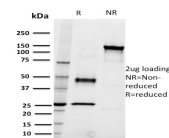


CD22 / BL-CAM Monoclonal Mouse Antibody (RFB4)



Product Description

Recognizes a protein of 130-140 kDa, identified as CD22 (also known as BL-CAM). CD22 expression is restricted to normal and neoplastic B cells and is absent from other haemopoietic cell types. In B-cell ontogeny, CD22 is first expressed in the cytoplasm of pro-B and pre-B cells, and on the surface as B cells mature to become IgD. It is not expressed by plasma cells. CD22 is found highly expressed in follicular mantle and marginal zone B-cells, and while germinal center B-cells are relatively weak. CD22 is a member of the immunoglobulin superfamily and serves as an adhesion receptor for sialic acid-bearing ligands expressed on erythrocytes and all leukocyte classes. It also associates with tyrosine kinases and play a role in signal transduction and B-cell activation.

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 1594, Anti-BL-CAM|CD22 (RFB4)

Call us : [800-304-5357](tel:800-304-5357)

Product attributes

Antibody number	#1594
Antibody reactivity (target)	BL-CAM, CD22
Antibody type	Primary
Host species	Mouse
Clonality	Monoclonal
Clone	RFB4
Isotype	IgG1, kappa
Molecular weight of antigen	130-140 kDa
Synonyms	B-lymphocyte cell adhesion molecule (BL-CAM); B-cell receptor CD22; CD22; Lectin 2; Lyb8; Sialic acid-binding Ig-like lectin 2 (Siglec-2); SIGLEC2; T-cell surface antigen Leu-14
Human gene symbol	CD22
Entrez gene ID	933
SwissProt	P20273
Unigene	579691
Immunogen	Human tonsil lymphocytes.
Verified antibody applications	Flow (verified)
Antibody target cellular localization	Plasma membrane
Species reactivity	Human
Expected antibody applications	Flow, surface (published for clone), Functional studies (published for clone)
Antibody application notes	ELISA: For coating, order antibody without BSA; Flow cytometry: 0.5-1 ug/million cells; Immunofluorescence: 0.5-1 ug/mL; 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	Ramos, Raji or Daudi cells. Human Tonsil, Lymph Node or Spleen.
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 (<i>Bos taurus</i>), or recombinant BSA produced in Chinese hamster ovary cells. Inquire for the specific lot.
Cell/tissue expression	B-cells, Lymphocytes

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.

- Clin Cancer Res (2005) 11(21): 7920-7928. (Flow, surface; functional studies)
- Biomolecules (2020) 10: 1140. (oncotoxins)