

## Data Sheet

### Siglec7, Fc Fusion, Avi-tag, Biotin-labeled

Human, Recombinant protein,  
C-terminal Fc-Avi-Tag, Biotin-labeled.

**Catalog #:** 100636

**Lot #:** 200227

**Conc.:** 0.42 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. Jandus, C. *et al.*, *JCI*. 2014;  
**124(4)**:1810-1820.
2. Orr, S.J. *et al.*, *JBC*. 2007; **282**:3418-  
3422.

**Description:** Human secreted Sialic Acid Binding Ig Like Lectin 7 (Siglec 7), also known as AIRM-1 and CD328, GenBank Accession No. NM\_014385, a.a. 19-353, fused at the C-terminus of the Fc portion of human IgG1, with C-terminal Avi-tag™, expressed in a HEK293 cell expression system and enzymatically biotinylated using Avi-tag™ technology. Biotinylation is confirmed to be  $\geq 90\%$ . MW = 66 kDa. This protein runs at a higher molecular weight due to glycosylation.

## Quality Assurance

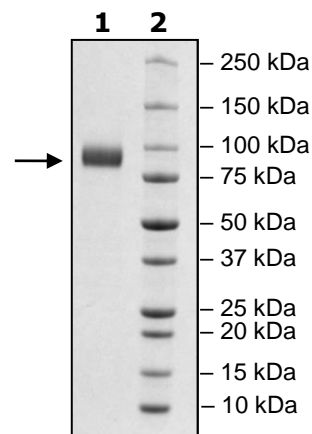
### 4-20% SDS-PAGE Coomassie staining

**Lane 1:**  
4 $\mu\text{g}$  Siglec7

**Lane 2:**  
Protein Marker

**MW:** 66 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 **1.858.202.1401** 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)