

## Data Sheet

### **B7-H7, Fc-Fusion, Avi-Tag, Biotin-Labeled**

Human, Recombinant, C-terminal Fc-Avi-Tag,  
Biotin-Labeled

**Catalog #:** 79366

**Lot #:** 150806

**Conc.:** 0.56 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. Protein may be diluted to  
 $\geq 100 \mu\text{g/ml}$  in PBS + glycerol and stored at  
 $-80^{\circ}\text{C}$ .

**References:**

1. Flajnik, M.F., *et al. Immunogenetics.*  
2012; **64(8)**: 571-590.
2. Schluter, F., *et al., Inorg. Chem.,*  
2011; **50(6)**: 2580-2589.

**Description:** Human B7-H7, also known as HHLA2, and HERV-H LTR-associating 2, GenBank Accession No. NM\_007072, a.a. 23-344, with a C-terminal Avi-Tag fused with the Fc portion of Human IgG1, expressed in a HEK293 cell expression system. This protein is enzymatically biotinylated using Avi-Tag™ technology. MW =66 kDa. This protein runs at a higher MW by SDS-PAGE due to glycosylation.

**Applications:** Useful for the study of protein-protein interactions and in vitro and in vivo protein function.

## **Quality Assurance**

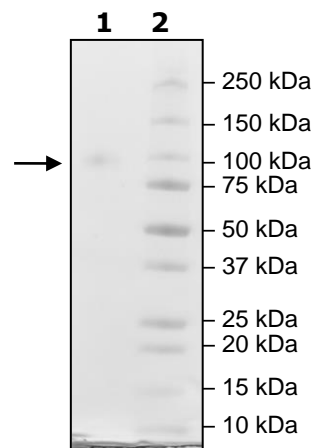
**4-20% SDS-PAGE  
Coomassie staining**

**Lane 1:**  
3  $\mu\text{g}$  B7-H7

**Lane 2:**  
Protein Marker

**MW:** 65.5 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 **1.858.829.3082** Fax **1.858.481.8694**

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 W, Ste B  
San Diego, CA 92121  
**Tel:** 1.858.829.3082  
**Fax:** 1.858.481.8694  
**Email:** [info@bpsbioscience.com](mailto:info@bpsbioscience.com)

### Sequence

IFPLAFFIYVPMNEQIVIGRLDEDIILPSSFERGVSEVVIHWKYQDSYKVHSYYKG  
SDHLESQDPRYANRTSLFYNEIQNGNASLFFRRVSLLEDEGIYTCYVGTAIQVITN  
KVVLKVGVFLLTPVMKYEKRNTNSFLICSVLSVYPRPIITWKMDNTPISENMEET  
GSLDSFSINSPLNITGSNSSYECTIENSLLKQTWTGRWTMKDGLHKMQSEHVSLS  
CQPVNDYFSPNQDFKVTWSRMKSGTFSVLAYYLSSSQNTIINESRFSWNKELINQ  
SDFSMNLMDLNLSDSGEYLCNISSDEYTLTITHTVHVEPSQETASHNIEGRMDPK  
SCDKTHTCPPCPAPELLGGPSVFLFPPKPKDTLMI SRTPEVTCVVVDVSHEDPEV  
KFNWYVDGVEVHNAKTKPREEQYNSTYRVVSVLTVLHQDWLNGKEYKCKVSNKAL  
PAPIEKTISKAKGQPREPQVYTLPPSRDELTKNQVSLTCLVKGFYPSDIAVEWES  
NGQPENNYKTTTPVLDSDGSFFLYSKLTVDKSRWQQGNVFSQVMSVHEALHNHYTQ  
KSLSLSPGKGGGLNDIFEAQKIEWHE

**Black:** B7-H7 (23-344)

**Green:** linker

**Red:** Fc of Human IgG1 (a.a. 100-330)

**Blue:** Avi-tag

OUR PRODUCTS ARE FOR RESEARCH USE ONLY. NOT FOR DIAGNOSTIC OR THERAPEUTIC USE.

To place your order, please contact us by Phone **1.858.829.3082** Fax **1.858.481.8694**

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)