

Data Sheet

CD48, Fc Fusion, Avi-Tag, HiP™

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

Catalog #: 72309

Lot #: 160419

Conc.: 1.33 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. Nakajima, H. *et al.*, *European Journal of Immunology*. 1999; **29(5)**: 1676-1683.
2. Brown, M.H., *et al.*, *Journal of Experimental Medicine*. 1998; **188(11)**: 2083.

Description:

Human CD48, Fc fusion protein, also known as Cluster of Differentiation 48, BLAST-1, or SLAMF2. GenBank Accession No. NM_001778, a.a. 27-220 expressed in a HEK293 cell expression system. This protein is fused at the C-terminus to the Fc region of human IgG1 using Avi-tag™ technology. MW = 51 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:

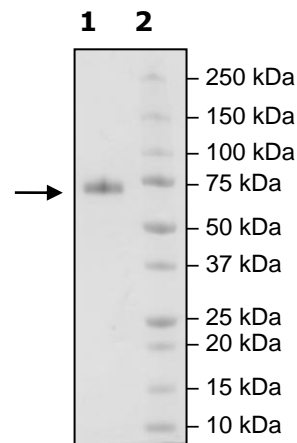
3 μg CD48

Lane 2:

Protein Marker

MW: 51 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**

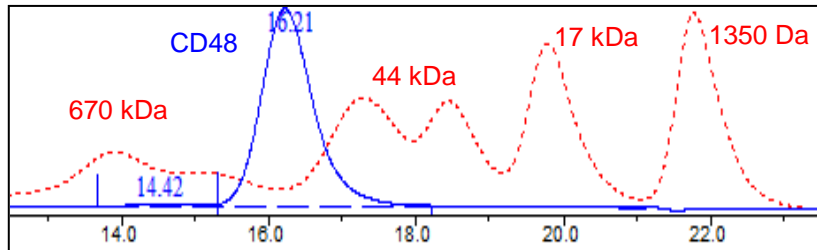
Or you can Email us at: info@bpsbioscience.com

Please visit our website at: www.bpsbioscience.com



6042 Cornerstone Court W, Ste B
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
Please visit our website at: www.bpsbioscience.com