

Data Sheet

Siglec7, Fc Fusion, Avi-tag

Human, Recombinant protein,
C-terminal Fc-Avi-Tag

Catalog #: 100589

Lot #: 200211

Conc.: 0.47 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.

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. MW = 66 kDa. This protein runs at a higher molecular weight due to glycosylation.

Quality Assurance

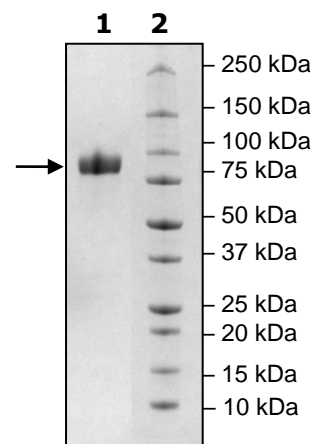
4-20% SDS-PAGE Coomassie staining

Lane 1:
4 μg 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

Please visit our website at: www.bpsbioscience.com