

Data Sheet

Siglec15, Fc Fusion Avi-tag

Human, Recombinant protein,
C-terminal Avi-tag

Catalog #: 100423

Lot #: 190710

Conc.: 0.97 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.

Description: Human secreted Sialic Acid Binding Ig Like Lectin 15 (Siglec15)-Fc fusion protein, also known as CD33 Antigen-Like 3 (CD33L3) and HsT1361, GenBank Accession No. NM_213602, a.a. 20-263, 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 = 55 kDa.

References:

1. Angata, T., *et al.*, *Glycobiology*. 2007 May 4;**17(8)**:838-846.
2. Hiruma, Y., *et al.*, *BBRC*. 2011 Jun 10;**409(3)**:424-429

Quality Assurance

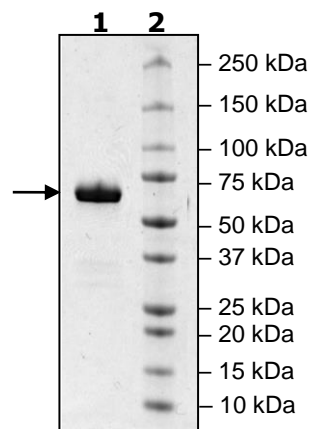
4-20% SDS-PAGE Coomassie staining

Lane 1:
4μg Siglec15

Lane 2:
Protein Marker

MW: 55 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

Please visit our website at: www.bpsbioscience.com