

VEGFR2 / Flk-1 / KDR3 / CD309 Monoclonal Rat Antibody (DC101)



Product Description

This MAb is specific to Mouse VEGFR2/FLK-1/CD309 and does not cross-react with FLK-2. VEGFR2 is a type I transmembrane glycoprotein. It is a member of the CSF-1/PDGF receptor family of type III tyrosine kinase receptors. Endothelial cells, embryonic tissues, and megakaryocytes mainly express VEGFR2. It plays an important role in the regulation of angiogenesis, vasculogenesis, and vascular permeability. The ligands of VEGFR2 include VEGF-A, VEGF-C, VEGF-D, and VEGF splice isoforms. Ligation of VEGFR2 with its ligands results in the receptor dimerization and auto-phosphorylation, stimulating endothelial cell proliferation and migration.

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 0531, Anti-CD309|Flk-1|KDR3|VEGF-R2 (DC101)

Product attributes

Call us: 800-304-5357

Product attributes				
Antibody number	#0531			
Antibody reactivity (target)	CD309, Flk-1, KDR3, VEGF-R2			
Antibody type	Primary			
Host species	Rat			
Clonality	Monoclonal			
Clone	DC101			
Isotype	IgG1, kappa			
Molecular weight	180-190 kDa			
Synonyms	CD309; Fetal liver kinase 1 (FLK-1); KDR; Kinase insert domain receptor (a type III receptor tyrosine kinase); KRD1; Ly73; Protein tyrosine kinase receptor FLK1; Vascular endothelial growth factor receptor 2			
Human gene symbol	Kdr (Mouse)			
Entrez gene ID	16542			
SwissProt	35918			
Unigene	285			
Immunogen	Recombinant full-length Mouse VEGFR2 protein			
Antibody target cellular localization	Plasma membrane			
Species reactivity	Mouse			
Expected antibody applications	IHC (frozen) (published for clone), Flow, surface (published for clone), Functional studies (published for clone)			
Positive control	Mouse epidermoid, renal cell carcinoma, pancreatic or glioblastoma cells.			
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 (Bos taurus), or recombinant BSA produced in Chinese hamster ovary cells. Inquire for the specific lot.			
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
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. Cancer Metastasis Rev (1998) 17(2):155-61. (functional studies)
- 2. J Exp Med (2001) 193(9): 1005-1014. (Flow)
- 3. FASEB J (2005) 19(14): 2005-2007. (IHC, frozen)

This datasheet was generated on November 30, 2025 at 01:24:58 AM. Visit product page to check for updated information before use. Product link: https://biotium.com/product/vegfr2-filk-1-kdr3-cd309-monoclonal-rat-antibody-dc101/