

Data Sheet

ULBP2, Avi-His-Tag

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

Catalog #: 100544

Lot#: 191106-1 **Conc.:** 0.60 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°C . Avoid
freeze/thaw cycles.

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 an N-terminal Avi-His-TagTM, 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:

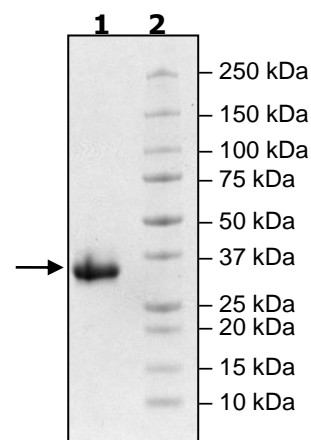
4 μg ULBP2

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

Please visit our website at: www.bpsbioscience.com