

PRODUCT INFORMATION



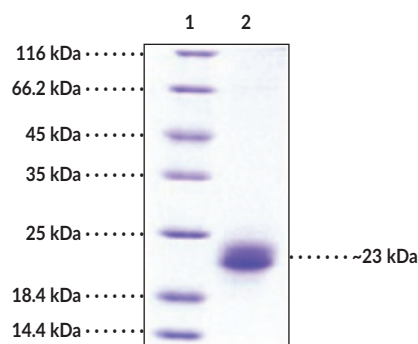
VEGF-C (human, recombinant)

Item No. 32054

Overview and Properties

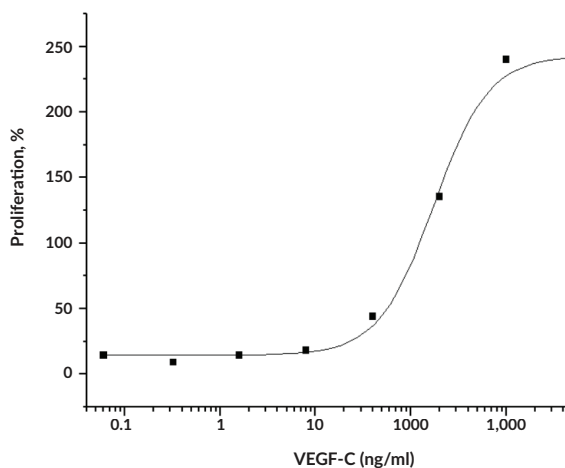
Synonyms: FLT4-L, FLT4 Ligand, Vascular Endothelial Growth Factor C
Source: Active recombinant human C-terminal His-tagged VEGF-C expressed in HEK293 cells
Amino Acids: 103-227
Uniprot No.: P49767
Molecular Weight: 15.5 kDa
Storage: -80°C (as supplied)
Stability: ≥1 year
Purity: ≥95% estimated by SDS-PAGE
Supplied in: Lyophilized from sterile PBS, pH 7.4, with 5% trehalose, 5% mannitol, and 0.01% Tween 80
Endotoxin Testing: <1.0 EU/μg, determined by the LAL endotoxin assay
Bioactivity: Measured by its binding ability in a binding assay. 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-C

SDS-PAGE Analysis of VEGF-C. This protein has a calculated molecular weight of 15.5 kDa. It has an apparent molecular weight of approximately 22 to 24 kDa by SDS-PAGE under reducing conditions due to glycosylation.



Cell Proliferation using HUVECs. The ED₅₀ for this effect is typically 0.1 - 0.5 μg/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, 09/01/2023

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-C is a member of the PDGF/VEGF family of growth factors that promotes lymphangiogenesis.¹ VEGF-C exists as a secreted homodimer composed of a central VEGF homology domain that contains the receptor binding region and is flanked by N- and C-terminal propeptides, which are proteolytically cleaved by furin and ADAMTS3 or plasmin to generate the mature, active protein.^{2,3} VEGF-C is produced by macrophages, fibroblasts, smooth muscle cells, and tumor cells.³ VEGFC expression is upregulated by stimulation with IL-1 β , TNF- α (Item Nos. 32020 | 32069), PDGF, TGF- β , or EGF (Item Nos. 32057 | 32025) and downregulated by dexamethasone (Item No. 11015).^{3,4} Binding of VEGF-C to its receptors, VEGFR2 and VEGFR3, which are expressed by endothelial cells, stimulates the migration, proliferation, and survival of endothelial cells and increases vascular permeability.³ VEGF-C induces angiogenesis in chick embryo chorioallantoic membranes and stimulates proliferation and migration of porcine aortic endothelial cells *in vitro*.⁵ Genome-wide deletion of *Vegfc* in mice induces embryonic edema and defects in lymphatic vascular development and is perinatal lethal.⁶ Increased VEGF-C tumor levels are positively correlated with poor prognosis, decreased overall survival, and resistance to hormone therapy in patients with breast cancer.⁷ VEGFC SNPs have been found in patients with non-small cell lung cancer (NSCLC), as well as colorectal or prostate cancer. Cayman's VEGF-C (human, recombinant) protein can be used for binding and cell-based assay applications. This protein consists of 136 amino acids, has a calculated molecular weight of 15.5 kDa, and a predicted N-terminus of Thr103 after signal peptide cleavage. By SDS-PAGE, under reducing conditions, the apparent molecular mass of the protein is approximately 22 to 24 kDa due to glycosylation.

References

1. Rauniyar, K., Jha, S.K., and Jeltsch, M. Biology of vascular endothelial growth factor C in the morphogenesis of lymphatic vessels. *Front. Bioeng. Biotechnol.* **6**, 7 (2018).
2. Jha, S.K., Rauniyar, K., Karpanen, T., *et al.* Efficient activation of the lymphangiogenic growth factor VEGF-C requires the C-terminal domain of VEGF-C and the N-terminal domain of CCBE1. *Sci. Rep.* **7**(1), 4916 (2017).
3. Narko, K., Enholm, B., Mäkinen, T., *et al.* Effect of inflammatory cytokines on the expression of the vascular endothelial growth factor-C. *Int. J. Exp. Path.* **80**(3), 109-112 (1999).
4. Enholm, B., Paavonen, K., Ristimäki, A., *et al.* Comparison of VEGF, VEGF-B, VEGF-C and Ang-1 mRNA regulation by serum, growth factors, oncoproteins and hypoxia. *Oncogene* **14**(20), 2475-283 (1997).
5. Cao, Y., Linden, P., Farnebo, J., *et al.* Vascular endothelial growth factor C induces angiogenesis *in vivo*. *Proc. Natl. Acad. Sci. USA* **95**(24), 14389-14394 (1998).
6. Karkkainen, M.J., Haiko, P., Sainio, K., *et al.* Vascular endothelial growth factor C is required for sprouting of the first lymphatic vessels from embryonic veins. *Nat. Immunol.* **5**(1), 74-80 (2004).
7. Jain, L., Vargo, C.A., Danesi, R., *et al.* The role of vascular endothelial growth factor SNPs as predictive and prognostic markers for major solid tumors. *Mol. Cancer Ther.* **8**(9), 2496-2508 (2009).

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