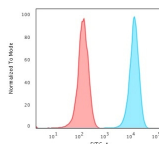


CD3e Monoclonal Mouse Antibody (UCHT1)



Product Description

Recognizes the epsilon-chain of CD3, which consists of five different polypeptide chains (designated as gamma, delta, epsilon, zeta, and eta) with MW ranging from 16-28 kDa. The CD3 complex is closely associated at the lymphocyte cell surface with the T cell antigen receptor (TCR). Reportedly, CD3 complex is involved in signal transduction to the T cell interior following antigen recognition. The CD3 antigen is first detectable in early thymocytes and probably represents one of the earliest signs of commitment to the T cell lineage. In cortical thymocytes, CD3 is predominantly intra-cytoplasmic. However, in medullary thymocytes, it appears on the T cell surface. CD3 antigen is a highly specific marker for T cells, and is present in majority of T cell neoplasms.

Also view our [CD3e Monoclonal Mouse Biotium Choice Antibody \(SK7\)](#). The 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 2473, Anti-CD3e (UCHT1)

Call us : [800-304-5357](tel:800-304-5357) Email: techsupport@biotium.com

Product attributes

Antibody number	#2473
Shelf life	Guaranteed for at least 24 months from date of receipt when stored as recommended
Antibody reactivity (target)	CD3e
Antibody type	Primary
Host species	Mouse
Clonality	Monoclonal
Clone	UCHT1
Isotype	IgG1, kappa
Molecular weight	20 kDa
Synonyms	CD3 epsilon; CD3 TCR complex; T cell antigen receptor complex epsilon subunit of T3; T-cell surface antigen T3/Leu-4 epsilon chain; T-cell surface glycoprotein CD3 epsilon chain; T3E; TCORE; TIT3 complex
Human gene symbol	CD3E
Entrez gene ID	916
SwissProt	P07766
Unigene	3003
Verified antibody applications	Flow (surface) (verified), IF (verified)
Immunogen	Human infant thymocytes and peripheral blood lymphocytes from a Sezary Syndrome donor
Expected antibody applications	ELISA (published for clone), Functional studies (published for clone), IP (published for clone), WB (published for clone)
Antibody target cellular localization	Plasma membrane
Species reactivity	Chimpanzee, Human
Positive control	Human peripheral blood. Jurkat cells. Tonsil or lymph node.
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
Cell/tissue expression	T-cells
Tumor expression	Leukemia/lymphoma
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.

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. Blood (1988) 71(3): 603-612. (immunofluorescence, western)
2. J Immunol Meth (1988) 106(1): 127-133. (ELISA)
3. J Exp Med (1991) 173: 7-17. (immunoprecipitation)
4. J Immunol (2002) 168: 3235-3241. (functional studies)

This datasheet was generated on August 25, 2025 at 06:47:14 AM. Visit product page to check for updated information before use.

Product link: <https://biotium.com/product/cd3e-monoclonal-mouse-antibody-ucht1/>