

Data Sheet

CD24, Fc Fusion, Avi-tag, Biotin-labeled

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

Catalog #: 100528

Lot #: 191024

Conc.: 1.62 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.

References:

1. Chen, G-Y., *et al.*, *Science*. 2009
March; **323(5922)**:1722-1725.
2. Phillips, T.M., *et al.*, *JNCI*. 2006 Dec;
98(24):1777-1785.

Description: Human secreted CD24, also known as Signal Transducer CD24, Small Cell Lung Carcinoma Cluster 4 Antigen and CD24A, GenBank Accession No. NM_013230, a.a. 27-59(end), fused at the C-terminus of the Fc portion of human IgG1, with C-terminal Avi-tag, expressed in a HEK293 cell expression system and enzymatically biotin-labeled using Avi-tag™ technology. Biotinylation is confirmed to be $\geq 90\%$. MW = 32 kDa. This protein runs at a higher MW by SDS-PAGE due to glycosylation.

Quality Assurance

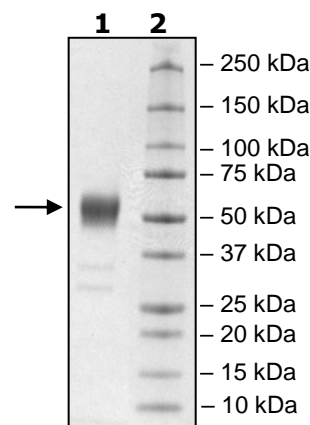
**4-20% SDS-PAGE
Coomassie staining**

Lane 1:
4 μg CD24

Lane 2:
Protein Marker

MW: 32 kDa + glycans

Purity: $\geq 90\%$



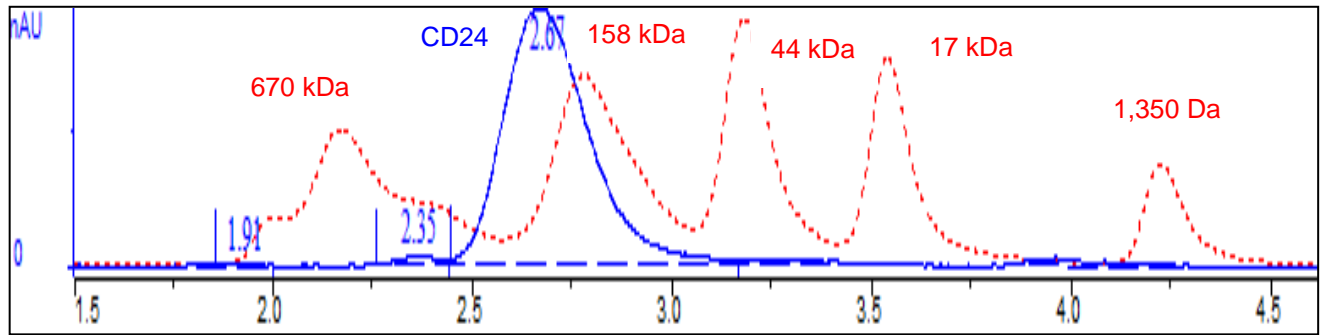
OUR PRODUCTS ARE FOR RESEARCH USE ONLY. NOT FOR DIAGNOSTIC OR THERAPEUTIC USE.

To place your order, please contact us by Phone **1.858.202.1401** 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: 1.858.202.1401
Fax: 1.858.481.8694
Email: info@bpsbioscience.com

Gel Filtration:



<10% Aggregation

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

To place your order, please contact us by Phone **1.858.202.1401** Fax **1.858.481.8694**
Or you can Email us at: info@bpsbioscience.com
Please visit our website at: www.bpsbioscience.com