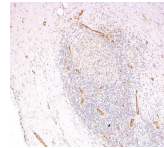


von Willebrand Factor / vWF Monoclonal Mouse Antibody (3E2D10 + VWF635)



Product Description

von Willebrand Factor (vWF) is a multimeric glycoprotein that is found in endothelial cells, plasma and platelets. It acts as a carrier protein for Factor VIII and promotes platelet adhesion and aggregation. vWF undergoes a variety of posttranslational modifications that influence the affinity and availability for Factor VIII, including cleavage of the propeptide and formation of N-terminal disulfide bonds. This antibody helps to establish the endothelial nature of some lesions of disputed histogenesis, e.g. Kaposi's sarcoma and cardiac myxoma. It is widely used for differentiating vascular lesions from those of other tissue differentiation within a panel of other vascular markers although not all tumors of endothelial differentiation contain this antigen.

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 0646, Anti-von Willebrand Factor (3E2D10 VWF635)

Product attributes

Antibody number	#0646
Antibody reactivity (target)	von Willebrand Factor, vWF
Antibody type	Primary
Host species	Mouse
Clonality	Monoclonal
Clone	3E2D10 + VWF635
Isotype	IgG's
Molecular weight	250 kDa
Synonyms	Coagulation Factor VIII, Factor VIII Related Antigen, F8VWF, von Willebrand Antigen 2, von Willebrand Disease (vWD)
Human gene symbol	VWF
Entrez gene ID	7450
SwissProt	P04275
Unigene	440848
Immunogen	Recombinant human vWF fragment
Antibody target cellular	Secreted (extracellular), Vesicular
Verified antibody applications	IHC (FFPE) (verified)
Species reactivity	Human
Antibody application notes	Higher concentration may be required for direct detection using primary antibody conjugates than for indirect detection with secondary antibody, Immunoprecipitation: 0.5-1 ug/500 ug protein lysate, Immunofluorescence: 0.5-1 ug/mL, Immunohistology (formalin), Staining of formalin-fixed tissues requires boiling tissue sections in 10 mM citrate buffer, pH 6.0, for 10-20 min followed by cooling at RT for 20 minutes, Flow Cytometry 0.5-1 ug/million cells/0.1 mL, Western blotting 0.5-1.0 ug/mL, Optimal dilution for a specific application should be determined by user
Positive control	HUVEC cells or Tonsil
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
Antibody research areas	Hematology
Cell/tissue expression	Endothelial cells

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
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, or LI-COR Bioscience.