

## Data Sheet

### ULBP3, Avi-His-Tag

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

**Catalog #:** 100506

**Lot#:** 191119-1

**Conc.:** 0.70 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 ULBP2, also known as UL16-binding protein 3, ALCAN-gamma, and NKG2D ligand 3, GenBank Accession No. NM\_024518.3, a.a. 30-217 with a C-terminal Avi-His-Tag™, expressed in a HEK293 cell expression system. MW=25 kDa. This protein runs at a higher molecular weight due to glycosylation.

#### 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.

---

### Quality Assurance

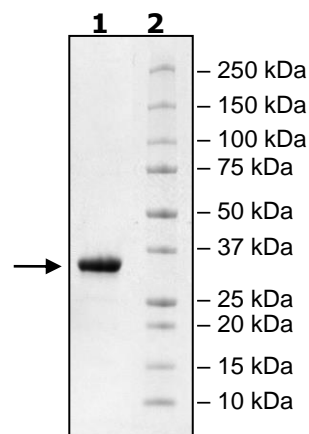
#### 4-20% SDS-PAGE Coomassie staining

**Lane 1:**  
4 $\mu\text{g}$  ULBP3

**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)