

## Data Sheet

### **Neuropilin-1, Avi-His-Tag, Biotin-labeled**

Human, Recombinant, C-terminal Avi-His-tag,  
Biotin-labeled

**Catalog #:** 100911

**Lot #:** 201029-2      **Conc.:** 1.40 mg/ml

**Description:** Human Neuropilin 1 (NRP1), also known as CD304 and VEGF165R, GenBank Accession No. NM\_001024628.3, a.a. 22-644(end) with C-terminal Avi-His-tag, expressed in a HEK293 expression system. MW=75 kDa. This protein is enzymatically biotinylated using Avi-Tag™ technology. Biotinylation is confirmed to be ≥90%.

**Formulated in:** 8 mM phosphate pH 7.4, 110 mM NaCl, 2.2 mM KCl, and 20% glycerol.

**Stability:** At least 6 months at –80°C. Avoid freeze/thaw cycles.

#### **References:**

1. Cantuti-Castelvetri, L. *et al.*, *BioRxiv*. doi.org/10.1101/2020.06.07.137802 (2020).
2. Weissenstein, B., *et al.*, *J. Regan Biol Med.* **2(4)**, 1-2 (2020).

## **Quality Assurance**

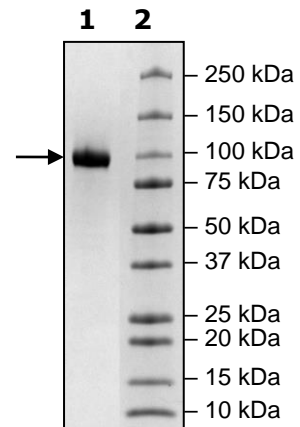
### **4-20% SDS-PAGE Coomassie staining**

**Lane 1:**  
4µg NRP1

**Lane 2:**  
Protein Marker

**MW:** 75 kDa +glycans

**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 **1.858.259.5734**

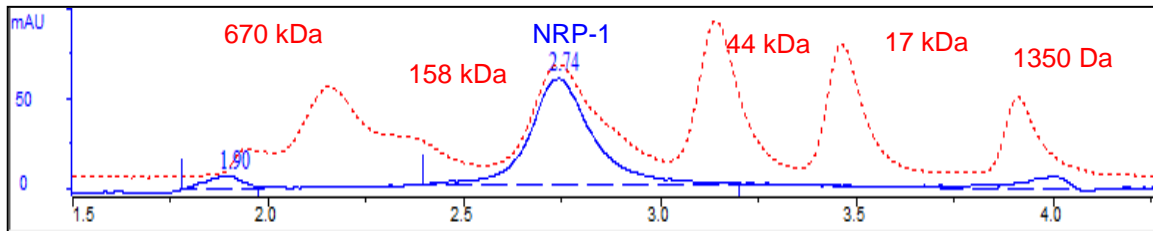
Or you can Email us at: [info@bpsbioscience.com](mailto:info@bpsbioscience.com)

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



6042 Cornerstone Court West, Suite B  
San Diego, CA 92121  
Tel: 1.858.202.1401  
Fax: 1.858.481.8694  
Email: [info@bpsbioscience.com](mailto:info@bpsbioscience.com)

### Gel Filtration Trace



<10% Aggregation

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 **1.858.259.5734**  
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)