

Data Sheet

ULBP1, 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°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

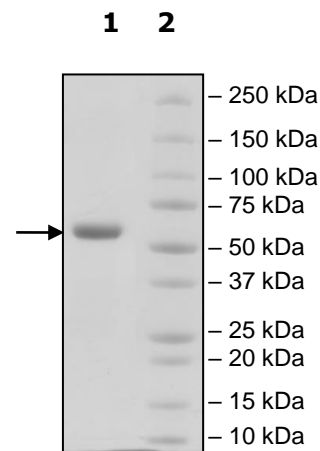
Lane 2:

Protein Marker

MW:

51 kDa + glycans

Purity: ≥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