## CD44v9 Recombinant Monoclonal Rabbit Antibody (CD44v9/2344R)



## **Product Description**

This antibody recognizes an epitope encoded by exon v4 on the variant portion of human CD44. The CD44 molecule belongs to a family of cellular adhesion molecules found on a wide range of normal and malignant cells in epithelial, mesothelial and hemopoiesis tissues. CD44 is a single gene with 20 exons, of which 10 are normally expressed to encode the basic CD44 (H-CAM) molecule. The additional 10 exons (v1 to v10) are only expressed by alternative splicing of the nuclear RNA. The expression of specific cell adhesion molecule CD44 splice variants has been reported to be associated with metastasis in certain human malignancies.

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 <a href="mailto:order@biotium.com">order@biotium.com</a> to inquire about stock status and lead times before placing your order.

Catalog number key for antibody number 2344, Anti-CD44v9 (CD44v9/2344R)

## Duadinat attailanta

Antibody research areas

Product origin

Antibody number #2344  Antibody reactivity (target) CD44v9  Antibody type Primary  Host species Rabbit  Clonality Recombinant Monoclonal  Clone CD44v9/2344R  Isotype IgG  Molecular weight 80-95 kDa  Synonyms LHR; BA-1; chondroitin sulfate proteoglycan 8 (CSPG8); Extracellular Matrix Receptor III (ECM III); GP90 Lymph Homing Adhesion Receptor; HOAH; HCELL; hematopoid E- and L-selectin ligand; Heparan Sulfate Proteoglycan; Antigen; HAS; HUTCH I; Hyaluronate Receptor; Indian b group; Inlu Related p80 Glycoprotein; Ly 24; MDU2; MDI MIC4; MUTCH I; Phagocytic Glycoprotein 1 (PGP-1)  Human gene symbol CD44	
Antibody type Primary  Host species Rabbit  Clonality Recombinant Monoclonal  Clone CD44v9/2344R  Isotype IgG  Molecular weight 80-95 kDa  Synonyms  LHR; BA-1; chondroitin sulfate proteoglycan 8 (CSPG8); Extracellular Matrix Receptor III (ECM III); GP90 Lymphc Homing Adhesion Receptor; HCAM; HCELL; hematopolic E- and L-selectin ligand; Heparan Sulfate Proteoglycan; Antigen; HAS; HUTCH I; Hyaluronate Receptor; Indian b group; Inlu Related p80 Glycoprotein; Ly 24; MDU2; MDI MIC4; MUTCH I; Phagocytic Glycoprotein 1 (PGP-1)	
Host species Rabbit  Clonality Recombinant Monoclonal  Clone CD44v9/2344R  Isotype IgG  Molecular weight 80-95 kDa  Synonyms  LHR; BA-1; chondroitin sulfate proteoglycan 8 (CSPG8); Extracellular Matrix Receptor III (ECM III); GP90 Lymphc Homing Adhesion Receptor; HCAM; HCELL; hematopolic E- and L-selectin ligand; Heparan Sulfate Proteoglycan; Antigen; HAS; HUTCH I; Hyaluronate Receptor; Indian b group; Inlu Related p80 Glycoprotein; Ly 24; MDU2; MDI MIC4; MUTCH I; Phagocytic Glycoprotein 1 (PGP-1)	
Clone CD44v9/2344R  Isotype IgG  Molecular weight 80-95 kDa  Synonyms LHR; BA-1; chondroitin sulfate proteoglycan 8 (CSPG8); Extracellular Matrix Receptor III (ECM III); GP90 Lymphc Homing Adhesion Receptor; HCAM; HCELL; hematopoid E- and L-selectin ligand; Heparan Sulfate Proteoglycan; Antigen; HAS; HUTCH I; Hyaluronate Receptor; Indian b group; Inlu Related p80 Glycoprotein; Ly 24; MDU2; MDI MIC4; MUTCH I; Phagocytic Glycoprotein 1 (PGP-1)	
Clone CD44v9/2344R  Isotype IgG  Molecular weight 80-95 kDa  Synonyms LHR; BA-1; chondroitin sulfate proteoglycan 8 (CSPG8); Extracellular Matrix Receptor III (ECM III); GP90 Lymphc Homing Adhesion Receptor; HCAM; HCELL; hematopoid E- and L-selectin ligand; Heparan Sulfate Proteoglycan; Antigen; HAS; HUTCH I; Hyaluronate Receptor; Indian b group; Inlu Related p80 Glycoprotein; Ly 24; MDU2; MDI MIC4; MUTCH I; Phagocytic Glycoprotein 1 (PGP-1)	
