	Flow cytometry: 0.5-1 ug/million cells; Immunofluorescence: 0.5-1 ug/mL; Optimal dilution for a specific application should be determined., Higher concentration may be required for direct detection using primary antibody conjugates than for indirect detection with secondary antibody		
Positive control	Ramos, Daudi or HuT78 cells. Tonsil or lymph node		
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.		

Email: btinfo@biotium.com

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. Infection Immunity (2002) 70(5): 2492-2501. (functional studies)
- 2. Microbiol Immunol (2000) 44(12): 1011-1017. (Flow)
- 3. Transplant International (2011) 24(10), 1018-1026. (Flow, IF, WB)

This datasheet was generated on November 4, 2025 at 03:00:39 PM. Visit product page to check for updated information before use. Product link: <a href="https://biotium.com/product/hla-drb-monoclonal-mouse-antibody-bra30/">https://biotium.com/product/hla-drb-monoclonal-mouse-antibody-bra30/</a>