

## Data Sheet

### VSIG3, Fc-fusion (IgG1), Avi-Tag

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

**Catalog #:** 79491

**Lot #:** 180410

**Conc.:** 1.14 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. Protein may be diluted to  
 $\geq 100 \mu\text{g/ml}$  in PBS + glycerol and stored at  
 $-80^{\circ}\text{C}$ .

**References:**

1. Kato, L., *et al.*, *International Journal of Oncology*. 2003; **23**: 525-531.
2. Watanbe, T., *et al.*, *Cancer Science*. 2005; **10**: 17858-17863.

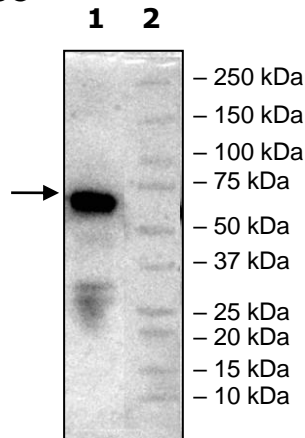
**Description:**

Human VSIG3, also known as V-set and immunoglobulin domain containing 3, IgSF11, GenBank Accession No. NM\_152538.3, a.a. 22-244, with C-terminal Avi-Tag™ fused to Fc region of Human IgG1, expressed in a HEK293 cell expression system. MW = 52 kDa. This protein runs at a higher MW by SDS-PAGE due to glycosylation.

### Quality Assurance

**Western Blot**

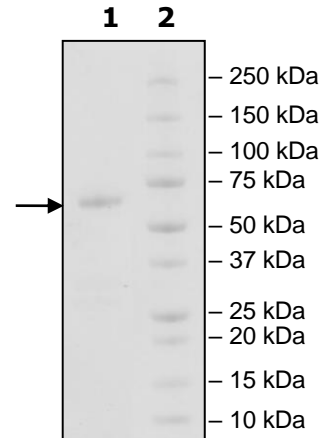
**Lane 1:**  
 3 $\mu\text{g}$  VSIG3  
**Lane 2:**  
 Protein Marker



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

**Lane 1:**  
 4 $\mu\text{g}$  VSIG3  
**Lane 2:**  
 Protein Marker

**MW:** 52 kDa  
**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: [support@bpsbioscience.com](mailto:support@bpsbioscience.com)  
 Please visit our website at: [www.bpsbioscience.com](http://www.bpsbioscience.com)