Isotype IgG  Molecular weight 80-95 kDa  Synonyms  LHR; BA-1; chondroitin sulfate proteoglycan 8 (CSPG8); Extracellular Matrix Receptor III (ECM III); GP90 Lymphc Homing Adhesion Receptor; HCAM; HCELL; hematopoid E- and L-selectin ligand; Heparan Sulfate Proteoglycan; Antigen; HAS; HUTCH I; Hyaluronate Receptor; Indian b group; Inlu Related p80 Glycoprotein; Ly 24; MDU2; MDI MIC4; MUTCH I; Phagocytic Glycoprotein 1 (PGP-1)	
Molecular weight  80-95 kDa  LHR; BA-1; chondroitin sulfate proteoglycan 8 (CSPG8); Extracellular Matrix Receptor III (ECM III); GP90 Lymphc Homing Adhesion Receptor; HCAM; HCELL; hematopoio E- and L-selectin ligand; Heparan Sulfate Proteoglycan; Antigen; HAS; HUTCH I; Hyaluronate Receptor; Indian b group; Inlu Related p80 Glycoprotein; Ly 24; MDU2; MDI MIC4; MUTCH I; Phagocytic Glycoprotein 1 (PGP-1)	
Synonyms  LHR; BA-1; chondroitin sulfate proteoglycan 8 (CSPG8); Extracellular Matrix Receptor III (ECM III); GP90 Lympho Homing Adhesion Receptor; HC3M; HCELL; hematopoi E- and L-selectin ligand; Heparan Sulfate Proteoglycan; Antigen; HAS; HUTCH I; Hyaluronate Receptor; Indian b group; Inlu Related p80 Glycoprotein; Ly 24; MDU2; MDI MIC4; MUTCH I; Phagocytic Glycoprotein 1 (PGP-1)	
Extracellular Matrix Receptor III (ECM III); GP90 Lymph; Homing Adhesion Receptor; HCAM; HCELL; hematopoi E- and L-selectin ligand; Heparan Sulfate Proteoglycan; Antigen; HAS; HUTCH I; Hyaluronate Receptor; Indian b group; Inlu Related p80 Glycoprotein; Ly 24; MDU2; MDI MIC4; MUTCH I; Phagocytic Glycoprotein 1 (PGP-1)	
Human gene symbol CD44	ocyte etic cell Hermes blood
- •	
Entrez gene ID 960	
SwissProt P16070	
<b>Unigene</b> 502328	
Immunogen Recombinant fragment corresponding to the v9 domain of human CD44	of
Antibody target cellular Plasma membrane localization	
Verified antibody applications IHC (FFPE) (verified)	
Species reactivity Human	
Antibody application notes  Higher concentration may be required for direct detection primary antibody conjugates than for indirect detection we secondary antibody. Immunohistology (formalin): 1-2 ug/30 minutes at RT, Stalning of formalin-fixed tissues requibolling tissue sections in 10 mM citrate buffer, pH 6.0, forminutes followed by cooling at RT for 20 minutes, Optima dilution for a specific application should be determined by	vith /mL for ires r 10-20 al
Positive control Cervix or Tongue Squamous Cell Carcinoma	
Shipping condition Room temperature	
Storage Conditions Store at 2 to 8 °C, Protect fluorescent conjugates from lie Note: store BSA-free antibodies at -10 to -35 °C	ght,
Shelf life Guaranteed for at least 24 months from date of receipt w stored as recommended	hen
Regulatory status For research use only (RUO)	
Antibody/conjugate formulation Conjugates: 0.1 mg/mL in PBS/0.1% BSA/0.05% azide, conjugates: 0.1 mg/mL in PBS/0.05% BSA, Purified: 0.2 in PBS/0.05% BSA/0.05% BSA/0.0	LIDD

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.

Call us: 800-304-5357 Email: techsupport@biotium.com

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.

This datasheet was generated on September 17, 2025 at 07:24:33 PM. Visit product page to check for updated information before use. Product link: <a href="https://biotium.com/product/cd44v9-recombinant-monoclonal-rabbit-antibody-cd44v9-2344r/">https://biotium.com/product/cd44v9-recombinant-monoclonal-rabbit-antibody-cd44v9-2344r/</a>