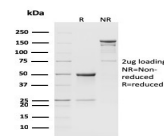


CD8b (Mouse) Monoclonal Rat Antibody (H35-17.2)



Product Description

This rat monoclonal antibody reacts with the mouse CD8 beta molecule. The CD8 beta chain associates with the CD8 alpha chain to form the CD8 alpha/beta heterodimer expressed on the surface of a majority of thymocytes, and on peripheral cytotoxic alpha beta TCR T cells. This monoclonal antibody identifies cytotoxic/suppressor T-cells that interact with MHC class I bearing targets. CD8 is thought to play a role in the process of T-cell mediated killing.

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 2003, Anti-CD8b, (H35-17.2)

Product attributes

Antibody number	#2003
Antibody reactivity (target)	CD8b,
Antibody type	Primary
Host species	Rat
Clonality	Monoclonal
Clone	H35-17.2
Isotype	IgG2b, kappa
Molecular weight	37 kDa
Synonyms	CD8b antigen; CD8 antigen beta polypeptide 1 (p37); CD8 beta; CD8B1; Leu2; Ly3; LYT3; P37; T lymphocyte surface glycoprotein beta chain; T-cell surface glycoprotein CD8 beta chain
Human gene symbol	Cd8b1 (Mouse)
Entrez gene ID	12526
SwissProt	P10300
Unigene	3331
Immunogen	Five day mixed leukocyte culture population, C57BL/6 anti-BALB/c
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 or Thymus.
Shipping condition	Room temperature
Storage Conditions	Note: store BSA-free antibodies at -10 to -35 °C, Store at 2 to 8 °C, Protect fluorescent conjugates from light
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	T-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. Infection Immunity (1994) 62(5): 1820-1829. (functional studies)
2. Meth Mol Biol vol 1559. Humana Press, New York, NY. https://doi.org/10.1007/978-1-4939-6786-5_2 (Flow, surface)
- 3.