

## Data Sheet

### Nkp30, Fc Fusion, Avi-tag

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

**Catalog #:** 100546

**Lot #:** 191112

**Conc.:** 1.70 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  $-80^{\circ}\text{C}$ . *Avoid  
freeze/thaw cycles.*

**Description:** Human secreted Nkp30, also known as Natural Cytotoxicity Triggering Receptor 3, NCTR3, Lymphocyte Antigen 117, Activating NK-A1 Receptor, CD337 Antigen, and MALS, GenBank Accession No. NM\_147130, a.a. 19-135, fused at the C-terminus of the Fc portion of human IgG1, with C-terminal Avi-tag™, expressed in a HEK293 cell expression system. MW = 42 kDa.

#### References:

1. Arnon, T.I., *et al.*, *Eu. J. Immunol.* 2001 Sep; **31(9)**:2680-2689.
2. Ferlazzo, G., *et al.*, *JEM.* 2002 Feb; **195(3)**:343

---

## Quality Assurance

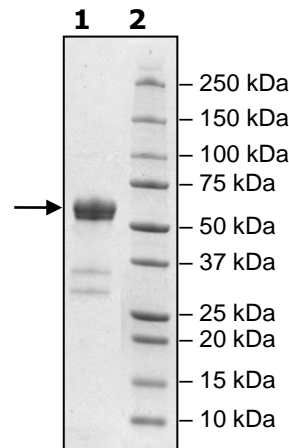
#### 4-20% SDS-PAGE Coomassie staining

**Lane 1:**  
4 $\mu\text{g}$ Nkp30

**Lane 2:**  
Protein Marker

**MW:** 42 kDa

**Purity:**  $\geq 86\%$



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)