

## Data Sheet

### Mouse CD155, His-tag, Biotin-labeled

Mouse, recombinant, C-terminal, His-tag, C-terminal Avitag™

**Catalog #:** 71168

**Lot #:** 150521

**Conc.:** 0.63 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.

**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:**

Mouse CD155, also known as Poliovirus receptor and Nectin-like protein 5, GenBank Accession No. NM\_027514, a.a. 29-348 with a C-terminal His-tag and Avitag™. This protein is expressed in a HEK293 cell expression system and enzymatically biotin-labeled using Avitag™ technology. MW = 37.5 kDa. This protein runs at a higher MW by SDS-PAGE due to glycosylation.

**Application:**

Useful for studying protein binding and for screening small molecules and antibodies.

### Quality Assurance

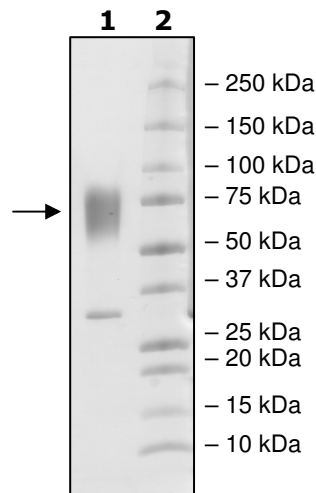
**4-20% SDS-PAGE  
Coomassie staining**

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

**Lane 2:**  
Protein Marker

**MW:** 37.5 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**

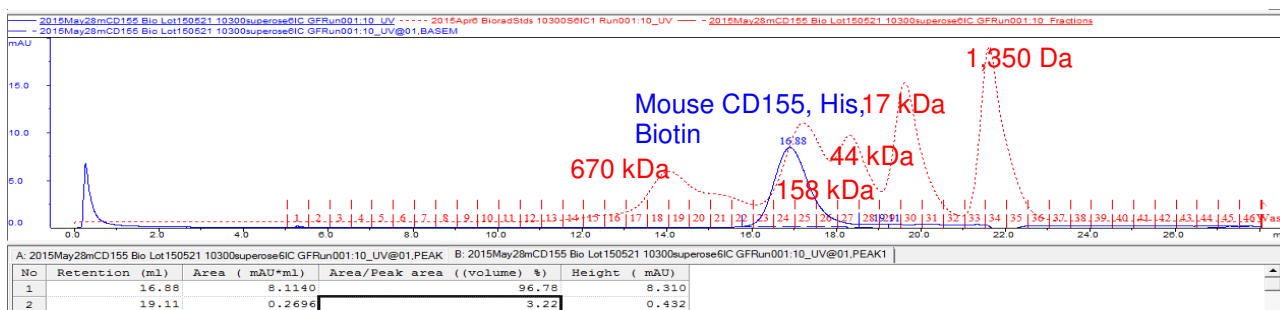
Or you can Email us at: [info@bpsbioscience.com](mailto:info@bpsbioscience.com)

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



6044 Cornerstone Court W, Ste E  
 San Diego, CA 92121  
**Tel:** 1.858.829.3082  
**Fax:** 1.858.481.8694  
**Email:** [info@bpsbioscience.com](mailto:info@bpsbioscience.com)

## Gel Filtration



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)