

## Data Sheet

### **NKG2D, Avi-Tag, Fc fusion**

Human, recombinant, N-terminal Avi-Fc-Tag  
**Catalog #:** 100252

**Lot #:** 190305      **Conc.:** 0.92 mg/ml

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

**Stability:** Stable for at least 12 months at -80°C. Avoid freeze/thaw cycles.

**Description:** Human NKG2d, also known as, Killer Cell Lectin Like Receptor K1, KLRK1, and NK Cell Receptor D. GenBank Accession No. NM\_007360.4, a.a. 38-216, with an N-terminal Avi-Tag™ fused to the Fc portion of human IgG1 and expressed in a HEK293 cell expression system. MW=53 kDa.

#### **References:**

1. Bauer, S, *et al.*, *Science*. 1999. Jul; **285(5428):**727-729.
2. Wu, J., *et al.*, *Science*. 1999. Jul; **285(5428):** 730-732

---

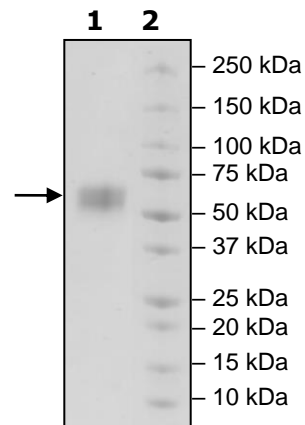
## **Quality Assurance**

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

**Lane 1:**  
4µg NKG2D

**Lane 2:**  
Protein Marker

**MW:** 44 kDa  
**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)