

## Data Sheet

### CD33, Fc Fusion, Avi-tag, Biotin Labeled

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

**Catalog #:** 100482

**Lot #:** 190815                      **Conc.:** 0.82 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. 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

**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 with 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. MW = 57 kDa. Biotinylation is confirmed to be  $\geq 90\%$ .

## Quality Assurance

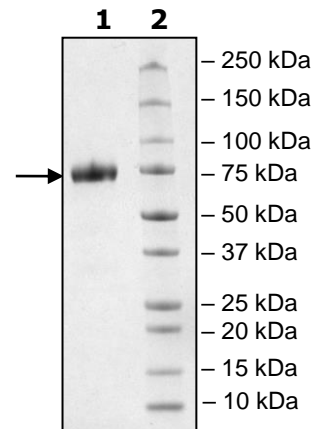
#### 4-20% SDS-PAGE Coomassie staining

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

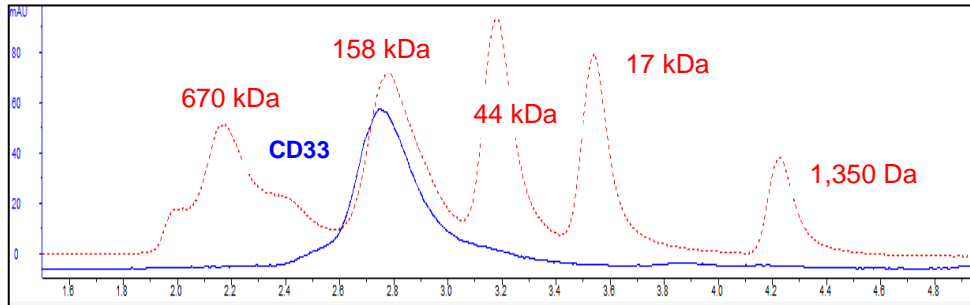
Or you can Email us at: [info@bpsbioscience.com](mailto:info@bpsbioscience.com)

Please visit our website at: [www.bpsbioscience.com](http://www.bpsbioscience.com)



6042 Cornerstone Court W, Ste B  
San Diego, CA 92121  
**Tel:** 1.858.202.1401  
**Fax:** 1.858.481.8694  
**Email:** info@bpsbioscience.com

**Gel Filtration Curve:**



<10%  
Aggregation

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)