

Data Sheet

B7-H5 Fc Fusion (IgG1), Avi-Tag, Biotin-labeled

Human, Recombinant, Fc fusion,
C-terminal Avi-Tag, Biotin labeled.

Catalog #: 71327

Lot #: 170330

Conc.: 2.91 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. 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 secreted B7-H5 Fc fusion protein, also known as VISTA and VSIR, with C-terminal Avi-Tag. GenBank Accession No. NM_022153, a.a 33-194 expressed in a HEK293 cell expression system. This protein is fused at the C-terminus to the Fc region of human IgG1 and is enzymatically biotinylated using Avi-tag™ technology. MW = 47 kDa. This protein runs at a higher M.W. by SDS-PAGE due to glycosylation.

Application: Useful for the study of protein-protein interaction and in vitro and in vivo protein function studies.

Quality Assurance

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

Lane 1:

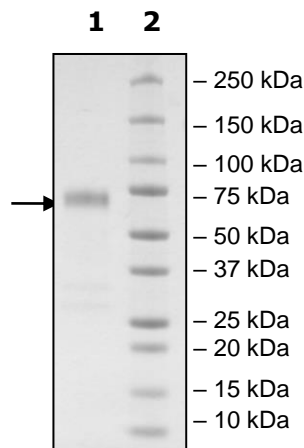
3 μg B7H5

Lane 2:

Protein Marker

MW: 47 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