

PRODUCT INFORMATION



VEGF-A 164 variant (mouse, recombinant)

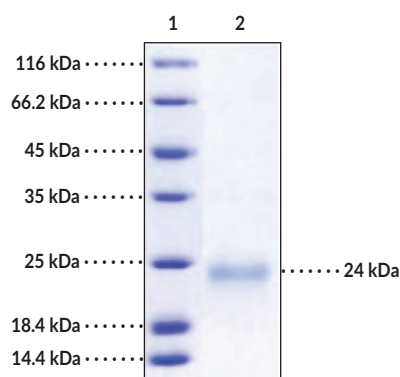
Item No. 32068

Overview and Properties

Synonym:	Vascular Endothelial Growth Factor A 164 variant
Source:	Active recombinant mouse VEGF-A 164 variant expressed in insect cells (baculovirus)
Amino Acids:	27-190
Uniprot No.:	Q00731-2
Molecular Weight:	19.4 kDa
Storage:	-80°C (as supplied)
Stability:	≥1 year
Purity:	≥90% estimated by SDS-PAGE
Supplied in:	Lyophilized from sterile 20 mM HAC-sodium acetate, pH 6.5, with 500 mM sodium chloride
Endotoxin Testing:	<1.0 EU/μg, determined by the LAL endotoxin assay
Bioactivity:	See figures for details

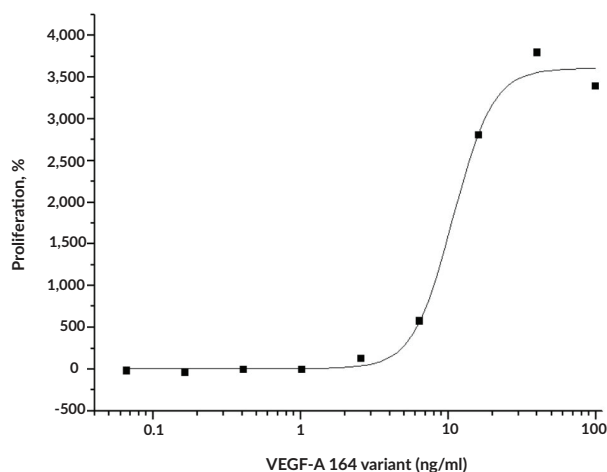
Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Images



Lane 1: MW Markers
Lane 2: VEGF-A 164 variant

SDS-PAGE Analysis of VEGF-A 164 variant.
This protein has a calculated molecular weight of 19.4 kDa. By SDS-PAGE, under reducing conditions, the molecular weight of the protein is 24 kDa due to apparent post-translational modifications.



Cell proliferation assay using human umbilical vein endothelial cells (HUVECs). The ED₅₀ value for this effect is typically 5-22 ng/ml.

WARNING
THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

SAFETY DATA
This material should be considered hazardous until further information becomes available. Do not ingest, inhale, get in eyes, on skin, or on clothing. Wash thoroughly after handling. Before use, the user must review the [complete](#) Safety Data Sheet, which has been sent via email to your institution.

WARRANTY AND LIMITATION OF REMEDY
Buyer agrees to purchase the material subject to Cayman's Terms and Conditions. Complete Terms and Conditions including Warranty and Limitation of Liability information can be found on our website.

Copyright Cayman Chemical Company, 03/09/2021

CAYMAN CHEMICAL
1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA
PHONE: [800] 364-9897
[734] 971-3335
FAX: [734] 971-3640
CUSTSERV@CAYMANCHEM.COM
WWW.CAYMANCHEM.COM

PRODUCT INFORMATION



Description

VEGF-A is a member of the PDGF/VEGF family of growth factors that regulates development, proliferation, and maintenance of the vascular system.¹ Alternative splicing of *Vegfa* pre-mRNA leads to the formation of VEGF-A 164 variant (VEGF-164), an isoform of VEGF-A that contains two heparin binding domains that interact with VEGFR1 and VEGFR2, the VEGFR co-receptors neuropilin-1 (NRP1) and NRP2, and matrix-associated glycosaminoglycans.^{2,3} VEGF-164 exists as a secreted or matrix-bound homodimer and is expressed in the brain, bone, eye, heart, lung, kidney and liver.^{4,5} Transgenic overexpression of VEGF-164 in the podocytes of mice induces complete collapse of the glomerular capillary network and is perinatal lethal.⁶ Inhibition of VEGF-164 with a neutralizing aptamer inhibits leukocyte adhesion and neovascularization in the retina in a rat model of hyperoxia-induced proliferative retinopathy.⁷ Serum levels of the human homolog VEGF-165 are increased in patients with melanoma and positively correlated with metastasis.⁸ VEGF-165 levels are also increased in the serum, synovial fluid, and inflamed joints of patients with rheumatoid arthritis.⁹ Cayman's VEGF-A 164 variant (mouse, recombinant) protein can be used for cell-based assay applications. This protein consists of 164 amino acids, has a calculated molecular weight of 19.4 kDa, and a predicted N-terminus of Ala27 after signal peptide cleavage. By SDS-PAGE, under reducing conditions, the molecular weight of the protein is 24 kDa due to apparent post-translational modifications.

References

1. Shibuya, M. Vascular endothelial growth factor (VEGF) and its receptor (VEGFR) signaling in angiogenesis: A crucial target for anti- and pro-angiogenic therapies. *Genes Cancer* **2**(12), 1097-1105 (2011).
2. Holmes, D.I.R. and Zachary, I. The vascular endothelial growth factor (VEGF) family: Angiogenic factors in health and disease. *Genome Biol.* **6**(2), 209 (2005).
3. Krilleke, D., DeErkenez, A., Schubert, W., et al. Molecular mapping and functional characterization of the VEGF164 heparin-binding domain. *J. Biol. Chem.* **282**(38), 28045-28056 (2007).
4. Carmeliet, P., Ng, Y.-S., Nuyens, D., et al. Impaired myocardial angiogenesis and ischemic cardiomyopathy in mice lacking the vascular endothelial growth factor isoforms VEGF₁₆₄ and VEGF₁₈₈. *Nat. Med.* **5**(5), 495-502 (1999).
5. Yamamoto, H., Rundqvist, H., Branco, C., et al. Autocrine VEGF isoforms differentially regulate endothelial cell behavior. *Front. Cell. Dev. Biol.* **4**, 99 (2016).
6. Eremina, V., Sood, M., Haigh, J., et al. Glomerular-specific alterations of VEGF-A expression lead to distinct congenital and acquired renal diseases. *J. Clin. Invest.* **111**(5), 707-716 (2003).
7. Ishida, S., Usui, T., Yamashiro, K., et al. VEGF164-mediated inflammation is required for pathological, but not physiological, ischemia-induced retinal neovascularization. *J. Exp. Med.* **198**(3), 483-489 (2003).
8. Osella-Abate, S., Quaglino, P., Savoia, P., et al. VEGF-165 serum levels and tyrosinase expression in melanoma patients: Correlation with the clinical course. *Melanoma Res.* **12**(4), 325-334 (2002).
9. Yoo, S.-A., Kwok, S.-K., and Kim, W.-U. Proinflammatory role of vascular endothelial growth factor in the pathogenesis of rheumatoid arthritis: Prospects for therapeutic intervention. *Mediators Inflamm.* 129873 (2008).