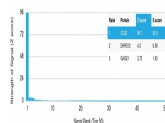


CD22 / BL-CAM Monoclonal Mouse Antibody (BLCAM/1796)



Product Description

This antibody 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 1796, Anti-CD22 / BL-CAM (BLCAM/1796)**

Product attributes

Antibody number	#1796
Antibody reactivity (target)	BL-CAM, CD22
Antibody type	Primary
Host species	Mouse
Clonality	Monoclonal
Clone	BLCAM/1796
Isotype	IgG2b, kappa
Molecular weight	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	Recombinant fragment of human CD22 protein (around aa 52-178) (exact sequence is proprietary)
Verified antibody applications	Flow (verified), IF (verified), WB (verified)
Antibody target cellular localization	Plasma membrane
Species reactivity	Human
Positive control	Raji or Ramos cells (IF/FACS). Raji cell lysate (WB). Human tonsil, lymph node or spleen (IHC).
Shipping condition	Room temperature
Storage Conditions	Store at 2 to 8 °C. Protect fluorescent conjugates from light. Note: store BSA-free antibodies at -10 to -35 °C
Shelf life	Guaranteed for at least 24 months from date of receipt when stored as recommended
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
Validated in protein array	Monospecific
Antibody research areas	Immunology
Cell/tissue expression	B-cells, Lymphocytes

Antibody # prefix	Conjugation	Ex/Em (nm)	Laser line	Detection channel
BNC04	CF [®] 405S	404/431	405	DAPI (microscopy), AF405
BNC88	CF [®] 488A	490/515	488	GFP, FITC
BNC68	CF [®] 568	562/583	532, 561	RFP, TRITC
BNC94	CF [®] 594	593/614	561	Texas Red [®]
BNC40	CF [®] 640R	642/662	633-640	Cy [®] 5
BNC47	CF [®] 647	650/665	633-640	Cy [®] 5
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

Dye Features
[CF[®]405S Features](#)
[CF[®]488A Features](#)
[CF[®]568 Features](#)
[CF[®]594 Features](#)
[CF[®]640R Features](#)
[CF[®]647 Features](#)

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 LI-COR Bioscience.