

## Data Sheet

### **SIRP- $\alpha$ (CD172a), Avi-His-Tag, Biotin (Monkey)**

Monkey, Recombinant, C-terminal Avi-His-tag,  
Biotin-Labeled

**Catalog #:** 79303

**Lot #:** 170921      **Conc.:** 1.49 mg/ml

**Formulated in:** 8mM phosphate, 110mM NaCl, 2.2mM KCl, pH7.4, 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. Hatherley, D., *et al.*, *Mol. Cell*, 2008, **31**:266-77
2. Baba, N., *et al.*, *J. Exp. Med.*, 2013, **210(6)**:1251-63

**Description:** Monkey Signal regulatory protein-alpha (SIRP- $\alpha$ ), also known as Tyrosine-protein phosphatase non-receptor type substrate 1, CD172a, Bit, and MyD-1, GenBank Accession No. NM\_001284750, a.a. 31-370, with C-terminal Avi-His-tag, expressed in a HEK293 expression system. MW=42 kDa. The protein is enzymatically biotinylated using AviTag™ technology. This protein is heavily glycosylated, resulting in apparent higher molecular weight.

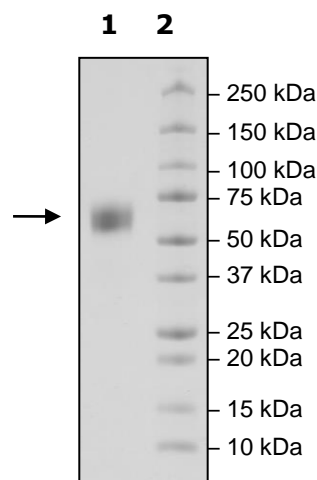
---

### **Quality Assurance**

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

**Lane 1:**  
4 $\mu\text{g}$  Sirpa(Monkey)  
**Lane 2:**  
Protein Marker

**MW:** 42 kDa + glycans  
**Purity:**  $\geq 90\%$



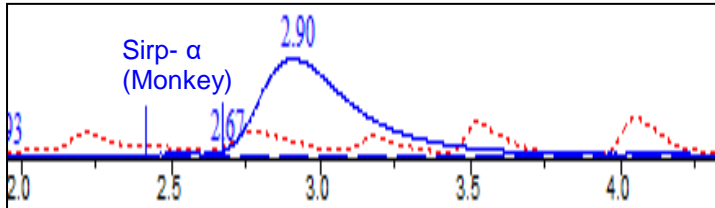
OUR PRODUCTS ARE FOR RESEARCH USE ONLY. NOT FOR DIAGNOSTIC OR THERAPEUTIC USE.

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)



6042 Cornerstone Court West, Suite B  
San Diego, CA 92121  
Tel: 1.858.829.3082  
Fax: 1.858.481.8694  
Email: [info@bpsbioscience.com](mailto:info@bpsbioscience.com)

### Gel Filtration Curve



<10% Aggregation

### Protein Sequence

EEELQVIQPEKSVSVAAGDSATLNCTVSSLI PVGPIQWFRGAGPGRELIYNLKEGH  
FPRVTAVSDPTKRNNMDFSIRISNITPADAGTY YCVKFRKGS PDVELKSGAGTELS  
VRAKPSAPVVSGPAVRATAEHTVSFTCESHGFS PRDITLKWFKNGNELSDVQTNVD  
PAGKSVSYSIRSTARVLLTRRDVHSQV ICEVAHVTLQGDPLRGTANLSEAIRVPPF  
LEVTQQSMRADNQVNVTCQVTKFY PQRQLQLTWLENGNVS RTEMASALPENKDGTYN  
WTSWLLVNVSAHRDDVKLTCQVEHDGQ PAVNKSFSVKVSAHPKEQGSNTAAENTGT  
NERNIEGRMDGGGLNDIFEAQKIEWHEHHHHHHHHHH

Black: Monkey Sirpa (aa 31-370)

Green: Linker

Red: Avi

Blue: 10xHis

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)