

Data Sheet

CD33, Fc Fusion, Avi-tag

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

Catalog #: 100155

Lot #: 190724

Conc.: 2.23 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 CD33, also known as Sialic Acid Binding Ig Like Lectin 3, Gp67 and Myeloid Cell Surface Antigen CD33, GenBank Accession No. NM_001772, a.a. 1-259, 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 = 57 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

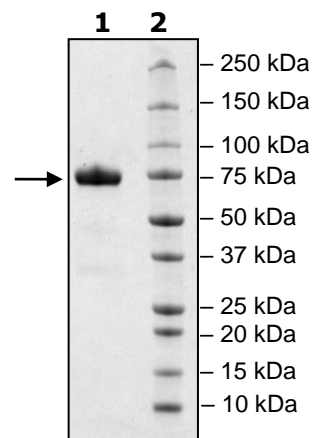
4-20% SDS-PAGE Coomassie staining

Lane 1:
4 μg CD33

Lane 2:
Protein Marker

MW: 57 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