

## Recombinant Human EG-VEGF

EG-VEGF is an angiogenic growth factor expressed in the ovaries, testis, adrenal, and placental tissues, and is one of main non-glycosylated proteins in the body. EG-VEGF expression in the cell correlates with vascularity in polycystic ovarian syndrome, a leading cause of infertility.

### Materials Provided

<b>Catalog Number:</b>	<b>92-1022</b>
<b>Quantity Supplied:</b>	1 vial(s), 5 µg per vial
<b>Reconstitution Buffer:</b>	1 mL

### Description

DiscoverX control ligands have been functionally tested and validated for optimal use with all cell line targets and Assay Ready kits.

### Product Information

<b>Molecular Weight:</b>	9600 g/mol
<b>Source:</b>	E. coli
<b>Purity:</b>	≥ 98% by SDS-PAGE and HPLC
<b>Endotoxin Level:</b>	< 0.1 ng/µg (<1EU/µg).
<b>Formulation:</b>	Sterile filtered through a 0.2 micron filter. Lyophilized with no additives
<b>Storage:</b>	≤ -20°C Please avoid multiple freeze/thaw cycles.

### Reagent Preparation

To avoid condensation, equilibrate the vial to ambient temperature before opening.

<b>Stock Concentration:</b>	20 µM
<b>Reconstitution Volume:</b>	0.026 mL
<b>Reconstitution Solvent:</b>	Reconstitution Buffer

***For Research Use Only***

Ordering: +1.510.979.1415 option 4 or e-mail [CustomerServiceDRX@eurofins.com](mailto:CustomerServiceDRX@eurofins.com)  
Technical support: +1.510.979.1415 option 5 or e-mail [DRX\\_SupportUS@eurofinsUS.com](mailto:DRX_SupportUS@eurofinsUS.com)  
General product information: [www.discoverx.com](http://www.discoverx.com)