

Data Sheet

Human CD226, Fc fusion, Biotin-labeled

Human, recombinant, Fc fusion protein,
C-terminal AviTag™

Catalog #: 71253

Lot #: 160314-G

Conc.: 1.04 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
≥ 100 µg/ml in PBS + glycerol and stored at -
80°C.

References:

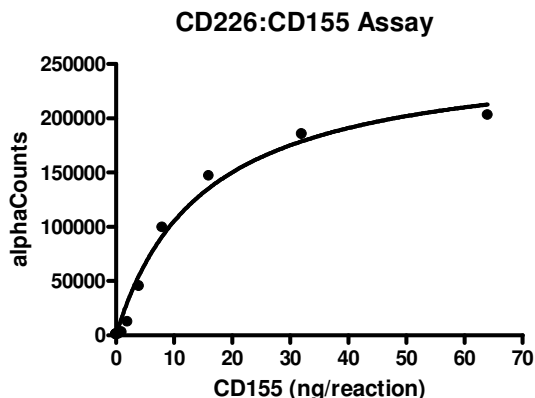
1. Bottino, C., *et al.*, *J. Exp. Med.* 2003;
198(4): 557-567.
2. Hafler, J., *et al.*, *Genes Immun.* 2009;
10: 5-10.

Description: Human CD226 antigen also
known as DNAX accessory molecule 1
(DNAM-1), GenBank Accession No.
NM_006566, a.a. 19-247 fused to Fc region
of human IgG1, this protein is enzymatically
biotinylated using AviTag™ technology and
expressed in a HEK293 cell expression
system. MW = 54.7 kDa (monomer). This
protein runs at a higher MW by SDS-PAGE
due to glycosylation.

Assay Conditions: CD226, biotin binding
to CD155, His-tag was confirmed according
to the protocol for BPS Cat. 72052. Briefly,
6 ng of CD226-biotin was mixed with
varying amounts of CD155, His-tag in
CD226 assay buffer at a final volume of 10
µl and allowed to incubate for 1 hr at room
temp. Nickle acceptor beads were added for
30 min, followed by Strep-conjugated
acceptor beads for 30 min and reading
alpha-counts.

Application: Useful for studying protein
binding and for screening small molecules
and antibodies.

Quality Assurance

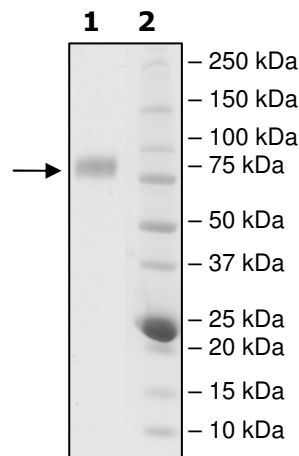


**4-20% SDS-PAGE
Coomassie staining**

Lane 1:
2 µg CD226

Lane 2:
Protein Marker

MW: 54.7 kDa+glycans
Purity: ≥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**

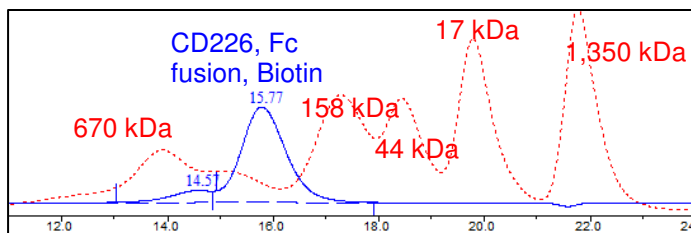
Or you can Email us at: info@bpsbioscience.com

Please visit our website at: www.bpsbioscience.com



6044 Cornerstone Court W, Ste E
San Diego, CA 92121
Tel: 1.858.829.3082
Fax: 1.858.481.8694
Email: info@bpsbioscience.com

Gel Filtration Curve



<10% aggregation

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)