



6405 Mira Mesa Blvd, Ste 100  
San Diego, CA 92121  
Tel: 1.858.202.1401  
Fax: 1.858.481.8694  
Email: support@bpsbioscience.com

## **Rat VEGF-A Recombinant**

**Catalog #:** 91008

**Lot #:** 220905

**Description:** Recombinant rat vascular endothelial growth factor (VEGF) expressed in *E. coli* cells. Rat recombinant VEGF contains 165 amino acids and has a molecular weight of 38.7 kDa. VEGF is an important growth factor involved in both vasculogenesis and angiogenesis.

**Formulation:** Recombinant rat VEGF lyophilized from 10 mM Na Phosphate, pH 7.5.

**Reconstitution:** It is recommended to reconstitute the lyophilized Vascular Endothelial Growth Factor in sterile 18M $\Omega$ -cm H<sub>2</sub>O not less than 100 $\mu$ g/ml, which can then be further diluted to other aqueous solutions.

**Storage:** Stable at room temperature for 3 weeks, should be stored desiccated below -18°C. Upon reconstitution VEGF should be stored at 4°C between 2-7 days and for future use below -18°C. For long term storage it is recommended to add a carrier protein (0.1% HSA or BSA). Avoid freeze/thaw cycles.

**Purity:** >97%, as determined by SDS-PAGE and HPLC.

**Biological Activity:** Determined by the dose-dependent stimulation of the proliferation of human umbilical vein endothelial cells (HUVEC) using a concentration range of 4.0 – 8.0 ng/ml.

### **Protein Sequence:**

MAPTTEGEQKAHEVVVKFMDVYQRSY  
CRPIETLVDIFQEYPDEIEYIFKPSCVPL  
MRCAGCCNDEALECVPTSESNVTMQI  
MRIKPHQSQHIGEMSFLQHSRCECRP  
KKDRTKPEKHCEPCSEERRKHLFVQDP  
QTCKCCKNTDSRCKARQLELNERTC  
RCDKPRR

### **References:**

1. Caires, K.C., *et al. Reproduction.* 2009; **138**: 667-677.
2. Linderholm, B.K., *et al. 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 **1.858.202.1401** Fax **1.858.481.8694**

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

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