

Data Sheet

B7-H6, Fc-Fusion-Avi-Tag (Human)

Human, Recombinant, C-terminal Fc-Fusion-Avi-Tag

Catalog #: 79383

Lot #: 171212-G2 **Conc.:** 0.98 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, J. Immunogenetics., 2012. Aug 64;8:571-590.
2. Wu, M.R., et al., Gene Therapy. 2015. Apr 22:675-684.

Description:

Human B7-H6, also known as NCR3LG1, B7 homolog 6, and natural killer cell cytotoxicity receptor. GenBank Accession No. NM_001202439, a.a. 25-262 with C-terminal Fc-Avi-Tag, expressed in a HEK293 cell expression system. MW = 52 kDa. This protein runs at a higher MW by SDS-PAGE due to glycosylation.

Quality Assurance

4-20% SDS-PAGE Coomassie staining

Lane 1:

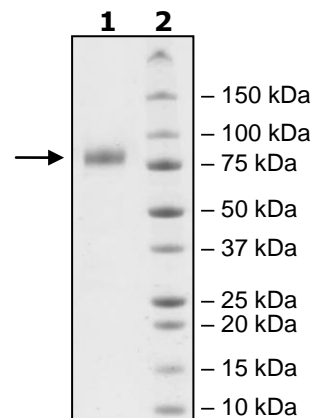
2 μg B7-H6

Lane 2:

Protein Marker

MW: 52 kDa + glycans

Purity: $\geq 90\%$



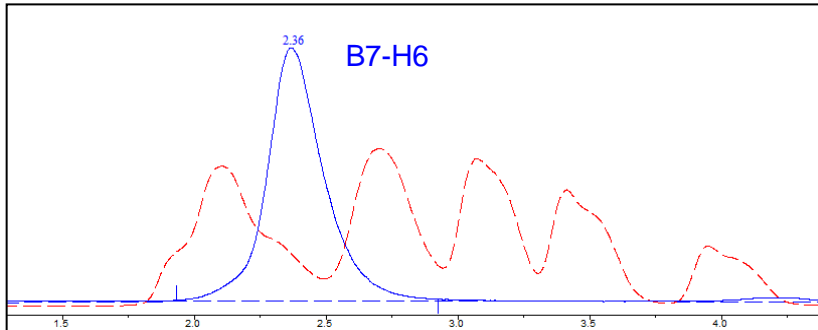
OUR PRODUCTS ARE FOR RESEARCH USE ONLY. NOT FOR DIAGNOSTIC OR THERAPEUTIC USE.

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

Please visit our website at: www.bpsbioscience.com

Gel Filtration Curve



<10% Aggregation

Sequence

DLKVEMMAGGTQITPLNDNVTIFCNIFYSQPLNITSMGITWFWKSLTFDKE
 VKVFEFFGDHQEAFRPGAIVSPWRLKSGDASLRLPGIQLEEAGEYRCEV
 VVTPLKAQGTVQLEVVASRLLLDQVGMKENEDKYMCESSGFYPEAI
 NITWEKQTQKFPHPHIEISEDVITGPTIKNMDGTFNVTSCCLKLNSSQEDPGT
 VYQCVVRHASLHTPLRSNFTLTAARHSLSETEKTDNFS**IEGRMDPKSCD**
KTHTCPPCPAPELLGGPSVFLFPPKPKDTLMISRTPEVTCVVVDVSHEDP
EVKFNWYVDGVEVHNAKTKPREEQYNSTYRVVSVLTVLHQDWLNGKEY
KCKVSNKALPAPIEKTISKAKGQPREPQVYTLPPSRDELTKNQVSLTCLV
KGFYPSDIAVEWESNGQPENNYKTTTPVLDSDGSFFLYSKLTVDKSRWQ
QGNVFSCSVMHEALHNHYTQKSLSLSPGKGGGLNDIFEAQKIEWHE

Black: B7-H6 (25-262)

Green: linker

Red: Fc

Blue: Avi

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