

Data Sheet

CD44

Human, recombinant, Fc fusion protein

Catalog #: 71187

Lot #: 140624-D

Conc.: 1.27 mg/ml

Formulated in: 40 mM Tris-HCl, pH 8.0,
110 mM NaCl, 2.2 mM KCl, and 20%
glycerol.

Stability: >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 .*

Endotoxin: <1 EU/ μg

References:

1. Prince, M.E., *et al.*, *Proc. Natl. Acad. Sci. (USA)*. 2006; **104(3)**: 973-978.
2. Lei, D., *et al.*, *Clin. Canc. Res.* 2008; **14**: 6751-6760.

Description:

Human CD44, also known as epican, extracellular matrix receptor III, GP90 lymphocyte homing/adhesion receptor, HUTCH-I, or phagocytic glyco-protein I (PGP-1), GenBank Accession No. NM_000610, a.a. 21-220, fused with Fc region of human IgG, expressed in a HEK293 cell expression system. MW = 48.7 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:

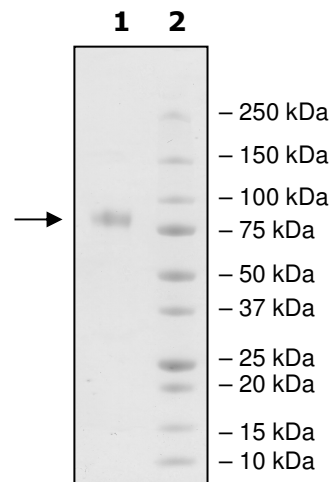
2 μg CD44

Lane 2:

Protein Marker

MW: 48.7 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