



6044 Cornerstone Court W, Ste E
San Diego, CA 92121
Tel: 1.858.829.3082
Fax: 1.858.481.8694
Email: info@bpsbioscience.com

Human VEGF121

Catalog #: 91005
Lot #: 140925

Description: Recombinant human vascular endothelial growth factor (VEGF)121 expressed in CHO cells. Human VEGF exists in multiple molecular variants and are named based on the number of amino acids they contain. VEGF121 is the smallest and most soluble form of VEGF.

Source: Optimized DNA sequence encoding human VEGF121 mature chain was expressed in CHO cells.

Formulation: Lyophilized powder

Reconstitution: Briefly centrifuge the vial, followed by reconstitution in distilled water to a concentration of ≥ 0.1 mg/ml. This solution can then be further diluted into other buffers containing carrier protein (0.1% BSA).

Storage: The lyophilized protein is stable for at least 2 years from date of receipt when stored at -20°C . Upon reconstitution, store in working aliquots at $2 - 8^{\circ}\text{C}$ for up to one month, or at -20°C for up to six months, in the presence of a carrier protein. Avoid repeated freeze/thaw cycles.

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

Endotoxin Level: Endotoxin level was found to be < 0.1 ng/ μg (1EU/ μg), using the LAL gel clot method.

Protein Sequence:

APMAEGGGQNHHEVVKFMDVYQRSY
CHPIETLVDIFQEYPDEIEYIFKPSCVPL
MRCGGCCNDEGLECVPTESNITMQI
MRIKPHQQQHIGEMSFLQHNKCECRP
KKDRARQENPCGPCSERRKHLFQDP
QTCKCCKNTDSRCKARQLELNERTC
RCDKPRR

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