CD8A Monoclonal Mouse Antibody (UCHT4)



Product Description

CD8 is a cell surface receptor expressed either as a heterodimer with the CD8 beta chain (CD8 alpha/beta) or as a homodimer (CD8 alpha/alpha). A majority of thymocytes and a subpopulation of mature T cells and NK cells express CD8a. CD8 binds to MHC class 1 and through its association with protein tyrosine kinase p56lck plays a role in T cell development and activation of mature T cells. For mature T-cells, CD4 and CD8 are mutually exclusive, so anti-CD8, generally used in conjunction with anti-CD4. It is a useful marker for distinguishing helper/inducer T-lymphocytes, and most peripheral T-cell lymphomas are CD4 /CD8-. Anaplastic large cell lymphoma is usually CD4 and CD8-, and in T-lymphoblastic lymphoma/leukemia, CD4 and CD8 are often co-expressed. CD8 is also found in littoral cell angioma of the spleen. This MAb binds to the CD8 alpha chain, at an epitope distinct from that recognized by UCHT14.

Also view our CD8a Monoclonal Mouse Biotium Choice Antibody (SK1). This antibody belongs to the Biotium Choice list of select antibodies that have been developed and validated in-house for optimal performance. The antibody has been validated in flow cytometry.

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 1628, Anti-CD8A (UCHT4)

Draduat attributes

Tumor expression

Call us: 800-304-5357

Product attributes				
Antibody number	#1628			
Antibody reactivity (target)	CD8A			
Antibody type	Primary			
Host species	Mouse			
Clonality	Monoclonal			
Clone	UCHT4			
Isotype	IgG2a, kappa			
Molecular weight	32 kDa			
Synonyms	CD8 antigen; alpha polypeptide (p32); T8/Leu-2 T-lymphocyte differentiation antigen; Ly3; LYT3; MAL; T-cell surface glycoprotein CD8 alpha chain			
Human gene symbol	CD8A			
Entrez gene ID	925			
SwissProt	P01732			
Unigene	85258			
Immunogen	Thymocytes followed by Sezary T cells.			
Antibody target cellular localization	Plasma membrane			
Verified antibody applications	Flow (surface) (verified)			
Species reactivity	Human			
Expected antibody applications	Functional studies (published for clone), IF (published for clone), IP (published for clone)			
Antibody application notes	Flow cytometry: 1-2 ug/million cells; 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				
FUSILIVE CUITIOI	Jurkat, HuT78 or hPBL. Tonsil.			
Shipping condition	Jurkat, HuT78 or hPBL. Tonsil. Room temperature			
Shipping condition	Room temperature Note: store BSA-free antibodies at -10 to -35°C, Store at 2 to 8			
Shipping condition Storage Conditions	Room temperature Note: store BSA-free antibodies at -10 to -35°C, Store at 2 to 8 °C, Protect fluorescent conjugates from light			
Shipping condition Storage Conditions Regulatory status Antibody/conjugate	Room temperature Note: store BSA-free antibodies at -10 to -35 °C, Store at 2 to 8 °C, Protect fluorescent conjugates from light For research use only (RUO) 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, Purified: 0.2 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			
Shipping condition Storage Conditions Regulatory status Antibody/conjugate formulation	Room temperature Note: store BSA-free antibodies at -10 to -35° C, Store at 2 to 8 °C, Protect fluorescent conjugates from light For research use only (RUO) 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 Guaranteed for at least 24 months from date of receipt when			

Leukemia/lymphoma

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,	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. Eur J Immunol (2005) 35: 2896-2908. (functional studies)
- 2. J Biol Chem (2007) 282(18): 13884-13894. (immunoprecipitation) 3. Traffic (2015) 16(12): 1330-1341. (immunofluorescence)