

## Data Sheet

### CD155, Fc-Avi-Tag

Human, recombinant, C-terminal Fc-Avi-Tag  
**Catalog #:** 79063  
**Lot #:** 170601                      **Conc.:** 1.35 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.

#### 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 CD155, also known as Poliovirus receptor and Nectin-like protein 5, GenBank Accession No. NM\_006505, a.a. 27-343, 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 = 63 kDa.

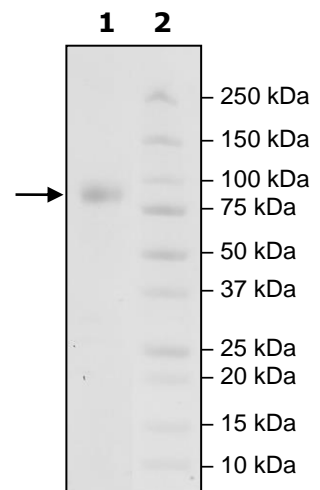
### Quality Assurance

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

**Lane 1:**  
3 $\mu\text{g}$  CD155  
**Lane 2:**  
Protein Marker

**MW:** 63 kDa

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