

## Data Sheet

### CD96, Fc-Fusion, Avi-Tag

Human, Recombinant, C-terminal Avi-tag

**Catalog #:** 71265

**Lot #:** 170130     **Conc.:** 0.68 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. Hosen, N. *et al.*, *PNAS*. 2007; **104**(26): 11008-11013.
2. Fuchs, A. *et al.*, *Journal of Immunology*. 2004; **172**(7): 3994-3998.

**Description:**

Human CD96, also known as cluster of differentiation 96 and Tactile, GenBank Accession No. NM\_198196, a.a. 22-519. This protein is fused at the C-terminus to the Fc region of Human IgG1 and an Avi-tag™. MW = 84 kDa. This protein runs at a higher MW by SDS-PAGE due to glycosylation.

**Application:**

Useful as a positive control for western blot analysis.

## Quality Assurance

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

**Lane 1:**

3 $\mu\text{g}$  CD96

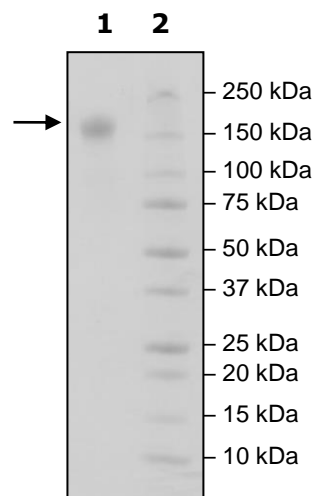
**Lane 2:**

Protein Marker

**MW:**

83.9kDa + 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)