

## Data Sheet

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

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

**Catalog #:** 100870**Lot #:** 200923**Conc.:** 2.67 mg/ml**Formulated in:** 8 mM phosphate pH 7.4,  
110 mM NaCl, 2.2 mM KCl, and 20%  
glycerol.**Stability:** At least 6 months at  $-80^{\circ}\text{C}$ . Avoid  
freeze/thaw cycles.**References:**

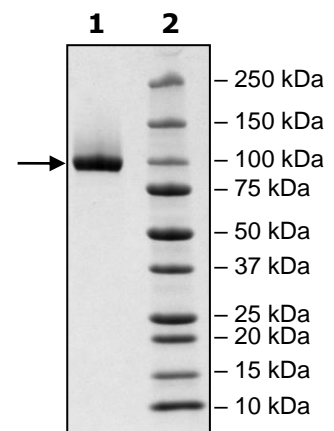
1. Cantuti-Castelvetri, L. *et al.*, *BioRxiv*.  
doi.org/10.1101/2020.06.07.137802 (2020).
2. Weissenstein, B., *et al.*, *J. Regan Biol  
Med.* **2(4)**, 1-2 (2020).

**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.

## **Quality Assurance**

**4-20% SDS-PAGE  
Coomassie staining****Lane 1:**4 $\mu\text{g}$  NRP1**Lane 2:**

Protein Marker

**MW:** 75 kDa**Purity:**  $\geq 90\%$ 

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)