

## Data Sheet

### SLAMF7, Avi-His-Tag, Biotin-Labeled

Human, Recombinant, C-terminal Avi-His-Tag, Biotin-Labeled

**Catalog #:** 100272

**Lot#:** 190214-1

**Conc.:** 1.12 mg/ml

**Description:** Human SLAMF7, also known as 19A, CD319, CRACC, CS1, and SLAM family member 7, GenBank Accession No. NM\_021181.5, a.a. 23-226 with C-terminal Avi-His-Tag, expressed in a HEK293 expression system. This protein is enzymatically biotinylated using Avi-Tag™ technology. Biotinylation is confirmed to be >90%. MW = 25 kDa.

**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.

#### References:

1. J. van de Donk, N. W. C. *et al.*, *Blood*. 2016; **127**: 681-695.
2. Huaijian, G. *et al.*, *MCB*. 2015; **35(1)**: 41-51.

---

## Quality Assurance

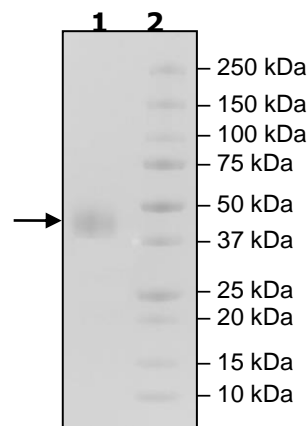
### 4-20% SDS-PAGE Coomassie staining

**Lane 1:**  
4 µg SLAMF7

**Lane 2:**  
Protein Marker

**MW:**  
25 kDa+glycans

**Purity:** ≥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**

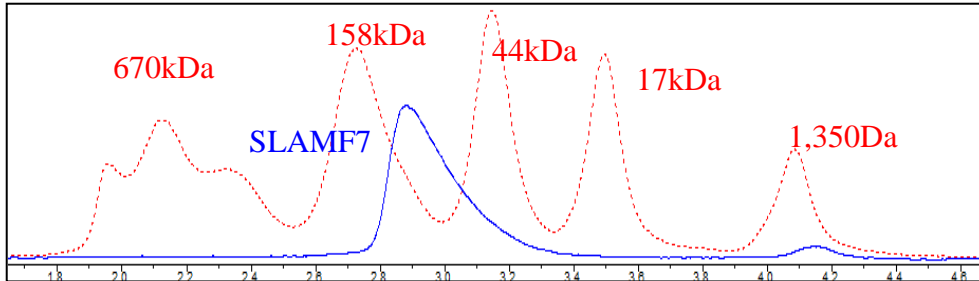
Or you can Email us at: [info@bpsbioscience.com](mailto:info@bpsbioscience.com)

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



6042 Cornerstone Court West, Suite B  
San Diego, CA 92121  
Tel: 1.858.202.1401  
Fax: 1.858.481.8694  
Email: [info@bpsbioscience.com](mailto:info@bpsbioscience.com)

**Gel Filtration Curve**



<10% Aggregation

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)