

MECA-79 Monoclonal Rat Antibody (MECA-79)

Product Description

MECA-79 recognizes a carbohydrate epitope shared with a group of sulfation decorated sialomucins, including sulfated ligands for CD62L (CD34, GlyCAM-1, Sgp200, and a subset of MAdCAM-1). This set of antigens has been referred to as peripheral node addressin (PNAd) with the molecular mass 50-250 kD. It has been identified that GlcNAc-6-SO4 sulfation contributes to MECA-79 binding and the core 1 beta1, 3-N-acetylglucosaminyltransferase is required for the generation of the MECA-79 epitope. MECA-79 is expressed on high endothelial venules (HEV) of lymphoid tissues, chronic inflamed tissues and rheumatoid synovia. The interaction of PNAd with CD62L receptor is involved in tethering and rolling of lymphocytes along HEV in lymphoid tissues. MECA-79 antibody reacts with PNAd from mouse, human and many other species and blocks L-selectin-dependent lymphocyte adhesion in vitro and in vivo.

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 1599, Anti-MECA-79 (MECA-79)

Product attributes

Antibody number	#1599
Antibody reactivity (target)	MECA-79
Antibody type	Primary
Host species	Rat
Clonality	Monoclonal
Clone	MECA-79
Isotype	IgM, kappa
Molecular weight of antigen	90-110 kDa
Synonyms	MECA-79; Peripheral Node Addressin (PNAd)
Entrez gene ID	Not Known
SwissProt	Not Known
Unigene	Not Known
Immunogen	Mouse lymph node stromal cells
Antibody target cellular localization	Plasma membrane
Species reactivity	Human, Mouse, Rat
Expected antibody applications	IHC (frozen) (published for clone), ELISA (published for clone), Flow, surface (published for clone), Functional studies (published for clone), IHC (FFPE) (published for clone), IF (published for clone)
Positive control	Human or mouse tonsil.
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.

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. J Cell Biol (1988) 107: 1853-1862. (IF; IHC, frozen; functional studies)
2. J Pathol (2005) 206: 260-268. (IHC, frozen)
3. Immunology (2006) 119: 461-469. (Flow; IHC, frozen)
4. Int J Clin Exp Pathol (2014) 7(5): 1967-1976. (IHC, FFPE)

This datasheet was generated on June 17, 2026 at 03:07:15 AM. Visit product page to check for updated information before use.

Product link: <https://biotium.com/product/meca-79-monoclonal-rat-antibody-meca-79/>