

Data Sheet

B7-H7, Fc-Fusion, Avi-Tag

Human, Recombinant, C-terminal Fc-Avi-Tag
Catalog #: 79365
Lot #: 150805 **Conc.:** 0.2 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. Protein may be diluted to $\geq 100 \mu\text{g/ml}$ in PBS + glycerol and stored at -80°C .

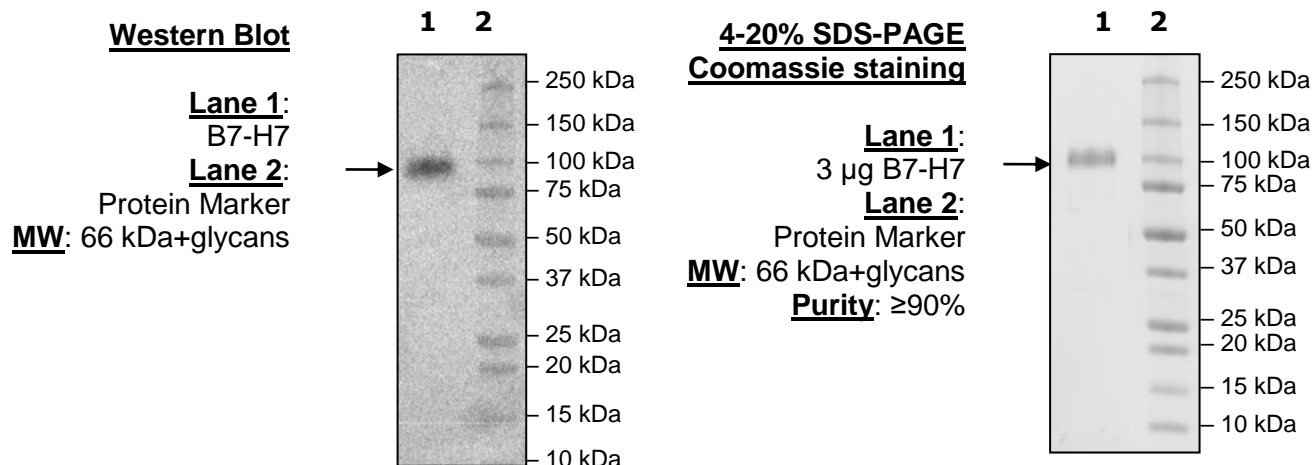
References:

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, 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. MW =66 kDa.

Applications: Useful as a positive control for Western blot studies and in protein binding studies.

Quality Assurance



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
 Please visit our website at: 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

Sequence

IFPLAFFIYVPMNEQIVIGRLDEDIILPSSFERGVSEVVIHWKYQDSYKVHSYYKG
SDHLESQDPRYANRTSLFYNEIQNGNASLFFRRVSLLEDEGIYTCYVGTAIQVITN
KVVLKVGVFELTPVMKYEKRNTNSFLICSVLSVYPRPIITWKMDNTPISENMEET
GSLDSFSINSPLNITGSNSSYECTIENSLLKQTWTGRWTMKDGLHKMQSEHVSLS
CQPVNDYFSPNQDFKVTWSRMKSGTFSVLAYYLSSQNTIINESRFSWNKELINQ
SDFSMNLMDLNLSDSGEYLCNISSDEYTLTITHTVHVEPSQETASHNIEGRMDPK
SCDKTHTCPPCPAPELLGGPSVFLFPPKPKDTLMI SRTPEVTCVVVDVSHEDPEV
KFNWYVDGVEVHNAKTKPREEQYNSTYRVVSVLTVLHQDWLNGKEYKCKVSNKAL
PAPIEKTISKAKGQPREPQVYTLPPSRDELTKNQVSLTCLVKGFYPSDIAVEWES
NGQPENNYKTTTPVLDSDGSFFLYSKLTVDKSRWQQGNVFSVMSVHEALHNHYTQ
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

Please visit our website at: www.bpsbioscience.com