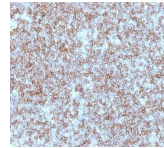


# MHC-II Monoclonal Mouse Antibody (Bra-14)



## Product Description

Reacts with a common epitope of human major histocompatibility (MHC) class II antigens, HLA-DR and DP. Human MHC class II antigens are transmembrane glycoproteins composed of an alpha chain (36 kDa) and a beta chain (27 kDa). They are expressed primarily on antigen presenting cells such as B lymphocytes, monocytes, macrophages, and thymic epithelial cells and are also present on activated T lymphocytes. Human MHC class II genes are located in the HLA-D region that encodes at least six and ten chain genes. Three loci, DR, DQ and DP, encode the major expressed products of the human class II region. The human MHC class II molecules bind intracellularly processed peptides and present them to T-helper cells. They, therefore, have a critical role in the initiation of the immune response. It has been shown that some autoimmune diseases are associated with certain class II alleles.

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 1129, Anti-MHC-II (Bra-14)**

## Product attributes

Antibody number	#1129
Antibody reactivity (target)	MHC-II
Antibody type	Primary
Host species	Mouse
Clonality	Monoclonal
Clone	Bra-14
Isotype	IgG3, kappa
Molecular weight	36 kDa (? chain) and 27 kDa (? chain)
Synonyms	Ia3; Ia-3; H-2Ea; H2-Ea; A1323765; I-Ealpha; E-alpha-f; MHC-H2-Ea
Human gene symbol	HLA-DPB1
Entrez gene ID	3115 (HLA-DP) & 3122 (HLA-DR)
SwissProt	P04440 (HLA-DP) & P01903 (HLA-DR)
Unigene	347270 (HLA-DP) & 520048 (HLA-DR)
Immunogen	Human REH cells
Antibody target cellular location	Plasma membrane
Verified antibody applications	IHC (FPPE) (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-fixed 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 minutes, Flow Cytometry 0.5-1 ug/million cells/0.1 mL, Optimal dilution for a specific application should be determined by user
Positive control	Raji 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
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	Immunology, MHC antigens

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, of LI-COR Bioscience.