

Data Sheet

B7-H5, Avi-His-Tag, Biotin-Labeled

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

Catalog #: 100359

Lot #: 190425-2

Conc.: 1.67 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°C . *Avoid
freeze/thaw cycles.*

Description: Human B7-H5 also known as VISTA, GI24, Dies1, PD-1H, SISP1, DD1alpha, and VSIR, with C-terminal Avi-His-Tag, GenBank Accession No. NM_022153, a.a 33-194, expressed in a HEK293 cell expression system. This protein is enzymatically biotinylated using Avi-Tag™ technology. Biotinylation is confirmed to be $\geq 90\%$. MW = 21 kDa. This protein runs at a higher molecular weight due to glycosylation.

References:

1. Zhu, Y., et al. *Nature Communications*. 2013 Feb 27; **4**.
2. Jung, K., et al., *Immune Network*. 2013, Oct. 13;**6**:184-193.

Quality Assurance

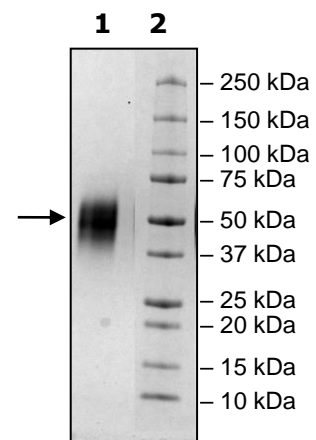
4-20% SDS-PAGE Coomassie staining

Lane 1:
4 μg VISTA

Lane 2:
Protein Marker

MW: 21 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**

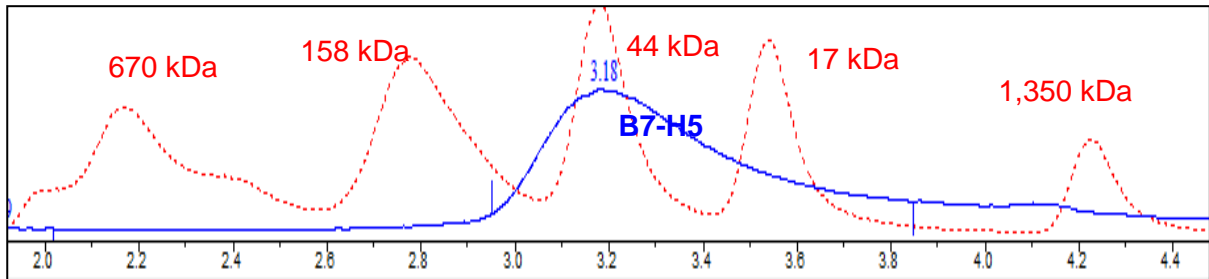
Or you can Email us at: info@bpsbioscience.com

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