

## RayBiotech, Inc.

150 Technology Parkway, Suite#136  
Norcross, GA 30092  
Tel: 770-729-2992, 1-888-494-8555  
Fax: 770-206-2393, 1-888-547-0580  
Email: info@raybiotech.com

# Certificate of Analysis and Data Sheet

## Recombinant Mouse Vacular Endothelial Growth Factor

<b>Source:</b> E.Coli	<b>Catalog No.</b> IP-01-336
--------------------------	---------------------------------

### Background:

Vascular endothelial growth factor-A was originally isolated from tumor cells and referred to as Tumor Angiogenesis Factor or Vascular Permeability Factor. Although expressed at high levels in certain tumor-derived cells it is produced by a wide variety of cell types. In addition to stimulating vascular growth and vascular permeability it may play a role in stimulating Vasolidation via nitric oxide-dependent pathways. Alternative splicing of the mRNA for VEGF-A results in several isoforms of the protein being produced. Rat and bovine VEGF are one amino acid shorter than the human factor, and the bovine and human sequences show a homology of 95 percent. In contrast to other factors mitogenic for endothelial cells such as FGF-1, FGF-2 and PDGF, VEGF is synthesized as a precursor containing a typical hydrophobic secretory signal sequence of 26 amino acids. Glycosylation is not required for efficient secretion of VEGF.

### Description :

Recombinant Murine VEGF produced in E.Coli is a double, non-glycosylated, polypeptide chain containing 165 amino acids and having a molecular mass of 39035 Dalton.

Recombinant VEGF is purified by proprietary techniques.

### Physical Appearance:

Sterile Filtered White Lyophilized (freeze-dried) powder.

### Formulation:

Recombinant VEGF was lyophilized from a concentrated (1mg/ml) solution with no additives.

### Solubility:

It is recommended to reconstitute the lyophilized VEGF in sterile 18M-cm H<sub>2</sub>O not less than 100µg/ml, which can then be further diluted to other aqueous solutions.

### Stability:

Lyophilized VEGF although stable at room temperature for 3 weeks, should be stored desiccated below -18 C. Upon reconstitution murine 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).

Please prevent freeze-thaw cycles.

### Purity:

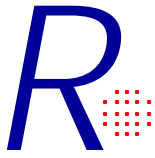
Greater than 98.0% as determined by:

- (a) Analysis by RP-HPLC.
- (b) Anion-exchange FPLC.
- (c) Analysis by reducing and non-reducing SDS-PAGE Silver Stained gel.

### Amino acid sequence:

The sequence of the first five N-terminal amino acids was determined and was found to be Met-Ala-Pro-Thr-Thr.

### Dimers and aggregates:



## RayBiotech, Inc.

150 Technology Parkway, Suite#136  
Norcross, GA 30092  
Tel: 770-729-2992, 1-888-494-8555  
Fax: 770-206-2393, 1-888-547-0580  
Email: [info@raybiotech.com](mailto:info@raybiotech.com)

---

Less than 1% as determined by silver-stained SDS-PAGE gel analysis.

### **Biological Activity:**

The biological activity is determined by the dose-dependent stimulation of the proliferation of human umbilical vein endothelial cells (HUVEC) using a concentration range of 1.0-5.0 ng/ml.

### **Endotoxin:**

Less than 0.1 ng/μg (IEU/μg) of VEGF.

### **Protein content:**

Protein quantitation was carried out by two independent methods:

1. UV spectroscopy at 280 nm using the absorbency value of 0.2875 as the extinction coefficient for a 0.1% (1mg/ml) solution. This value is calculated by the PC GENE computer analysis program of protein sequences (IntelliGenetics).
2. Analysis by RP-HPLC, using a calibrated solution of VEGF as a Reference Standard.

### **Usage:**

This material is offered for research, laboratory or further evaluation purposes.