

## Data Sheet

### CD113, Fc-Fusion, Avi-Tag

Human, Recombinant, C-terminal Fc-Fusion-Avi-Tag

**Catalog #:** 72510

**Lot #:** 170619

**Conc.:** 1.43 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. Protein may be diluted to  $\geq 100 \mu\text{g/ml}$  in PBS + glycerol and stored at  $-80^{\circ}\text{C}$ .

#### References:

1. Bottino, C., *et al.*, *J. Exp. Med.* 2003; 198(4): 557-567.
2. Pende, D., *et al.*, *Blood.* 2005; 105(5): 2066-2073.

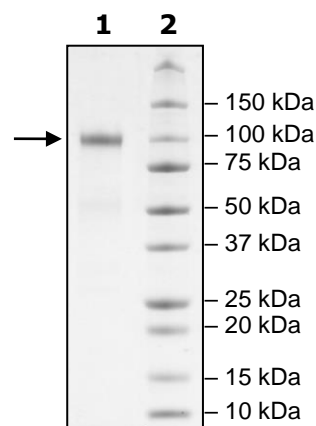
#### Description:

Human secreted CD113, also known as Poliovirus receptor-related 3, PVRL3, and nectin-3, GenBank Accession No. NM\_015480.2, a.a. 58-400 expressed in a HEK293 cell expression system, with C-terminal Fc-Fusion-Avi-tag. MW = 66 kDa. This protein runs at a higher MW by SDS-PAGE due to glycosylation.

## Quality Assurance

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

**Lane 1:**  
2  $\mu\text{g}$  CD113  
**Lane 2:**  
Protein Marker  
**MW:** 66 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 **1.858.829.3082** 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)