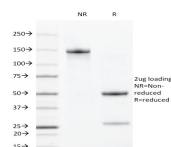


# VEGFR2 / Flk-1 / KDR3 / CD309 Monoclonal Rat Antibody (DC101)



## Product Description

This MAbs 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](mailto: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

<b>Antibody number</b>	#0531
<b>Antibody reactivity (target)</b>	CD309, Flk-1, KDR3, VEGF-R2
<b>Antibody type</b>	Primary
<b>Host species</b>	Rat
<b>Clonality</b>	Monoclonal
<b>Clone</b>	DC101
<b>Isotype</b>	IgG1, kappa
<b>Molecular weight</b>	180-190 kDa
<b>Synonyms</b>	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
<b>Human gene symbol</b>	Kdr (Mouse)
<b>Entrez gene ID</b>	16542
<b>SwissProt</b>	35918
<b>Unigene</b>	285
<b>Immunogen</b>	Recombinant full-length Mouse VEGFR2 protein
<b>Antibody target cellular localization</b>	Plasma membrane
<b>Species reactivity</b>	Mouse
<b>Expected antibody applications</b>	IHC (frozen) (published for clone), Flow, surface (published for clone), Functional studies (published for clone)
<b>Positive control</b>	Mouse epidermoid, renal cell carcinoma, pancreatic or glioblastoma cells.
<b>Shipping condition</b>	Room temperature
<b>Storage Conditions</b>	Note: store BSA-free antibodies at -10 to -35 °C, Store at 2 to 8 °C, Protect fluorescent conjugates from light
<b>Regulatory status</b>	For research use only (RUO)
<b>Antibody/conjugate formulation</b>	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
<b>Shelf life</b>	Guaranteed for at least 24 months from date of receipt when stored as recommended
<b>Product origin</b>	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.
<b>Cell/tissue expression</b>	Endothelial cells

Antibody # prefix	Conjugation	Ex/Em (nm)	Laser line	Detection channel	Dye Features
BNC04	CF®405S	404/431	405	DAPI (microscopy), AF405	<a href="#">CF®405S Features</a>
BNC88	CF®488A	490/515	488	GFP, FITC	<a href="#">CF®488A Features</a>
BNC68	CF®568	562/583	532, 561	RFP, TRITC	<a href="#">CF®568 Features</a>
BNC94	CF®594	593/614	561	Texas Red®	<a href="#">CF®594 Features</a>
BNC40	CF®640R	642/662	633-640	Cy®5	<a href="#">CF®640R Features</a>
BNC47	CF®647	650/665	633-640	Cy®5	<a href="#">CF®647 Features</a>
BNC74	CF®740	742/767	633-685	775/50	<a href="#">CF®740 Features</a>
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 March 16, 2025 at 12:49:55 PM. Visit product page to check for updated information before use. Product link: <https://biotium.com/product/vegfr2-flk-1-kdr3-cd309-monoclonal-rat-antibody-dc101/>