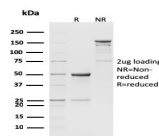


# CD8b (Mouse) Monoclonal Rat Antibody (H35-17.2)



## 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](mailto: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

<b>Antibody number</b>	#2003
<b>Antibody reactivity (target)</b>	CD8b,
<b>Antibody type</b>	Primary
<b>Host species</b>	Rat
<b>Clonality</b>	Monoclonal
<b>Clone</b>	H35-17.2
<b>Isotype</b>	IgG2b, kappa
<b>Molecular weight of antigen</b>	37 kDa
<b>Synonyms</b>	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
<b>Human gene symbol</b>	Cd8b1 (Mouse)
<b>Entrez gene ID</b>	12526
<b>SwissProt</b>	P10300
<b>UniGene</b>	3331
<b>Immunogen</b>	Five day mixed leukocyte culture population, C57BL/6 anti-BALB/c
<b>Antibody target cellular localization</b>	Plasma membrane
<b>Species reactivity</b>	Mouse
<b>Expected antibody applications</b>	Flow, surface (published for clone), Functional studies (published for clone)
<b>Positive control</b>	MPBMC's or Thymus.
<b>Shipping condition</b>	Room temperature
<b>Storage Conditions</b>	Note: store BSA-free antibodies at -10 to -35 °C, Store at 2 to 8 °C, Protect fluorescent conjugates from light
<b>Regulatory status</b>	For research use only (RUO)
<b>Antibody/conjugate formulation</b>	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
<b>Shelf life</b>	Guaranteed for at least 24 months from date of receipt when stored as recommended
<b>Product origin</b>	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.
<b>Cell/tissue expression</b>	T-cells

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>
BNC74	CF®740	742/767	633-685	775/50	<a href="#">CF®740 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, 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.

- Infection Immunity (1994) 62(5): 1820-1829. (functional studies)
- Meth Mol Biol vol 1559. Humana Press, New York, NY. [https://doi.org/10.1007/978-1-4939-6786-5\\_2](https://doi.org/10.1007/978-1-4939-6786-5_2) (Flow, surface)
-