

Data Sheet

CD244, Fc Fusion, Avi-Tag

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

Catalog #: 72305

Lot #: 160419

Conc.: 0.59 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. Eissmann, P. *et al.*, *Blood*. 2004; **105**:
4722-4729.
2. Watzl, C. *et al.*, *Journal of Immunology*.
2000; **165(7)**: 3545-3548.

Description:

Human CD244, Fc fusion protein, also
known as Natural Killer Cell Receptor 2B4.
GenBank Accession No. NM_016382, a.a.
22-229 expressed in a HEK293 cell
expression system. This protein is fused at
the C-terminus to the Fc region of human
IgG1. MW = 52 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:

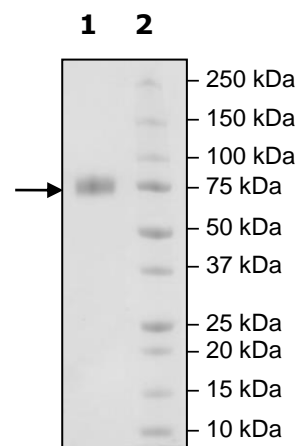
3 μg CD244

Lane 2:

Protein Marker

MW: 52 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