

Data Sheet

CD44, Fc fusion, Biotin-labeled

Human, recombinant, Fc fusion, biotin labeled

Catalog #: 71133

Lot #: 150312

Conc.: 2.84 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. Aruffo, A., *et al.*, *Cell*. 1990; **61(7)**: 1303-1313.
2. Sreaton, G.R., *et al.*, *Proc. Natl. Acad. Sci. (USA)*. 1992; **89(24)**: 12160-12164.

Description:

Human secreted CD44, also known as epican, extracellular matrix receptor III, GP90 lymphocyte homing/adhesion receptor, HUTCH-I, or phagocytic glycoprotein I (PGP-1), GenBank Accession No. NM_000610, a.a. 21-220 expressed in a HEK293 cell expression system, fused at the C-terminus to the Fc portion of human IgG1, and enzymatically biotin-labeled using Avi-tag™ technology. MW = 50.6 kDa (monomer). This protein runs at a higher M.W. by SDS-PAGE due to glycosylation.

Application:

Useful for the study of enzyme kinetics, screening inhibitors, and selectivity profiling.

Quality Assurance

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

Lane 1:

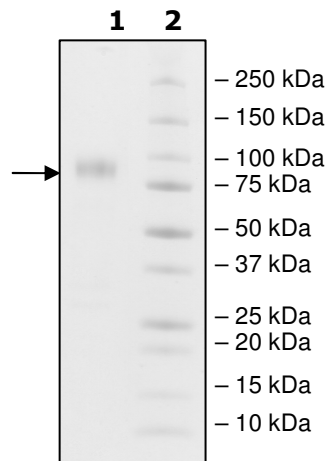
3 μg CD44

Lane 2:

Protein Marker

MW: 50.6 kDa+glycans

Purity: $\geq 90\%$



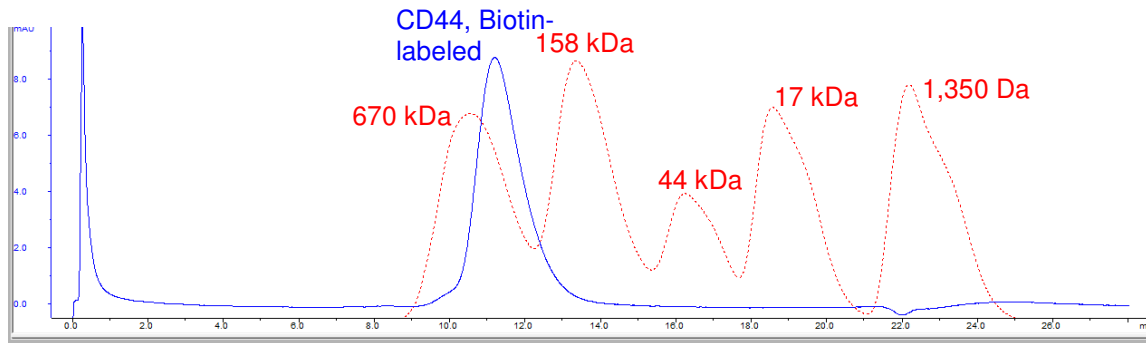
OUR PRODUCTS ARE FOR RESEARCH USE ONLY. NOT FOR DIAGNOSTIC OR THERAPEUTIC USE.

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

Gel Filtration Curve



Sequence

QIDLNITCRFAGVFHVEKNGRYSISRTEAADLCKAFNSTLPTMAQMEKALSIGFETCRY
 GFIEGHVVIPRIHPNSICAANNTGVYILTSNTSQYDYCFNASAPPEEDCTSVTDLPNAF
 DGPITITIVNRDGTTRYVQKGEYRTNPEDIYPSNPTDDDVS SSSERSSTSGGYIFYTFS
 TVHPIPEDSPWITDSTDRIPIEGRMDPKSCDKTHTCPPCPAPELLGGPSVFLFPPKPK
 DTLMISRTPEVTCVVVDVSHEDPEVKFNWYVDGVEVHNAKTKPREEQYNSTYRVVSVL
 TVLHQDWLNGKEYKCKVSNKALPAPIEKTISKAKGQPREPQVYTLPPSRDELTKNQVSL
 TCLVKGFYPSDIAVEWESNGQPENNYKTTTPVLDSDGSFFLYSKLTVDKSRWQQGNV
 FSCSVMHEALHNHYTQKSLSLSPGKGGGLNDIFEAQKIEWHE

Black: CD44 (a.a. 21-220)
 Green: linker
 Red: Fc of human IgG1 (a.a. 100-330)
 Blue: Avi-tag™ sequence

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