

## Data Sheet

### **NKp46, Fc-fusion (IgG1), Avi-Tag, Biotin-labeled**

Human, recombinant, C-terminal Fc-Avi-Tag,  
Biotin-labeled

**Catalog #:** 100466

**Lot #:** 190829

**Conc.:** 0.78 mg/ml

**Formulation:** 8 mM Phosphate, pH 7.4, 110 mM NaCl, 2.2 mM KCl, and 20% glycerol.

**Stability:** Stable for at least 6 months at -80°C. Avoid freeze/thaw cycles.

**References:**

1. Mandelboim, O., *et al.*, *Nature*. 2001 Feb; **409**: 1055-1060.
2. Mandelboim, O., *et al.*, *The Int'l J. of Biochem. & Cell Bio.* 2001 Jul; **33(12)**: 1147-1150

**Description:** Human NKp46, also known as CD335, Natural Cytotoxicity Triggering Receptor 1, Lymphocyte Antigen 94 Homolog, and Natural Killer Cell P46-Related Protein, GenBank Accession No. NM\_004829, a.a. 22-258, fused at the C-terminus with the Fc portion of human IgG1 followed by a C-terminal Avi-tag™ and expressed in a HEK293 cell expression system. MW=56 kDa. This protein runs at a higher MW by SDS-PAGE due to glycosylation. This protein is enzymatically biotinylated using Avi-Tag technology. Biotinylation is confirmed to be ≥90%.

### **Quality Assurance**

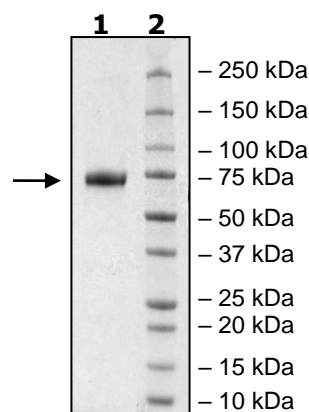
**4-20% SDS-PAGE  
Coomassie staining**

**Lane 1:**  
4µg NKp46

**Lane 2:**  
Protein Marker

**MW:** 56 kDa +glycans

**Purity:** ≥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](mailto:info@bpsbioscience.com)

Please visit our website at: [www.bpsbioscience.com](http://www.bpsbioscience.com)