

## Recombinant Human VEGF (121aa)

Recombinant human protein expressed in E. coli cells

**Catalog # V810-40N**

Lot # G3223-25

### Product Description

Recombinant Human VEGF (121aa) was expressed in E. coli cells. The protein accession number is [P15693](#)

### Gene Aliases

Vascular endothelial growth factor A, VEGF-A, Vascular permeability factor, VPF, VEGF-121aa, Recombinant Human Vascular Endothelial Growth Factor 121AA (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 Tris and NaCl, pH 8.0

### Reconstitution Protocol

A quick spin of the vial followed by reconstitution in distilled water to a concentration not less than 0.1 mg/mL. This solution can then be diluted into other buffers.

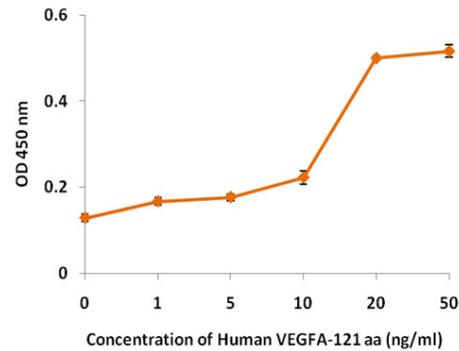
### Storage and Stability

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

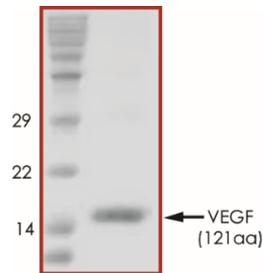
VEGF is a heavily glycosylated, homodimeric protein. The human factor occurs in several molecular variants of 121, 162, 145, 148, 165, 183, 189, 206 amino acids, arising by alternative splicing of the mRNA. The splice forms of VEGF differ in biological properties such as the receptor types, which they recognize and their interaction with heparan sulfate proteoglycans. VEGF121 and VEGF165 are soluble secreted forms of the factor. The interaction of VEGF with heparin-like molecules of the extracellular matrix is required for efficient receptor binding. Protamine sulfate and suramin are capable of replacing the receptor-bound factor. This recombinant construct represents the isoform 121 of VEGFA.

### Specific Activity



The ED50 as determined by the dose-dependent proliferation of human umbilical vein endothelial cells (HUVEC) was found to be 10 ng/mL

### Purity



Recombinant Human VEGF (121aa) resolved on a 15% SDS-PAGE gel under reducing conditions and stained with Coomassie Brilliant Blue G-250

Approx. MW **14.3 kDa**

## Recombinant Human VEGF (121aa)

Recombinant human protein expressed in E. coli cells

Catalog Number	V810-40N
Specific Lot Number	G3223-25
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](mailto:orders@signalchem.com)  
[www.signalchem.com](http://www.signalchem.com)

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