

## Data Sheet

### VEGF165, His-tag (Lyophilized)

Human, Recombinant, N-terminal His-tag  
**Catalog #:** 91001-1  
**Lot #:** 180605  
**Original Conc.:** 0.19 mg/ml

**Original Formulation:** White powder.  
 Lyophilized from 0.19 mg/ml solution in 10 mM Phosphate, pH 7.4, 138 mM NaCl and 2.75 mM KCl

**Reconstitution:** Supplied as a lyophilized solid. Reconstitute in sterile dH<sub>2</sub>O.

**Stability:** ≥12 months at -80°C (as a solid). Store reconstituted protein in aliquots at -80°C for up to 6 months. Avoid freeze/thaw cycles.

**Description:** Human VEGF165, also known as Vascular endothelial growth factor 165, and VEGF, a.a. 27-191 with N-terminal His-Tag. GenBank Accession number NM\_001025368, and expressed in a Baculovirus infected Sf9 cell expression system. MW = 20 kDa.

**Endotoxin Level:** <1 EU/μg.

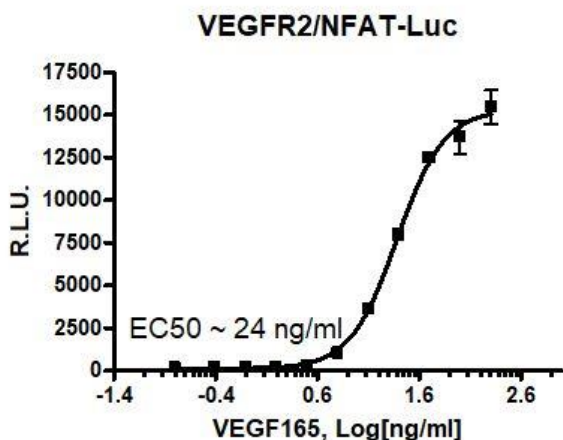
**Assay Conditions:** VEGF165 was tested using VEGFR2/NFAT-Luc HEK293 reporter cell line (BPS cat#79387).

**Application:** Useful for cell culture.

#### References:

1. Ruckman, J., et al. Journal of Biological Chemistry. 273, 20556-20567 (1998).
2. Soker, S., et al. Cellular Biochemistry. 85(2), 357-368 (2002).

## Quality Assurance

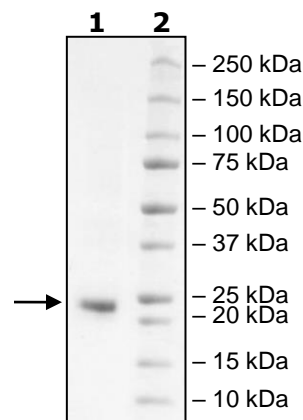


#### 4-20% SDS-PAGE Coomassie staining

**Lane 1:**  
2μg VEGF165

**Lane 2:**  
Protein Marker

**MW:** 20 kDa  
**Purity:** ≥90%



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

To place your order, please contact us by Phone **1.858.202.1401** Fax **1.858.481.8694**

Or you can Email us at: [info@bpsbioscience.com](mailto:info@bpsbioscience.com)

Please visit our website at: [www.bpsbioscience.com](http://www.bpsbioscience.com)