

CD15 / FUT4 / Lewis-X Recombinant Monoclonal Rabbit Antibody (FUT4/1478R)



Product Description

CD15 plays a role in mediating phagocytosis, bactericidal activity, and chemotaxis. It is present on >95% of granulocytes including neutrophils and eosinophils and to a lesser degree on monocytes. In addition, CD15 is expressed in Reed-Sternberg cells and some epithelial cells. CD15 antibody is very useful in the identification of Hodgkin's disease. CD15 is occasionally expressed in large cell lymphomas of both B and T phenotypes which otherwise have a guite distinct histological appearance.

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 1478, Anti-CD15|FUT4|Lewis-x (FUT4/1478R)

Product attributes

Call us: 800-304-5357

Product attributes				
Antibody number	#1478			
Antibody reactivity (target)	CD15, FUT4, Lewis-x			
Antibody type	Primary			
Host species	Rabbit			
Clonality	Recombinant Monoclonal			
Clone	FUT4/1478R			
Isotype	IgG, kappa			
Molecular weight	220 kDa			
Synonyms	CD162; CLA; Cutaneous lymphocyte associated antigen; P-selectin glycoprotein ligand 1; PSGL1; Selectin P ligand; SELPLG			
Human gene symbol	FUT4			
Entrez gene ID	2526			
SwissProt	P22083			
Unigene	654379			
Immunogen	Purified human neutrophils			
Verified antibody applications	Flow (verified), IF (verified), IHC (FFPE) (verified)			
Antibody target cellular localization	Plasma membrane, Perinuclear			
	Human, Mouse, Rat			
Species reactivity	Human, Mouse, Rat			
Species reactivity Antibody application notes	Human, Mouse, Rat Staining of formalin-fixed tissues requires boiling tissue sections in 10 mM Tris, 1 mM EDTA pH 9.0 for 10-20 min followed by cooling at RT for 20 min, Flow cytometry 2-4 ug/million cells, Immunofluorescence 2-4 ug/mL, Immunohistology (formalin) 5-10 ug/mL, Optimal dilution for a specific application should be determined by user			
•	Staining of formalin-fixed tissues requires boiling tissue sections in 10 mM Tris, 1 mM EDTA pH 9.0 for 10-20 min followed by cooling at RT for 20 min, Flow cytometry 2-4 ug/million cells, Immunofluorescence 2-4 ug/mL, Immunohistology (formalin) 5-10 ug/mL, Optimal dilution for a specific application should be			
Antibody application notes	Staining of formalin-fixed tissues requires boiling tissue sections in 10 mM Tris, 1 mM EDTA pH 9.0 for 10-20 min followed by cooling at RT for 20 min, Flow cytometry 2-4 ug/mlliminon cells, Immunofluorescence 2-4 ug/ml, Immunofistology (formalin) 5-10 ug/ml, Optimal dilution for a specific application should be determined by user			
Antibody application notes Positive control	Staining of formalin-fixed tissues requires boiling tissue sections in 10 mM Tris, 1 mM EDTA pH 9.0 for 10-20 min followed by cooling at RT for 20 min, Flow cytometry 2-4 ug/million cells, Immunofluorescence 2-4 ug/mL, Immunohistology (formalin) 5-10 ug/mL, Optimal dilution for a specific application should be determined by user U937 cells, Reed-Sternberg's cells in Hodgkin's lymphoma			
Antibody application notes Positive control Shipping condition	Staining of formalin-fixed tissues requires boiling tissue sections in 10 mM Tris, 1 mM EDTA pH 9.0 for 10-20 min followed by cooling at RT for 20 min, Flow cytometry 2-4 ug/million cells, Immunofluorescence 2-4 ug/mL, Immunohistology (formalin) 5-10 ug/mL, Optimal dilution for a specific application should be determined by user U937 cells, Reed-Sternberg's cells in Hodgkin's lymphoma Room temperature Store at 2 to 8 °C, Protect fluorescent conjugates from light,			
Antibody application notes Positive control Shipping condition Storage Conditions	Staining of formalin-fixed tissues requires boiling tissue sections in 10 mM Tris, 1 mM EDTA pH 9.0 for 10-20 min followed by cooling at RT for 20 min, Flow cytometry 2-4 ug/mllion cells, immunofluorescence 2-4 ug/ml., Immunofluorescence 2-4 ug/ml., Immunofluorescence 2-4 ug/ml., Optimal dilution for a specific application should be determined by user U937 cells, Reed-Sternberg's cells in Hodgkin's lymphoma Room temperature Store at 2 to 8 °C, Protect fluorescent conjugates from light, Note: store BSA-free antibodies at -10 to -35 °C Guaranteed for at least 24 months from date of receipt when			
