

## Data Sheet

### ULBP2, Avi-His-Tag,

### Biotin-labeled

Human, Recombinant, C-terminal Avi-His-Tag,  
Biotin-labeled

**Catalog #:** 100545

**Lot#:** 191107-1      **Conc.:** 0.83 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.

#### References:

1. Li, K. *et al.*, *Cancer Immunology*.  
2009; **58(5)**: 641-652.
2. Paschen, A. *et al.*, *Clinical Cancer  
Research*. 2009; **15(16)**: 5017-5019.

**Description:** Human ULBP2, also known as UL16 Binding Protein 2, Retinoic Acid Early Transcript 1H, NKG2D Ligand 2, and ALCAN-Alpha, GenBank Accession No. NM\_025217, a.a. 26-216 with a C-terminal Avi-His-Tag, expressed in a HEK293 cell expression system and enzymatically biotin-labeled using Avi-tag™ technology. Biotinylation is confirmed to be  $\geq 90\%$ . MW=25 kDa. This protein runs at a higher molecular weight due to glycosylation.

---

## **Quality Assurance**

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

**Lane 1:**

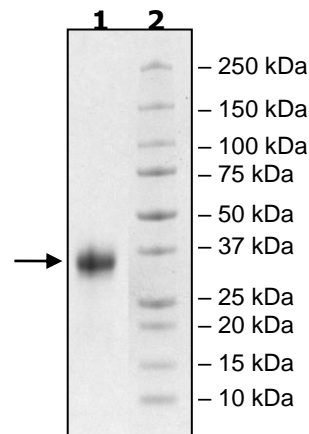
4 $\mu\text{g}$  ULBP2

**Lane 2:**

Protein Marker

**MW:** 25 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)