

Data Sheet

Nectin4, His-Avi-Tag

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

Catalog #: 100674

Lot#: 200407

Conc.: 3.41 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°C . Avoid freeze/thaw cycles.

References:

1. Mühlebach, M.D., *et al.*, *Nature*. 2011; **480**: 530-533
2. Noyce, R.S., *et al.*, *PLoS Pathog.* 2011; **7(8)**: e1002240.

Description: Human Nectin Cell Adhesion Molecule 4 (Nectin4), also known as Poliovirus Receptor-Related 4 (PVL4) and Ig superfamily receptor LNIR, GenBank Accession No. NM_030916, a.a. 32-349 with a C-terminal His-Avi-TagTM, expressed in a HEK293 cell expression system. MW=37 kDa. This protein runs at a higher M.W. by SDS-PAGE due to glycosylation.

Quality Assurance

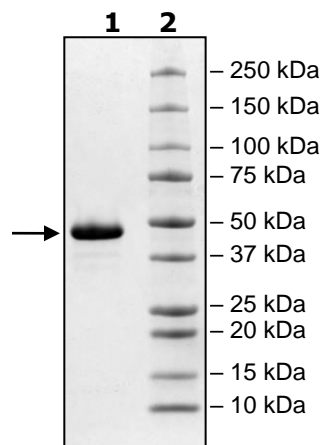
**4-20% SDS-PAGE
Coomassie staining**

Lane 1:
4 μg Nectin4

Lane 2:
Protein Marker

MW: 37 kDa + glycans

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 **1.858.481.8694**

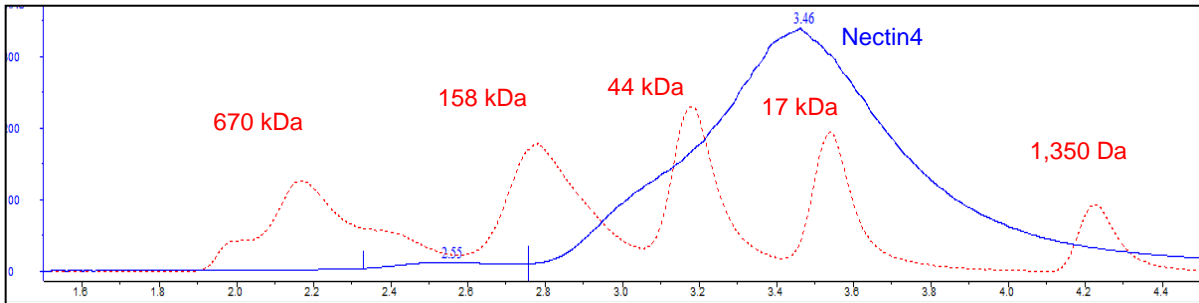
Or you can Email us at: info@bpsbioscience.com

Please visit our website at: www.bpsbioscience.com



6042 Cornerstone Court W, Ste B
San Diego, CA 92121
Tel: 1.858.202.1401
Fax: 1.858.481.8694
Email: info@bpsbioscience.com

Gel Filtration Curve



<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 **1.858.481.8694**
Or you can Email us at: info@bpsbioscience.com
Please visit our website at: www.bpsbioscience.com