

## Data Sheet

### SIRP- $\gamma$ (CD172g), Fc fusion, Biotin-labeled

Human, recombinant, Fc fusion with C-terminal Avitag™

**Catalog #:** 71236

**Lot #:** 150717

**Conc.:** 1.49 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. Barclay, AN and Brown, MH., *Nature Rev. Immunol.* 2006; **6**: 457-464.

**Description:** Signal regulatory protein-gamma (SIRP- $\gamma$ ), also known as signal regulatory protein beta 2, and CD172g, GenBank Accession No. NP\_018556, a.a. 29-360, fused at the C-terminus to the Fc portion of human IgG1 and C-terminal Avitag™, MW 67.6 kDa, expressed in a HEK293 expression system and enzymatically biotinylated using Avitag™ technology. Note: SIRP- $\gamma$  is heavily glycosylated, resulting in higher molecular weight. Biotinylation was confirmed to be  $>90\%$ .

**Application:** Useful for studying protein binding and screening small molecules.

## Quality Assurance

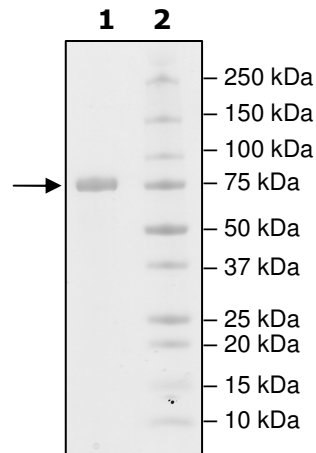
**4-20% SDS-PAGE  
Coomassie staining**

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

**Lane 2:**  
Protein Marker

**MW:** 67.6 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 **1.858.259.5734**

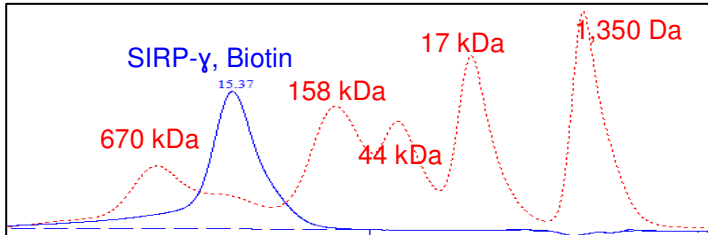
Or you can Email us at: [info@bpsbioscience.com](mailto:info@bpsbioscience.com)

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



6044 Cornerstone Court West, Suite 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 Trace



< 5% aggregated

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 **1.858.259.5734**  
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)