



6042 Cornerstone Court W, Ste B
San Diego, CA 92121
Tel: 1.858.829.3082
Fax: 1.858.481.8694
Email: info@bpsbioscience.com

Human VEGF165 (E. coli derived)

Catalog #: 79516
Lot #: 180507

Description: Recombinant human vascular endothelial growth factor (VEGF)165 expressed in E. coli cells. Human VEGF exists in multiple molecular variants and are named based on the number of amino acids they contain. VEGF165 is the most common form of VEGF. It is a 38.2 kDa disulfide-linked homodimeric protein consisting of two 165 amino acid polypeptides.

Source: Optimized DNA sequence encoding Human VEGF mature chain was expressed in Escherichia Coli

Formulation: Recombinant VEGF lyophilized from 0.2 µm filtered 20 mM PBS solution, pH 7.0.

Reconstitution: 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: The lyophilized protein is stable for at least 2 years from date of receipt at -20° 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..

Purity: >95%, as determined by SDS-PAGE and HPLC.

Endotoxin Level: Endotoxin content was assayed using a LAL gel clot method. Endotoxin level was found to be less than 0.1 ng/µg (1EU/µg).

Biologic Activity: The ED(50) was determined by the dose-dependent proliferation of human umbilical vein endothelial cells (HUVEC) using a concentration range of 1.0-10.0 ng/ml.

References:

1. *Reproduction*. 2009; **138**: 667-677.
2. *Ann. Onc.* 2009; **20**: 1639-1646.

OUR PRODUCTS ARE FOR RESEARCH USE ONLY. NOT FOR DIAGNOSTIC OR THERAPEUTIC USE.

To place your order, please contact us by Phone **1.858.829.3082** Fax **1.858.481.8694**

Or you can Email us at: info@bpsbioscience.com

Please visit our website at: www.bpsbioscience.com