

VEGF-C

Recombinant human protein expressed in E. coli cells

Catalog # V812-40N

Lot # S346-4

Product Description

Recombinant human VEGF-c was expressed in E. coli cells and purified by proprietary chromatographic techniques. The protein accession number is [NP_005420.1](#).

Gene Aliases

VEGFC; VRP; Flt4-L

Endotoxin Level

Less than 0.1 ng/µg (IEU/µg) of VEGF-C.

Formulation

Recombinant protein stored in 1 X PBS, 10% glycerol.

Storage and Stability

Stored at desiccated below -20°C. It is stable at 4°C between 2-7 days. For long term storage, aliquot target into smaller quantities after centrifugation and store at recommended temperature. For most favorable performance, avoid repeated handling and multiple freeze/thaw cycles.

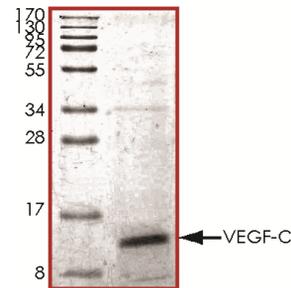
Scientific Background

VEGF-C is a member of the PDGF/VEGF family which is active in angiogenesis and endothelial cell growth, and can also affect the permeability of blood vessels. It is structurally and functionally similar to vascular endothelial growth factor D. VEGFC is an osmosensitive, hypertonicity-driven gene intimately involved in salt-induced hypertension and tonicity-responsive enhancer-binding protein (TONEBP)-VEGFC signaling in mononuclear phagocytes is a major determinant of extracellular volume and blood pressure homeostasis (1). VEGF-C is also essential for embryonic survival and lymphangiogenesis (2).

References

1. Machnik, A. et.al: Macrophages regulate salt-dependent volume and blood pressure by a vascular endothelial growth factor-C-dependent buffering mechanism. *Nature Med.* 15: 545-552, 2009.
2. Karkkainen, M. J. et.al: Vascular endothelial growth factor C is required for sprouting of the first lymphatic vessels from embryonic veins. *Nature Immun.* 5: 74-80, 2004.

Purity



The purity was determined to be **>90%** by densitometry analysis in reducing SDS-PAGE Coomassie blue gel. Approx. MW of the major band was **14 kDa (14,226 Daltons)** by MW calculation).

VEGF-C

Recombinant human protein expressed in E. coli cells

Catalog Number	V812-40N
Specific Lot Number	S346-4
Purity	>90%
Concentration	0.2 µg/µl
Stability	1 yr At -20°C from date of shipment
Storage & Shipping	Stored at desiccated below -20°C. It is stable at 4°C between 2-7 days. For long term storage, aliquot target into smaller quantities after centrifugation and store at recommended temperature. For most favorable performance, avoid repeated handling and multiple freeze/thaw cycles. Product shipped on dry ice.

To place your order, please contact us by phone 1-(604)-232-4600, fax 1-604-232-4601 or by email: orders@signalchem.com
www.signalchem.com

FOR IN VITRO RESEARCH PURPOSES ONLY. NOT INTENDED FOR USE IN HUMAN OR ANIMALS.