

## Data Sheet

### CD277, Fc-Fusion (IgG1) Avi-Tag, Biotin-Labeled

Human, Recombinant, Fc fusion protein, with C-terminal Avi-tag™, Biotin-Labeled

**Catalog #:** 100074

**Lot #:** 180510

**Conc.:** 1.22 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. Protein may be diluted to  $\geq 100 \mu\text{g/ml}$  in PBS + glycerol and stored at -80°C.

#### Description:

Human CD227, also known as BTN3A1, with C-terminal Avi-Tag™ fused to the Fc region of Human IgG1. GenBank Accession No. NM\_007048, a.a. 31-254, expressed in a HEK293 cell expression system and enzymatically biotin-labeled using Avi-tag™ technology. MW = 53 kDa. This protein runs at a higher MW by SDS-PAGE due to glycosylation.

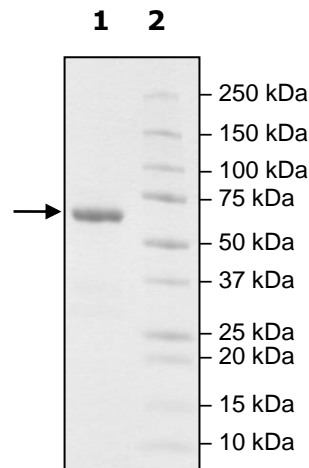
#### References:

1. Wykes, M., et al., *Journal of Leukocyte Biology*, 2002 Oct; **72(4)**: 692-701.
2. Mukherjee, P., et al., *Journal of Leukocyte Biology*. 2004 Jan; **77(1)**: 90-99.

## **Quality Assurance**

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

**Lane 1:**  
4 $\mu\text{g}$  CD277  
**Lane 2:**  
Protein Marker  
**MW:**  
53 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: [support@bpsbioscience.com](mailto:support@bpsbioscience.com)

Please visit our website at: [www.bpsbioscience.com](http://www.bpsbioscience.com)