

Data Sheet

Siglec9, Fc Fusion, Avi-tag, Biotin-labeled

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

Catalog #: 100559

Lot #: 200213

Conc.: 0.65 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. Ikehara, Y., *et al.*, *JBC*. 2004 Oct 8;**279**:43117-43125.
2. Zhiang, J.Q., *et al.*, *JBC*. 2000 Jul 21;**275**:22121-22126

Description: Human secreted Sialic Acid Binding Ig Like Lectin 9 (Siglec 9), also known as CD329, CDw329, FOAP-9, and OBBP-LIKE, GenBank Accession No. NM_014441.2, a.a. 18-348, 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

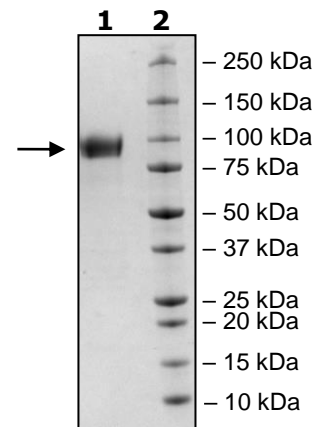
**4-20% SDS-PAGE
Coomassie staining**

Lane 1:
4 μg Siglec9

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