Antibody application notes Positive control Shipping condition Storage Conditions Shelf life	Staining of formalin-fixed tissues requires boiling tissue sections in 10 mM Tris, 1 mM EDTA pH 9.0 for 10-20 min followed by cooling at RT for 20 min, Flow cytometry 2-4 ug/million cells, immunofluorescence 2-4 ug/mL, Immunofluorescence 2-4 ug/mL, Immunofluorescence 2-4 ug/mL, optimal dilution for a specific application should be determined by user U937 cells, Reed-Sternberg's cells in Hodgkin's lymphoma Room temperature Store at 2 to 8 °C, Protect fluorescent conjugates from light, Note: store BSA-free antibodies at -10 to -35 °C Guaranteed for at least 24 months from date of receipt when stored as recommended			
Antibody application notes Positive control Shipping condition Storage Conditions Shelf life Regulatory status Antibody/conjugate	Staining of formalin-fixed tissues requires boiling tissue sections in 10 mM Tris, 1 mM EDTA pH 9.0 for 10-20 min followed by cooling at RT for 20 min, Flow cytometry 2-4 ugimillion cells, Immunofluorescence 2-4 ug/mL, Immunofluorescence 2-4 ug/mL, Immunofluorescence 2-4 ug/mL, Immunofluorescence 2-4 ug/mL, Optimal dilution for a specific application should be determined by user U937 cells, Reed-Sternberg's cells in Hodgkin's lymphoma Room temperature Store at 2 to 8 °C, Protect fluorescent conjugates from light, Note: store BSA-free antibodies at -10 to -35 °C Guaranteed for at least 24 months from date of receipt when stored as recommended 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; BSA-free: 1 mg/mL in PBS/0.05% BSA, Purified; BSA-free: 1 mg/mL in			
Antibody application notes Positive control Shipping condition Storage Conditions Shelf life Regulatory status Antibody/conjugate formulation	Staining of formalin-fixed tissues requires boiling tissue sections in 10 mM Tris, 1 mM EDTA pH 9.0 for 10-20 min followed by cooling at RT for 20 min, Flow cytometry 2-4 ug/mll.lino cells, Immunofluorescence 2-4 ug/ml., Optimal dilution for a specific application should be determined by user U937 cells, Reed-Sternberg's cells in Hodgkin's lymphoma Room temperature Store at 2 to 8 °C, Protect fluorescent conjugates from light, Note: store BSA-free antibodies at -10 to -35 °C Guaranteed for at least 24 months from date of receipt when stored as recommended For research use only (RUO) Conjugates: 0.1 mg/mL in PBS/0.05% BSA/0.05% azide, HRP conjugates: 0.1 mg/mL in PBS/0.05% BSA/10.05% azide, Purified: 0.2 mg/mL in PBS/0.05% BSA/10.05% azide, Purified. BSA-free: 1 mg/mL in PBS without azide			
Antibody application notes Positive control Shipping condition Storage Conditions Shelf life Regulatory status Antibody/conjugate formulation Antibody research areas	Staining of formalin-fixed tissues requires boiling tissue sections in 10 mM Tris, 1 mM EDTA pH 9.0 for 10-20 min followed by cooling at RT for 20 min, Flow cytometry 2-4 ug/million cells, Immunofluorescence 2-4 ug/mL, Immunohistology (formalin) 5-10 ug/mL, Optimal dilution for a specific application should be determined by user U937 cells, Reed-Sternberg's cells in Hodgkin's lymphoma Room temperature Store at 2 to 8 °C, Protect fluorescent conjugates from light, Note: store BSA-free antibodies at -10 to -35 °C Guaranteed for at least 24 months from date of receipt when stored as recommended 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 Immunology Product may contain either bovine serum albumin (BSA) from bovine serum (Bos taurus), or recombinant BSA produced in			

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.

This datasheet was generated on November 30, 2025 at 01:39:19 AM. Visit product page to check for updated information before use. Product link: https://biotium.com/product/recombinant-rabbit-monoclonal-anti-cd15-fut4-fut41478r/