

Data Sheet

B7-H5, Avi-His-Tag

Human, Recombinant, C-terminal Avi-His-Tag
Catalog #: 100358
Lot #: 190424-2 **Conc.:** 1.48 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. Zhu, Y., et al. *Nature Communications*.
2013 Feb 27; **4**.
2. Jung, K., et al., *Immune Network*. 2013,
Oct. 13; **6**:184-193.

Description: Human B7-H5 also known as VISTA, GI24, B7-H5, 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. MW = 21 kDa. This protein runs at a higher MW by SDS-PAGE due to glycosylation.

Quality Assurance

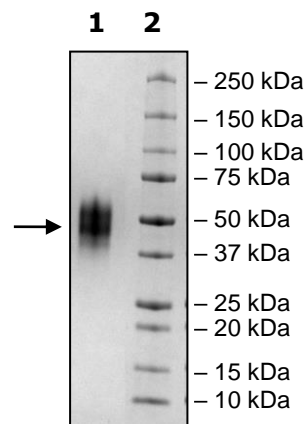
4-20% SDS-PAGE Coomassie staining

Lane 1:
4 μg B7-H5

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