

Data Sheet

CD244, Fc Fusion, Biotin-Labeled

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

Catalog #: 72306

Lot #: 160421

Conc.: 2.03 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:

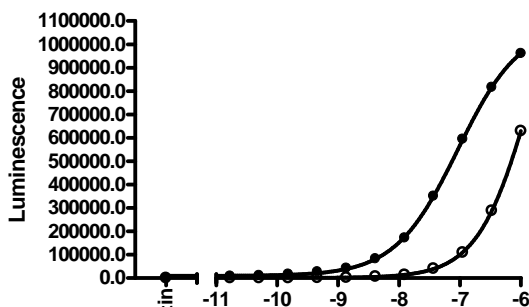
Human CD244, Fc fusion protein, also
 known as Natural Killer Cell Receptor 2B4.
 GenBank Accession No. NM_016382, a.a.
 22-229 expressed in a HEK293 cell
 expression system. This protein is fused at
 the C-terminus to the Fc region of human
 IgG1 and is enzymatically biotinylated using
 Avi-tag™ technology. 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

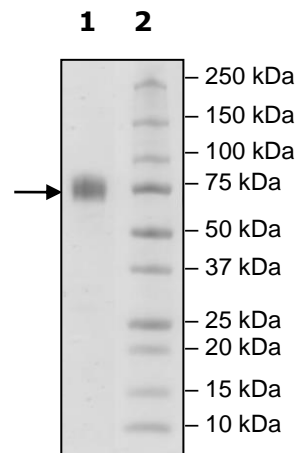
CD48: CD244-Biotin Binding Assay



- no coat
- 100 ng/well CD48

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

Lane 1:
 3μg 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