

Data Sheet

CD150, Fc Fusion, Biotin-Labeled

Human, Recombinant, Fc fusion protein,
C-terminal Avi- tag™

Catalog #: 72307

Lot #: 160421

Conc.: 0.94 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. Tatsuo, H. *et al.*, *Journal of Virology*.
2001; **75**: 5842-5850.
2. Sidorenko, S. *et al.*, *Nature Immunology*.
2003; **4**: 19-24.

Description:

Human CD150, Fc fusion protein, also
known as signaling lymphocytic activation
molecule (SLAM) or SLAMF1. GenBank
Accession No. NM_003037, a.a. 21-237
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 = 53 kDa. This
protein runs at a higher MW by SDS-PAGE
due to glycosylation.

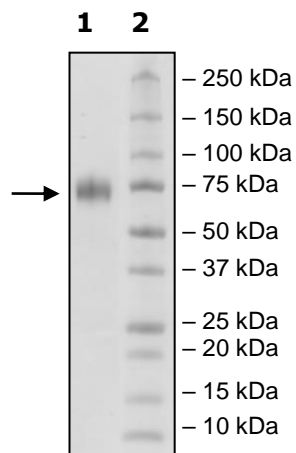
Application:

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

Quality Assurance

4-20% SDS-PAGE
Coomassie staining

Lane 1:
3µg CD150
Lane 2:
Protein Marker
MW: 53 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