

## Data Sheet

### ULBP1/NKG2DL1, Fc-Fusion (IgG1) Avi-Tag

Human, Recombinant, Fc-Avi-Tag,  
Biotin Labeled

**Catalog #:** 100253

**Lot #:** 181211

**Conc.:** 1.28 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 ULBP1, also known as RAET1I, N2DL-1, NKG2DL1, UL16 and binding protein 1. GenBank Accession No. NM\_025218, a.a. 26-215, 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 = 51 kDa. This protein runs at a higher molecular weight due to glycosylation.

#### References:

1. Rölle, A. *et al.*, *J Immunol.* 2003; **171**(2): 902-908.
2. Kubin, M., *et al.*, *European Journal of Immunology.* 2001; **31**(5): 1428-1437.

### Quality Assurance

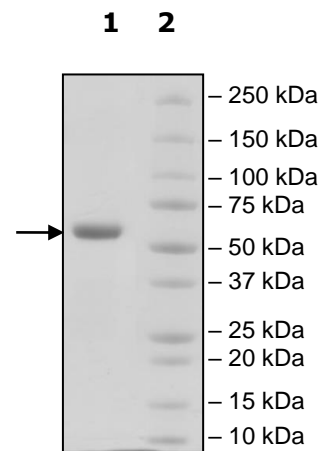
#### 4-20% SDS-PAGE Coomassie staining

**Lane 1:**  
4µg ULBP1

**Lane 2:**  
Protein Marker

**MW:**  
51 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 **1.858.202.1401** 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)