

Recombinant Human VEGF (165aa)

Recombinant human protein expressed in *E. coli* cells

Catalog # V815-40N

Lot # G3223-26

Product Description

Recombinant Human VEGF (165aa) was expressed in *E. coli* cells. The protein accession number is [P15692](#)

Gene Aliases

Vascular endothelial growth factor A, VEGF-A, Vascular permeability factor, VPF, VEGFA, VEGF-165aa, Recombinant Human Vascular Endothelial Growth Factor 165AA (VEGFA)

Endotoxin Level

<1.0 EU/µg of recombinant protein as determined by the LAL method

Formulation

Lyophilized from 0.2 µm filtered solution in sodium phosphate and NaCl (pH 6.5)

Reconstitution Protocol

A quick spin of the vial followed by reconstitution in distilled water to a concentration no less than 0.1 mg/mL. This solution can then be diluted into other buffers.

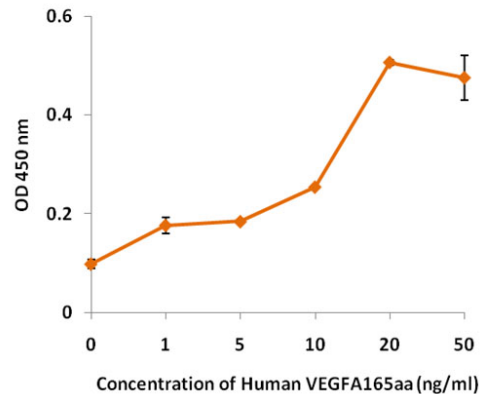
Storage and Stability

The lyophilized protein is stable for at least one year from date of receipt at -70°C. Upon reconstitution, this cytokine can be stored in working aliquots at 2° - 8°C for one month, or at -20°C for six months, with a carrier protein without detectable loss of activity. Avoid repeated freeze/thaw cycles.

Scientific Background

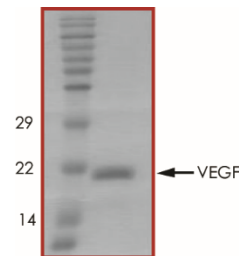
VEGF is a homodimeric glycoprotein that is actively involved in angiogenesis, vasculogenesis and endothelial cell growth. Owing to alternative splicing of the mRNA, VEGF occurs in several molecular variants of 121, 162, 145, 148, 165, 183, 189 and 206 amino acids. The various isoforms of VEGF differ in biological properties such as; (a) the receptors they recognize and (b) their interaction with heparan sulfate proteoglycans. VEGF-121 and VEGF-165 are soluble secreted forms of the factor while VEGF-189 and VEGF-206 are mostly bound to heparin-containing proteoglycans in the cell surface or in the basement membrane. VEGF-165 is the most common variant in most tissues and one with heparin binding properties. Recombinant VEGF-165aa is a non-glycosylated, disulfide linked homodimer.

Specific Activity



The ED50 as determined by the dose-dependent proliferation of human umbilical vein endothelial cells, using a concentration range of 1-100 ng/mL, was found to be <10ng/mL

Purity



Recombinant Human VEGF (165aa) resolved on a 15% SDS-PAGE gel under reducing conditions and stained with Coomassie Brilliant Blue G-250

Approx. MW **19.0 kDa**

Recombinant Human VEGF (165aa)

Recombinant human protein expressed in *E. coli* cells

Catalog Number	V815-40N
Specific Lot Number	G3223-26
Purity	95%
Stability	6 mos. at -20°C from date of shipment
Storage & Shipping	The lyophilized protein is stable for at least one year from date of receipt at -70°C. Upon reconstitution, this cytokine can be stored in working aliquots at 2° - 8°C for one month, or at -20°C for six months, with a carrier protein without detectable loss of activity. Avoid repeated freeze/thaw cycles.

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.